



Smart-Vue Pro Monitoring Solution

Web application for
Smart-Vue Pro Duo/Quattro Data Loggers

User Guide

331676H01 • Revision A • 5/27/2020

IMPORTANT Read this user guide. Failure to follow the instructions in this user guide can result in damage to the unit, injury to operating personnel and poor equipment performance.

CAUTION All internal adjustments and maintenance must be performed by qualified service personnel.

Material in this manual is for informational purposes only. The contents and the product it describes are subject to change without notice. Thermo Fisher Scientific makes no representations or warranties with respect to this manual. In no event shall Thermo be held liable for any damages, direct or incidental, arising from or related to the use of this manual.

© 2020 Thermo Scientific. All rights reserved. Thermo Scientific, the Thermo Scientific logo, Smart-View Pro Duo and Smart-View Pro Quatro are the exclusive property of Thermo Scientific. All other brands mentioned are the property of their respective owners. This is a non-contractual document. Features and specifications mentioned here are subject to change.

Contents

Safety Instructions	1
Introduction	2
About this Manual	2
Overview of Smart-Vue Pro	2
Getting Started	5
Connecting to Smart-Vue Pro	5
Creating your Smart-Vue Pro Company Account	6
Forgot your password?	8
Using Smart-Vue Pro	9
User Interface.....	9
Home Screen Overview	12
Plan View	12
Watch Mode	12
Users	15
Authentication Modes	15
Adding a New User	17
Editing a User.....	18
User Access Profiles (Roles)	22
Your Organization	24
Company Information	24
Organization with Sites and Departments	25
Adding Plan Images	28
Infrastructure Devices	30
Adding Infrastructure.....	30

Editing Infrastructure	31
Adding a LoRaWAN Receiver as Infrastructure	31
Data Loggers	32
Adding Data Loggers	32
Viewing the Data Logger List	34
Viewing Data Logger Details	34
Data Logger Actions	35
Editing Data Logger Details	36
Sensors	37
Viewing the Sensor List	37
Viewing Sensor Details	37
Calibration Parameters	38
4-20 mA, 0-5 V, 0-10 V Loop Sensors	39
Simulating a Sensor Alarm for Testing	40
Removing a Sensor	40
Disabling / Enabling a Sensor	41
Equipment	42
Adding Equipment	42
Viewing Equipment Details	44
Modifying / Managing Equipment	45
Deleting Equipment	45
Data Logging	46
How does it work?	46
Starting Data Logging	46
Stopping Data Logging	48
Viewing Sensor Data	49
Displaying Sensor Graph	49
List of Sensor Readings	51
Creating Reports	51

Alert Devices	53
Smart Siren	53
Smart Remote Contact	55
Alarms & Alerts	59
Handling Alarms	59
Acknowledging Alarms	60
Alarm Report Subscriptions	61
Alert Notifications	63
Alert Rules with Users	64
Time Slots	66
Getting Help	68
User Manuals	68
Viewing Smart-Vue Pro Version Release Notes	68
Appendix	69
Alarms	69
FAQ	69
WEEE Compliance	71
Contact Information	72

Safety Instructions



IMPORTANT NOTE: Do not use this product for protection or as part of an automated emergency system or as for any other application that involves protecting people and/or property. This product is designed for use in environments where children are not likely to be present. Customers and users of Thermo Scientific products are responsible for making sure that the product is fit for the intended usage. Do not open the product casing and do not disassemble or modify internal components in any manner. Thermo Scientific products do not contain any internal components that require user intervention or repair. If the device shows signs of improper operation, disconnect it immediately from its power source or remove the battery and contact Thermo Scientific technical services.

Introduction

This user guide describes how to use the Smart-Vue Pro web application to manage and monitor Smart-Vue Pro data logger featuring LoRaWAN™ long-range wireless connectivity.

Smart-Vue Pro data loggers collect data automatically from connected sensors and transmit the data to the Cloud. These data loggers are part of Thermo Scientific's latest generation of wireless IoT (Internet of Things) devices.

About this Manual

This user guide describes how to use the features offered by the Smart-Vue Pro application. Before using Smart-Vue Pro, LoRaWAN™ connectivity for your devices must be configured to use a private (on-premises) network with a Thermo Scientific LoRaWAN™ receiver. See Thermo Scientific LoRaWAN receiver user guide for more details.

Table 1. Terminology used in this doc.

Terminology	Description
Alarm	An alarm occurs when the system observes a sensor reading that is out-of-bounds, such as a temperature reading or humidity level or 4-20 mA / 0-5 V / 0-10 V reading that is too high or too low with respect to programmed range limits. The system can notify users when alarms occur by sending alerts.
Alert	An alert is a notification sent by the system to users when the system observes an alarm condition or potential problem. Alerts can be sent by e-mail, SMS or voice calls.
Equipment	The material or space that you are monitoring with a data logger.
LoRaWAN	Very-long-range wireless communication protocol available with public or private network connectivity. Wireless range can reach up to nearly 10 miles (16 km).
Cloud	Internet-based platform on which data from data loggers is stored and accessed via the Smart-Vue Pro application.

Overview of Smart-Vue Pro

The concept of Smart-Vue Pro is based on managing sensors that are assigned to entities called “equipment”. Equipment refers to any assets (such as refrigerators and freezers), rooms or locations in/on which sensors may be placed.

The overall process of using Smart-Vue Pro Web is divided into the following steps:

1. You must first order a license voucher to cover the number of sensors you will be using. Contact the Thermo Scientific sales representative for more information.
2. Log in to Smart-Vue Pro Web to create your company account, add users and set up your equipment.

The process for configuring Smart-Vue Pro is shown in the following figure:

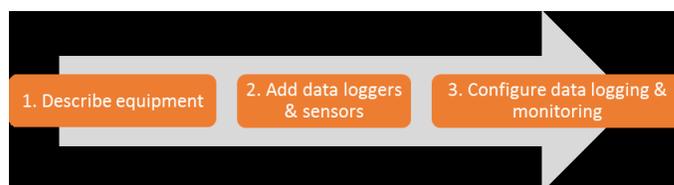


Figure 1. Process Overview

1. **Describe equipment:** The notion of equipment is the central principle for sensor monitoring with Smart-Vue Pro. This approach is based on the idea that you are more likely concerned about the assets you are monitoring than the tool used to accomplish the task. For example, you monitor equipment such as a cold room, refrigerator, deep freezer or incubator. You therefore must add this equipment to your system.
2. **Add data loggers and sensors:** Equipment is monitored by one or more data loggers with sensors that record temperature or humidity or 4-20 mA / 0-5 V / 0-10 V etc. at regular intervals. Sensors are added automatically when the data logger to which they are attached is added to the system.
3. **Configure data logging and monitoring:** As different pieces of equipment may have different needs in terms of monitoring, sensors must be configured to monitor different ranges at different reading frequencies with different alert strategies, etc.

Smart-Vue Pro Feature Highlights

Smart-Vue Pro enables you to visualize your Smart-Vue Pro data logger's sensor status and data with a management interface for configuration and system administration.

Key application features include:

User and license management

- Creating company and user accounts.
- Managing users, profiles and access rights.
- Managing activation keys (license voucher).
- Managing your company's different sites and departments.
- Compatibility with FDA 21 CFR Part 11 requirements.
 - Readings and settings stored in secure database
 - Password protection

Data

- Configuring sensors, receivers and infrastructure elements.
- Viewing data recorded by Thermo Scientific data loggers.
- Monitoring equipment and associated sensors.
- Managing data transmitted by Thermo Scientific Smart-Vue Pro data loggers.
- Viewing detailed sensor readings, history and more.
- Managing sensor calibrations and calibration reminders.

Alarms

- Configuring and sending alerts when alarms are detected.
- Managing various types of alarms: upper and lower limits, technical alarms in case of a sensor problem or wireless communication problem and more.
- Adapting alarm handling for days, nights and weekends according to user level and desired contact methods.

Sensor reading

- Viewing detailed sensor readings, history and more.
- Displaying sensor graphs.
- Creating reports for all key parameters, including configuration, status, events and alarms.

Prerequisites

Smart-Vue Pro is a web application designed to work with standard web browsers.

The minimum system requirements are as follows:

- Personal computer
- Internet connection
- Screen resolution 1280 x 900 pixels minimum
- Supported browsers (indicated version or higher):
 - Google Chrome 66
 - Mozilla Firefox 60
 - Microsoft Edge version 42
 - Safari 9 (iOS 13 and above)
 - Opera 53

Note: The Smart-Vue Pro web application does not support Internet Explorer.

Cloud Access Requirements

When using Smart-Vue Pro in Cloud mode, the following ports must be open:

Table 2. Ports

Purpose	URL	Ports	Protocol
For users: Smart-Vue Pro web application	smart-vuepro.thermoscientific.com	443	https
For infrastructure: LoRaWAN receiver connection	smartvueconnect.com	1700	UDP

These ports are used for two-way communication between the Smart-Vue Pro web application running in your browser and the server.

Getting Started

This section describes how to:

1. Create an account on the Smart-Vue Pro web application.
2. Login to Smart-Vue Pro.
3. Create equipment and add a data logger to monitor it.
4. Adjust monitoring settings and complete the data logging configuration.



CAUTION: This section assumes that you have purchased a license key that is available. Contact the Thermo Scientific sales representative for more information.

Connecting to Smart-Vue Pro

You must connect to the Smart-Vue Pro web platform to use the software.

1. Launch your internet browser application.
2. Enter `smart-vuepro.thermoscientific.com` in the address bar.

The application detects your browser language automatically and displays the login window either in English, French, German, Italian, Spanish or Portuguese. For all other languages, English is used:

Figure 2. Authentication window (login)

Login Authentication

Smart-Vue Pro supports several authentication methods depending on the type of solutions used. The application can either use its own user management platform or it can connect to an external authorization platform (on-premises solutions only).

To use the integrated Smart-Vue Pro authentication:

1. Enter your assigned user-name and password.
2. Click **Log in** to enter the application.

Passwords are case sensitive and encrypted. If you forget your password, click **Forgot password**.

As a security measure, the account is locked if you enter an incorrect password for three consecutive times. To unlock the account, reset the password. Proceed with the **Forgot password** procedure or contact an Application Manager to reset the password (which you must then complete).

To use an external authorization platform (such as LDAP) for Smart-Vue Pro Server (on-premise server) solution:

1. Enter your credentials in the form “authentication_mode/username” and enter the password.
2. Click **Log in** to enter the application.



CAUTION: For specific information on Authentication modes, see **Authentication Modes**. For more information on how to create user accounts, see **Users**.

Changing the Display Language:

The language used for Smart-Vue Pro depends on your browser configuration.

Once the Smart-Vue Pro is connected, you can change the application language at any time:

1. Click on your profile image (avatar) at the top right-hand corner of the screen.
2. Click **Profile** → **Options(3 vertical dots)** → **Edit localization settings** .
3. Click on the **Language** and select the desired language from the list. For the language change to be effective, you must logout and log back into the system.

By default, the user session will disconnect from Smart-Vue Pro automatically after one hour of inactivity. To reconnect to the platform, enter the login credentials in the authentication screen.



CAUTION: The Smart-Vue Pro application version is shown at the bottom of the login window. Move your mouse cursor over the version number for more information. This information may be requested from the technical support team.

Creating your Smart-Vue Pro Company Account

You must create a Smart Vue Pro account to interact and manage Smart-Vue Pro data loggers. You may add users to an existing account at any time (see **Authentication Modes** for details).



CAUTION: When you first create your company account in Smart-Vue Pro, you also create the first user. The user has Application Manager rights and can create additional users. See **Users** for more information.

To create your company account and the first Application Manager user:

1. Enter `smart-vuepro.thermoscientific.com` in the address bar of your web browser.
2. Click **New company?** as shown in Figure 3.

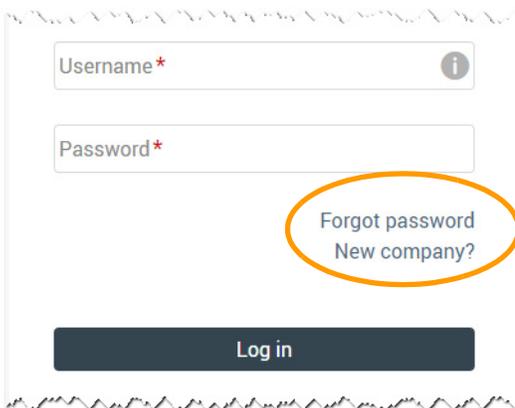


Figure 3. Authentication Window (login)

Company Information

Smart-Vue Pro includes a 2-step setup wizard for creating your company account and a first user profile. The first user profile created has Application Manager rights and can configure the entire system and add other users. The fields marked with * are required.

Figure 4. Filling in the company information

Fill the required fields (marked with *) as appropriate. You can change or update this information later if necessary:

Company name: Enter the name of your company.

Platform: Unique company identifier, filled in automatically based on the company name you enter.

Company address: City, Postal code, Country, International phone number.



CAUTION: Make sure you enter the telephone number in international format, with the prefix “[country code]”, such as: +14153817894. Do not include any extra digits or leading zeros.

The following settings apply as a default for your system. They can be overridden by each individual’s settings for local operation and display purposes, but this information is used

as a common reference for the company in case people have conflicting profile settings for shared actions.

Example: If you have e-mail alerts sent to people in different regions, English could be the default company language for the same alert message.

Language: Choose the desired default language for your company. You are free to choose your personal preference at any time.

Time zone: Select the time zone based on the primary geographical area for the company.

Date format: Choose the date format to be displayed in the application.

Hour format: Choose the time format to be displayed in the application. Readings can be displayed in the **12-hour (AM/PM)** or **24-hour** format.

Unit: Choose the unit of temperature (**degree celsius (°C)** or **Fahrenheit (°F)**).

Decimal separator: Select the character to use as a decimal separator in numerical values.

License key: Enter your license key in this field. Based on the number of measurement points (sensors) in your system, this key is provided when you subscribe to Smart-Vue Pro.

After completing all the company information, click **Next** to fill in the user information as described in the next section.

User Information

Enter information for the first user.



CAUTION: The e-mail address you enter here is the only information that you cannot change subsequently for your company account. We recommend using an e-mail address that is not strictly tied to a single person, such as “smart-vuepro@your_company.name”.

Figure 5. Entering user information

Fill in the required fields (marked with *) as appropriate:

First name / Last name: Enter the first name and last name.

Email: Enter the e-mail address that is used to login to the system. The e-mail alerts and confirmations by the system are sent to this email address. The e-mail address may only be used once into the system and may not be edited once entered.

Job title: Enter the job function within the company.

After completing the information, select the checkbox “**I accept the Terms of Use and the Privacy Policy Terms**” and click **Proceed**.

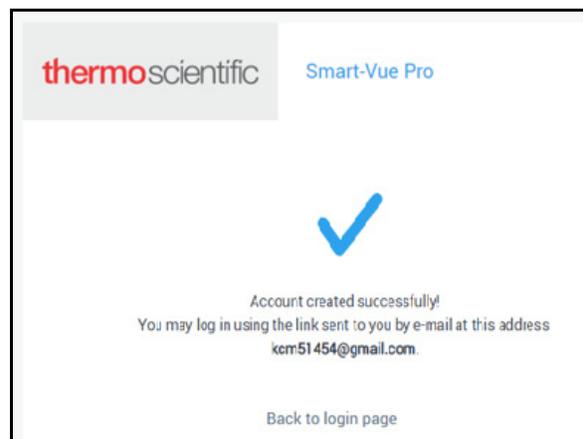


Figure 6. Smart-Vue Pro confirmation of account creation

Your account is created and a confirmation e-mail is sent to the address entered. Click on the link in the e-mail to activate the account.

First Connection to the Application



CAUTION: This section applies to systems using Smart-Vue Pro's integrated authentication (based on the user-name and password). For external authentication platforms, notably LDAP, use your regular corporate login credentials.

If your system uses Smart-Vue Pro's integrated authentication mode, you can change the initial password the first time you login to the application.

1. Check the confirmation e-mail in your inbox and click on the link provided. If you do not see the e-mail, check the spam or junk e-mail folders in case your application filtered it automatically.
2. After you click on the link, the following window opens to set your password:

Figure 7. Changing a password upon first connection

3. Enter a new password. Passwords must be at least 8 characters long and contain a combination of upper- and lower-case letters with at least one number and one special character.
4. Confirm the new password by re-typing in the field below.
5. Click **OK**. The password is reset successfully.

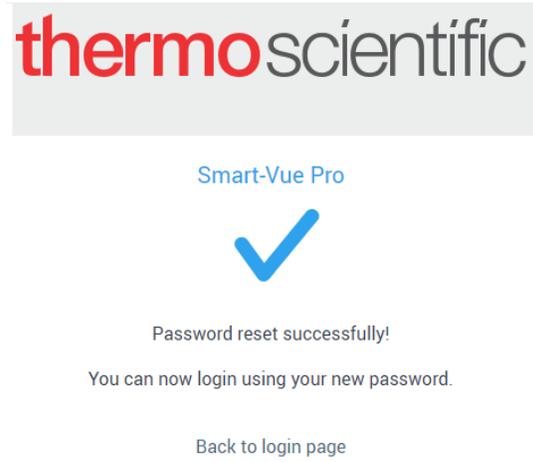


Figure 8. Successful password change

6. Click **Back to login page** to connect to the application using your e-mail address and the password you created.

Forgot your password?



CAUTION: This section describes password recovery for systems using Smart-Vue Pro's integrated authentication (based on the user's e-mail address and password). For external authentication platforms, notably LDAP, check with your system administrator to recover or reset a forgotten password.

If you are unable to access Smart-Vue Pro or if you cannot remember your user-name or password, the system can send password reset instructions to you by e-mail. Follow these instructions if you must reset your Smart-Vue Pro password:

1. Use your browser to open the address: smart-vuepro.thermoscientific.com
2. Click **Forgot password** on the login screen.
3. Enter your Smart-Vue Pro account email address and click **OK**.

Figure 9. Entering an e-mail address to receive a new password

4. Check your e-mail in-box and click on the link in the email you received to set a new password.

Using Smart-Vue Pro

User Interface

The Smart-Vue Pro interface consists of the following main zones:

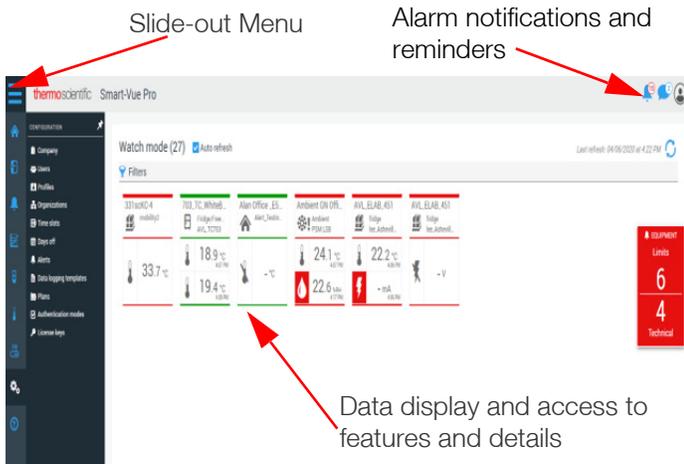


Figure 10. Overview of Smart-Vue Pro User Interface

The watch mode screen shown above provides direct access to key features in your monitoring system.

Main Menu

The main menu is located on the left-hand side of the screen that slides open or close when you hover the mouse on the menu icon . Click Menu to access Smart-Vue Pro features. Depending on the context, the menu item displays a statistics pane or a sub-menu with additional features or information (See **Figure 11**):

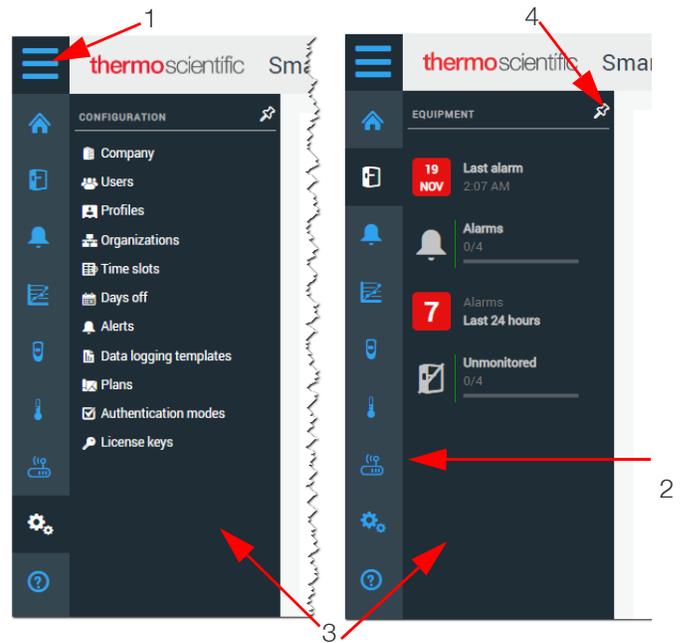


Figure 11. General layout of Smart-Vue Pro Main Menu

Table 3. Main menu features

S.No	Description
1	Menu slide open
2	Navigation menu:  Home  Equipment  Alarms  Reports  Data Logger  Sensors  Infrastructure  Configuration  Help

Table 3. (Cont.)Main menu features

S.No	Description
3	Sub-menu of additional features or Statistics pane depending on the context.
4	Toggle the thumb tack to Pin or Unpin the menu (keep it open or allow it to slide back automatically).

The background color changes as you move your mouse over each feature and when you select a feature.

Sub-Menu Indicators

Your system constantly monitors the health of all its components to take an action if necessary. As you slide the main menu on the left-hand side of the screen, sub-menus with key indicators are displayed. Some icon colors vary according to status, such as alarms.

Equipment: This information pertains specifically to the status of your equipment.

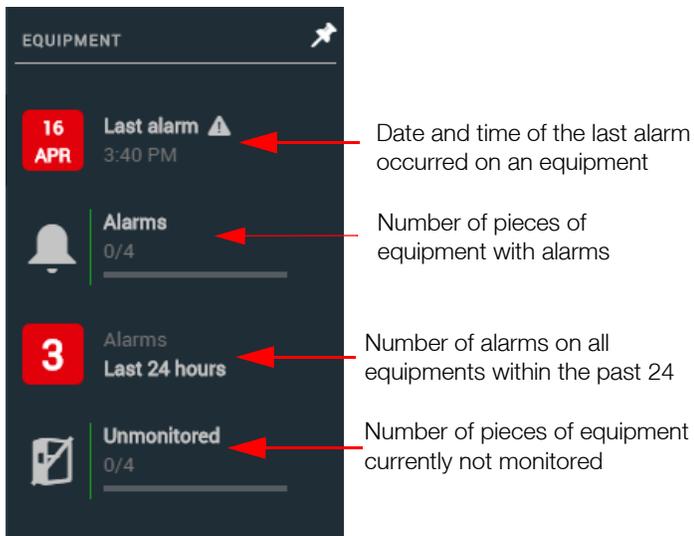


Figure 12. Equipment

Alarms: This information pertains specifically to alarms detected by the system.

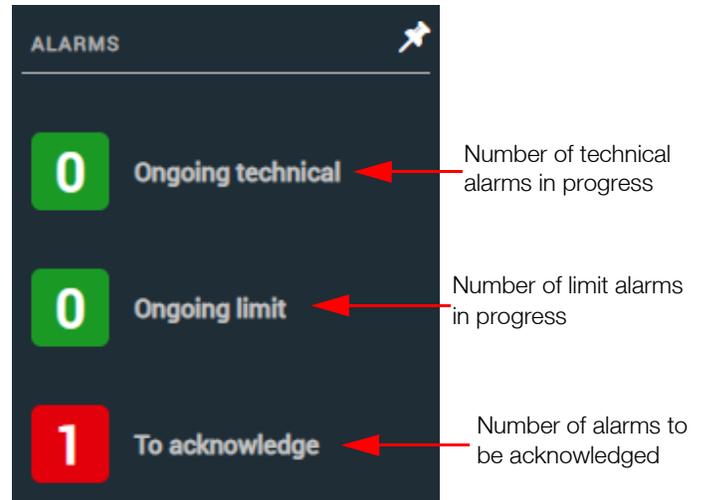


Figure 13. Alarms sub-menu indicators

Data loggers: This information pertains specifically to the status of your data loggers (i.e. not their sensors or equipment to which they are assigned).

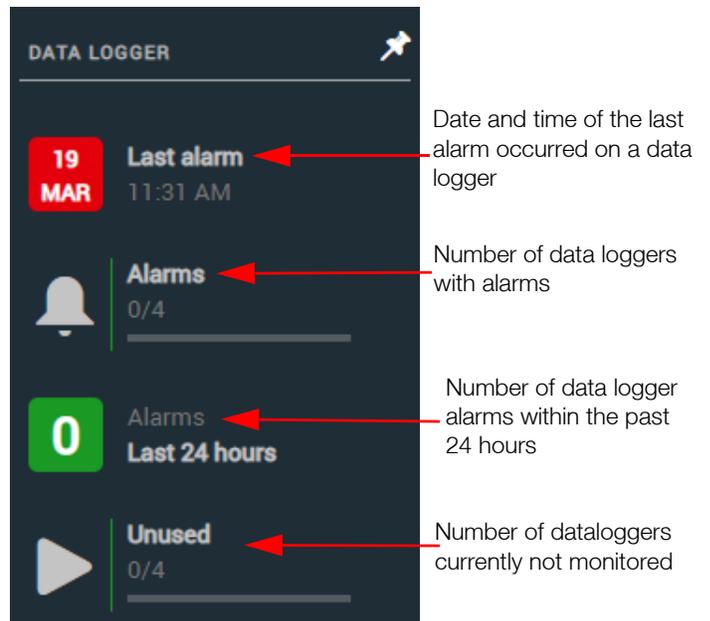


Figure 14. Data logger sub-menu indicators

Sensors: This information pertains specifically to the status of the sensors on your data loggers.

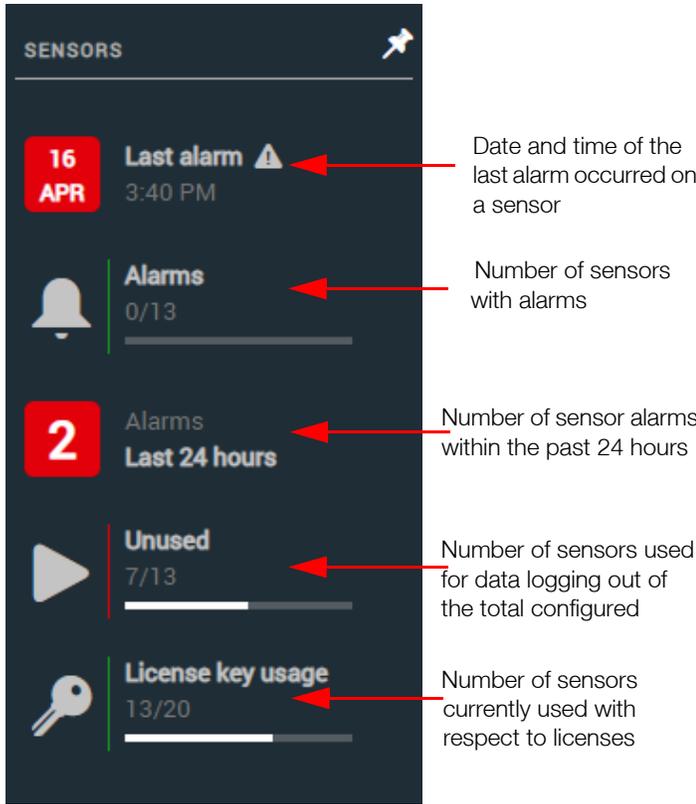


Figure 15. Sensor sub-menu indicators

Infrastructure: This information pertains specifically to the status of your infrastructure devices, such as receivers and smart remote contact or siren alert devices.

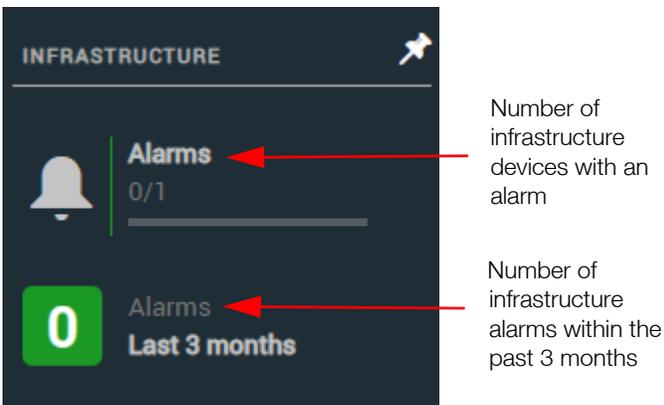


Figure 16. Infrastructure

Using Filters

Many Smart-Vue Pro screens include a **Filters** option at the top of the display. You may use various fields and search criteria to adapt lists displayed on screen or sort data in different ways. Filters can be very useful for finding specific information quickly or limiting long lists.

The **Filters** feature always works as described depending on the context.

To use **Filters**:

1. Click **Filters** (1) and use the drop-down menu to choose the information you want to display. The following examples show the filters options on the **Equipment** and **Sensors** screens:

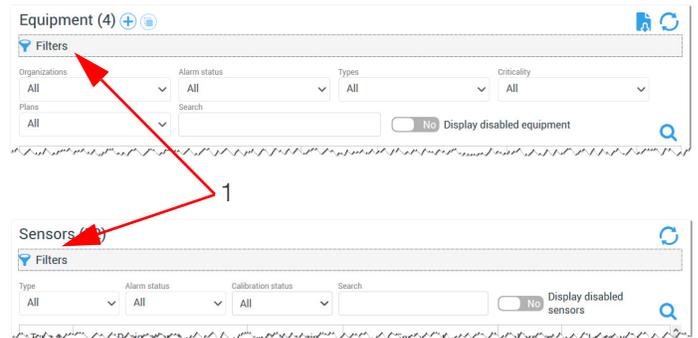


Figure 17. Filters

2. You may filter by entering text in the **Search** field.
3. Apply the filters by clicking the search icon (🔍). The list only displays items related to the criteria you specified.
4. To clear the active filter(s), simply uncheck the selected options and click on the search icon (🔍).

Home Screen Overview

The Smart-Vue Pro home screen contains the equipment monitoring dashboard which provides quick access to information about the system. The home screen features two display modes:

Plan View: Displays an image of the facility's floor plan to place the equipment you are monitoring according to its physical location (see **Placing Equipment on the Plan**).

Watch mode: Allows you to visually check the status of all the equipment and information such as the last recorded reading, sensor status, alarm status and more.

You may choose your default view using the **"Favorites"** bookmark, as shown below.

To pin your preferred view to your home screen, click on the desired bookmark (1) as shown:

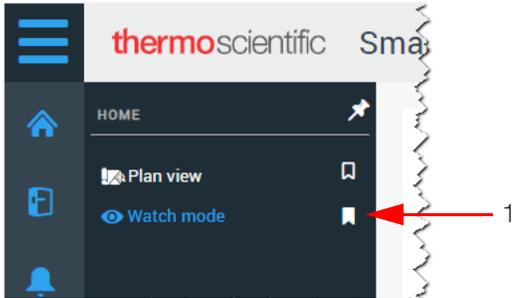


Figure 18. Customizing Smart-Vue Pro home screen

The selected view is opened automatically the next time you login to Smart-Vue Pro.

Plan View

Smart-Vue Pro enables you to load a floor plan for a quick visual indication of where your equipment and devices are located. The following is an example showing equipment at various locations on the floor plan (1). For more details, see **Adding Plan Images**.

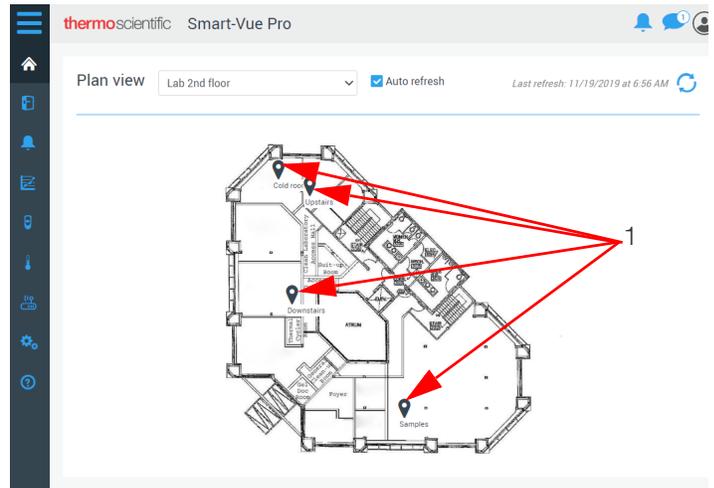


Figure 19. Viewing your sensors on a floor plan of

Watch Mode

The **Watch Mode** screen shows the status of all the equipments your profile allows you to view. The thumbnails on the dashboard provide an instant view of your equipment as well as the physical parameters recorded by your data loggers. Enable **Auto refresh** (1) to update the data on your dashboard automatically every five minutes.

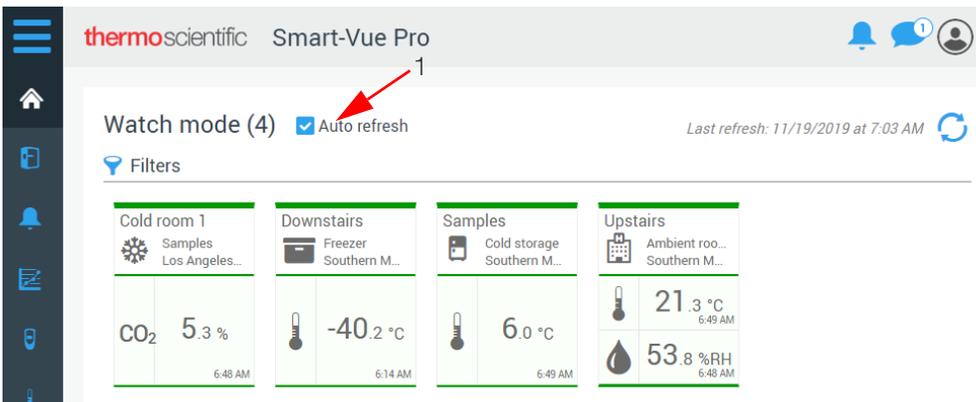


Figure 20. Overview of the Watch Mode



CAUTION: Actions and updates with respect to remote data loggers may not always be visible instantaneously. For example, data transfer is based on the configured transfer interval. You may observe a reading on a Smart-View Pro data logger screen that has not yet been transferred to the system.

The color of the thumbnail reflects sensor status:

- **Green:** Equipment with sensors enabled and functioning within programmed high and low alarm limits (if configured).
- **Red:** All alarms are indicated in red. This includes low limit alarms, high limit alarms, faulty sensors, and technical alarms (low battery or wireless transmission issues).
- Sensors that are not currently logging data are shown with a hyphen in place of the sensor value.

Each piece of equipment is represented as a tile showing the key details at a glance:

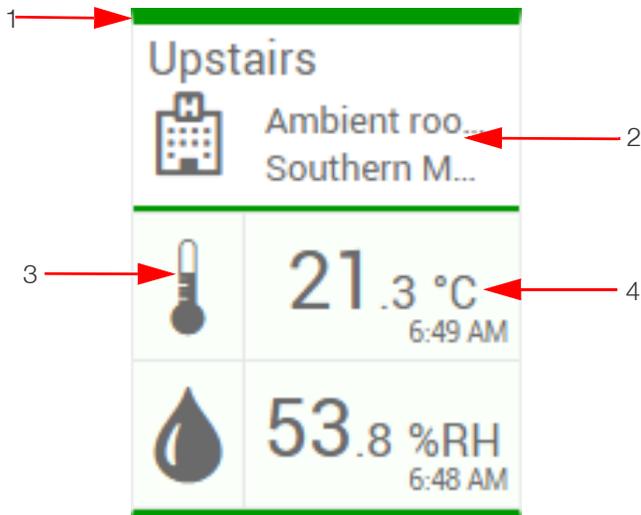


Figure 21. Equipment tile with key details

1. The color coding reflects the status of the overall equipment. Green indicates that there are no alarms; red indicates an alarm state.
2. Equipment name, type and associated organizational structure.
3. Measurement icon displayed in color according to the sensor status.
4. Latest reading with the color indicating sensor status.
 - The time of the latest reading is indicated if within the past 24 hours. Otherwise, click on the tile for more details.

- A “-” is shown in place of a value if the sensor could not be read.

The same color-coding is also used in the detailed pop-up pane described below. Click on a tile to open equipment details or sensor details:

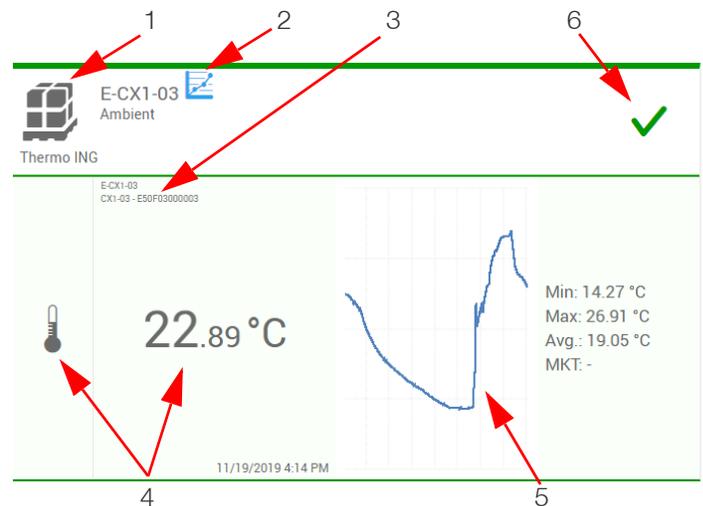


Figure 22. Summary display of selected equipment

Table 4. Equipment Details

S.No	Description
1	Equipment name, type and associated organizational structure.
2	Click on the graph icon to view readings recorded by the sensor(s) associated with this equipment (measurements and events) or to download the report in PDF, CSV or Excel format (see Viewing Sensor Data).
3	The first line (top) indicates data logging name and second line (bottom) indicates data logger name.
4	Visual indicator of the connected sensor and latest reading displayed in color according to status (with or without an alarm). The date and time of the latest reading are displayed as well as the date and time of the alarm (if present).
5	Displays the detailed sensor graph or download reading reports in PDF, Excel or Word format (see Viewing Sensor Data).

Table 4. (Cont.)Equipment Details

S.No	Description
6	<p>If there are no alarms, a green ✓ is displayed. Otherwise, the number of alarms is shown in red (low limit, high limit, technical alarm):</p>  <p>The number increments with the number of detected alarms.</p> <p>Click on the alarm counter to get more information, acknowledge one or more alarms, or generate an alarm report. This leads you directly to the alarm management page (see Alarms & Alerts).</p>

Click on the detailed equipment pane to close the equipment window.

If the system detects an alarm while you are checking your dashboard, an alarm panel appears on the right-hand side of the screen (1), with counters to indicate the number of pieces of equipment with alarms. This feature alerts if there are any critical incidents. Click on the alarm panel to open the alarm management window (see **Acknowledging Alarms**):

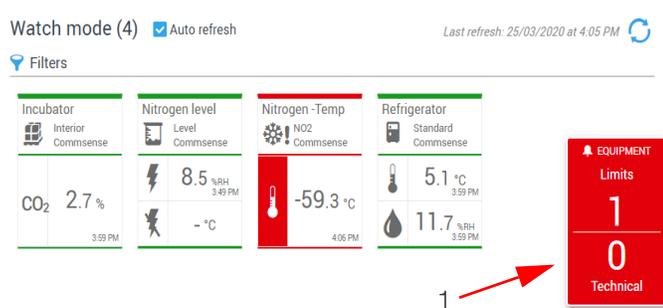


Figure 23. Pop-up for alarm notification

Users

Smart-Vue Pro includes a complete user management interface. When you create your company account in Smart-Vue Pro (described in **Creating your Smart-Vue Pro Company Account**), you also create the first user. The user defined has the role of an **“Application Manager”** who can create user profiles as well as manage all aspects of the system.

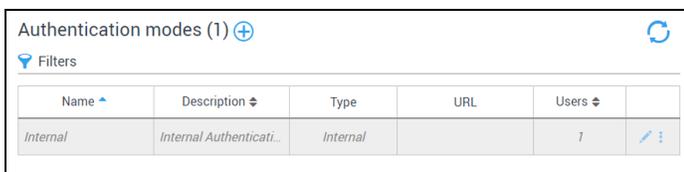
Authentication Modes

Smart-Vue Pro supports several user authentication modes:

- Smart-Vue Pro’s own **integrated user authentication platform** identifies users by their user-name and password.
- An **external authorization platform, namely LDAP** (Lightweight Directory Access Protocol) identifies users by their corporate, network or system login name and password. Generally used by large organizations, this option provides consistent and non-redundant user management while offering a broader range of user permissions and password control within corporate IT tools.

Default User Authentication with Integrated Platform

In a default system, only the Smart-Vue Pro internal authentication platform is used. To see the authentication mode(s) used by your system, click **Configuration**  **Authentication modes:**



Name	Description	Type	URL	Users	
Internal	Internal Authenticali...	Internal		1	

Figure 24. Default Smart-Vue Pro authentication platform

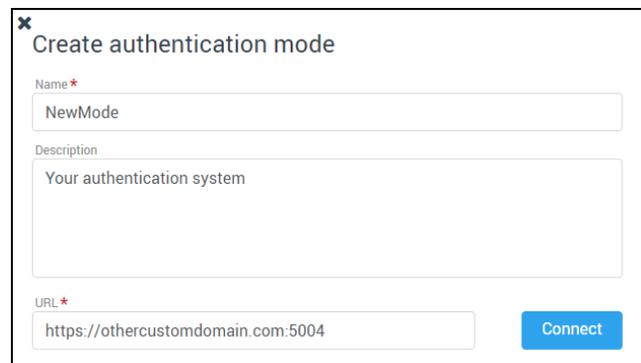


CAUTION: The internal authentication mode may not be disabled or edited in any way.

Adding an External Authentication Platform: LDAP

For your system to access an external LDAP authentication platform, you must add a new authentication mode to Smart-Vue Pro and configure it as described:

1. In the main menu, click **Configuration**  **Authentication modes:**
2. Then click  **(Add mode)**, which opens this screen:



Create authentication mode

Name *
NewMode

Description
Your authentication system

URL *
https://othercustomdomain.com:5004

Connect

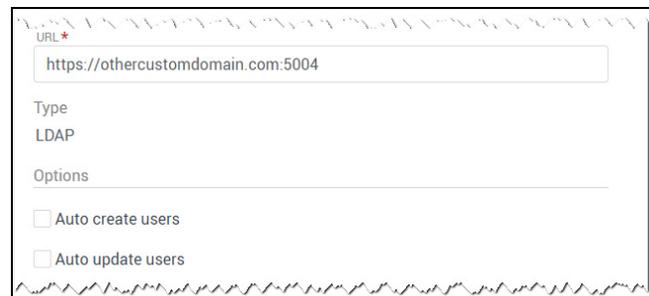
Figure 25. Adding an external authentication platform

Name: Assign a name for the authentication mode. This name is used as a reference by Smart-Vue Pro only and must be unique within your entire Smart-Vue Pro system. If the name is already in use, you will receive an error message.

Description: Enter optional information in this field.

URL: Enter the exact URL for connecting to your platform’s LDAP authentication agent.

3. Click **Connect**. If the URL is correct and connection to the LDAP authentication agent was established, additional information is displayed:



URL *
https://othercustomdomain.com:5004

Type
LDAP

Options

Auto create users

Auto update users

Figure 26. Automatic options with LDAP

In this example, the authentication agent was identified as LDAP, which has the following two options:

Auto-create users: When users who do not exist in the system try to login for the first time, the system creates them automatically. They can login to Smart-Vue Pro and continue with their regular credentials. Their user profile then shows which authentication mode is being used (i.e. the name assigned above in step (2)).

Auto-update users: With this option, information from user profiles is automatically updated based on information in the external directory.



CAUTION: The user's e-mail address must be:

1. Entered into the LDAP directory, otherwise the user will not be imported.
2. Unique in the system (the same address cannot be used more than once)

Editing an External Authentication Platform: LDAP

You may edit an authentication mode that you have added to the system (but not Smart-Vue Pro's native authentication):

1. Click on **Configuration**  → **Authentication modes**.
2. To change information, click on the Edit button () on the line containing the authentication mode you want to edit.

Edit authentication mode

Name *
LDAP

Description

URL *
http://localhost:5004

Type Hash
LDAP TGZGVm5YTU5HWkNIVXd2aU12cmtnYS9qMkV...

Options

Auto create users

Auto update users

Cancel Save

Figure 27. Editing an external authentication platform

If you change the URL to the authentication agent, the system checks to make sure that the new URL corresponds to the same agent as before. If not, an error occurs and you cannot save changes.

Options for External Authentication Platforms

Several options are available when you have configured one or more external authentication platforms.



CAUTION: You cannot edit or delete the system's internal authentication mechanism.

To access options for external authentication platforms (LDAP):

1. Click on **Configuration**  → **Authentication modes**.

Authentication modes (2) 

Filters

Name	Description	Type	URL	Users	
Internal	Internal Authenticati...	Internal		4	
LDAP		LDAP	http://localhost:5004	0	

Figure 28. Editing an external authentication platform

2. Click on the vertical dots () (1) to access options:

Synchronize: This option connects to your LDAP agent and fetches all the users configured as Users or Application Managers for Smart-Vue Pro and creates them automatically. This allows you to include users in alert rule definitions even if they have not yet connected to Smart-Vue Pro (which would create those users upon their first login if you enable the Auto-create option described previously). A message is displayed to confirm the operation, informs you if any required information is missing from any LDAP profile, or if the connection could not be established.

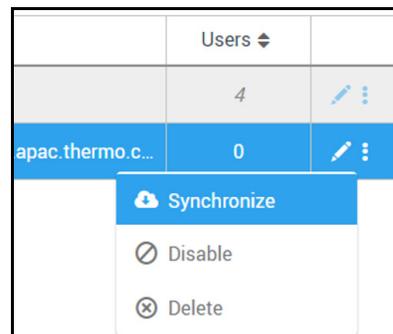


Figure 29. Synchronize

Note: If, **Auto-create user** and **Auto-update users** are not selected (see **Figure 26**), the synchronize button only updates the status of already-existing users (Enabled, Disabled, Expired).

Disable: Prevents associated users from connecting to the system.

Delete: You may only delete an existing authentication mode if there are no associated users.

Adding a New User

Login to Smart-Vue Pro as an Application Manager and follow the steps to add a new user:

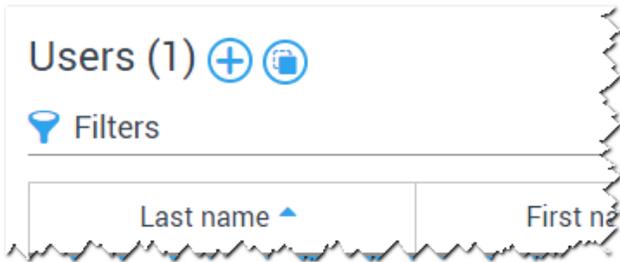


Figure 30. Add new user

- In the main menu, click on **Configuration** → **Users**.
You may either: Click on (Add user) or click on a user-name in the list and then (Create from selected) to use an existing user as a template for creating a new one.
- The new user identification window opens on the right side of the screen, with fields either empty or with information:

Figure 31. User details

- Enter the required information in the fields marked with a red asterisk (*):

First name / Last name: Enter the first and last names.

E-mail: Enter the e-mail address to receive e-mail alerts and confirmations sent by the system. Email id must be unique as per 21CFR.

User-name: If you are using Smart-Vue Pro’s integrated authentication, the user-name can either be the same as the e-mail address or it may be a name that you enter here.



CAUTION: When logging in, you must use the information in the **User-name** field.

If you are using an LDAP authentication agent, the user-name is your regular corporate system user-name (in the format: authentication_mode /username).

Authentication mode: This field enables you to choose whether to use Smart-Vue Pro’s internal authentication (based on user-name and a password) or an external LDAP platform using the regular corporate system credentials. The option specified when creating the company account is displayed here by default.

When you select LDAP, a confirmation button is displayed to test access directly from this screen.

Job title: Enter the role within the company.

International phone: Enter the telephone number in international format, with the prefix “[country code]”, such as: +14153817894. Do not include any extra digits or leading zeros.

Expiration date: When using Smart-Vue Pro’s native authentication mode, you may set a date at which this user account will expire. For external authentication platforms, you may not change the expiration date here.

Role: Assign the overall role within the Smart-Vue Pro solution by choosing one of the following options:

- Application manager:** Controls all aspects of the entire system, from managing licenses to adapting the individual profiles of other users.
- User:** This type of user starts with view only access.

Additional Rights such as acknowledging alarms and managing data logging are then added by an Application Manager at an organizational level (See **Organization with Sites and Departments**), offering flexibility.



CAUTION: When using external authentication platforms, you may not change the role directly via Smart-Vue Pro.

Authorize PIN code: This option allows you to determine whether the user is allowed to perform actions that require a PIN code:

- Acknowledging alarms via the Smart-Vue Pro data logger screen.
- Acknowledging telephone alerts.
- Accessing advanced settings menus on the Smart-Vue Pro data logger screen.

Yes Indicates that you can have a PIN code.

No Indicates that you will not have a PIN code.

PIN code: If you are authorized to have a PIN code, enter 4 digits into this field. For additional security, the system will complete the PIN code with 2 additional random digits.



CAUTION: To see the full 6-digit code, open your profile and select **PIN code** → **Show**. No other system users, including Application Managers, can see another user's full PIN code.

Language: You may change the application language at any time. Select your default language from the **Language** pull-down menu.

Time zone: Select the time zone based on the geographical area when first logged in.

Date format: Choose the date format to be displayed in the application.

Time format: Choose the time format to be displayed in the application. Readings can be displayed in 12-hour (AM/PM) or 24-hour format.

Unit: Choose the unit of temperature to be displayed in degrees Celsius (°C) or Fahrenheit (°F).

Decimal separator: Select the character (period or comma) to use as a decimal separator in numerical values in the Smart-Vue Pro display.

4. Click **Save** to create the user profile or **Cancel** to return to the previous screen.

Editing a User

To edit any information in your Smart-Vue Pro system:

1. Click **Configuration**  → **Users**.
2. Click on a line in the table to view or edit the profile.

3. The user identification window is displayed:

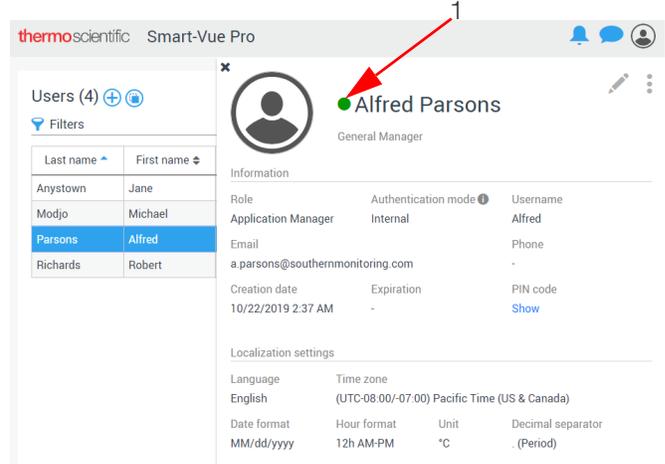


Figure 32. Details for selected user

4. The account status is indicated next to the name (1): indicates green if the status is OK; red if the account is expired or blocked; or grey if the account is deactivated.
5. To edit any information, click on the button  in the top right-hand corner of the window and edit the fields as needed.
6. Click **Save** to save your changes or **Cancel** to exit this screen without saving changes.



CAUTION: Changes made to an existing user:

1. The user's e-mail address cannot be changed, as it serves as a unique identifier for this user in the system.
2. If an external authentic mode (LDAP) is being used by the system, and the option "Auto-update users" is activated, then some user information presented here may not be modified. If you change the Authentication mode, test the user's account validity by clicking on the test button: .
3. For traceability reasons, it is not possible to delete a user profile from the database. A profile can only be disabled.
4. Any changes will take effect the next time the user logs in to the system.

Updating Contact Details



CAUTION: If you want the user to receive alert notifications by e-mail and/or SMS, you must first assign the user to a Call Group as described in **Alert Rules with Users**.

To edit contact details for a user associated with an alert rule:

1. Click **Configuration**  → **Users**.
2. Click on a line in the table to edit that user's profile.
3. The user details appear at the right of the screen.
4. Click **Options**  at the top right of the user profile and then click **Edit contact information**.

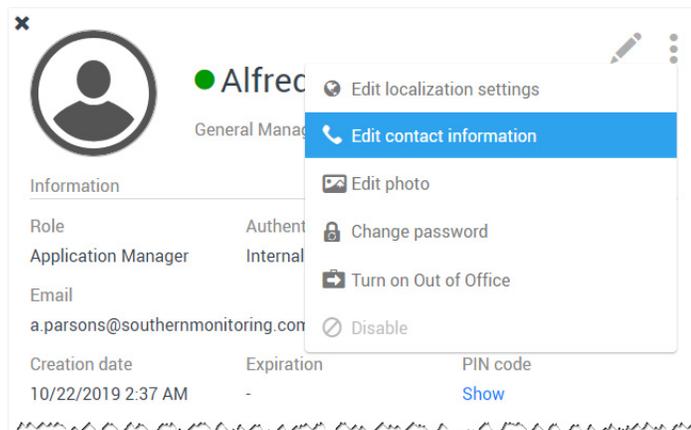


Figure 33. Accessing contact information

Enter the telephone numbers and/or e-mail addresses for the different types of alerts you have configured, as described in **Configuring Alert Rules**.

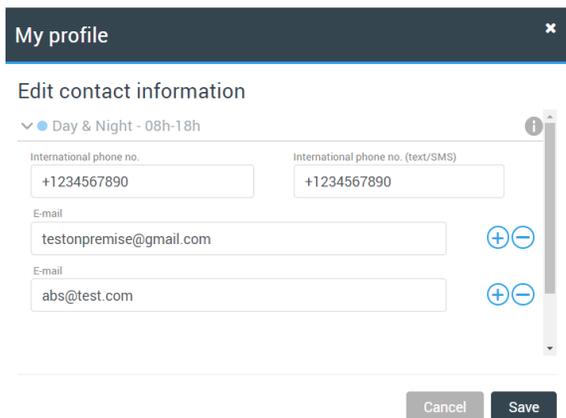


Figure 34. Editing contact information



CAUTION: These fields are independent from any external authentication (LDAP) platform information.

With Smart-View Pro, you may customize alert rules to meet your needs (such as daytime hours, night calls, shift work, etc.) and each alert rule can contact multiple users. Once you add a user to an alert rule, you may modify that user's contact information accordingly, such as to have different phone numbers or e-mail addresses at different times. For the various alert types, you may add up to 2 phone numbers and 3 email addresses by clicking on Add .



CAUTION: Smart-View Pro Alert notification via voice messaging works with all types of phones in many regions. SMS/text messages are supported on most mobile phones. Remember to enter the "+" sign and country code prefix for both your mobile and land-line phone numbers. You cannot disable a user account if that user is associated with an alert rule.

Changing a User Password on a System with Integrated Authentication



CAUTION: This section does not apply to users in systems using external authentication (LDAP). The audit trail is updated to reflect a password change, unless the new password and existing password are identical.

Who can change a password? Users can change their own passwords. Users with Application Manager rights can initiate a password reset for other people but cannot actually set the password for someone else. In all cases, users must set up their own passwords.

To change your own password:

1. Click on your user profile in the upper right-hand corner of the screen then click on **Profile**.
2. Your profile window opens directly.
3. Click **Options**  in the top right-hand corner of the window and select **Change password**.
4. Enter the new password and re-type it in the second field to confirm.
5. Click **OK** to save your changes or **Cancel** to return to the previous screen.



CAUTION: Passwords should be at least 8 characters long and contain a combination of upper-case and lower-case letters with at least one number and one special character.

Resetting a User Password

An application manager may launch the password reinitialization process for another user. That user will be forced to change their password upon the next login:

1. Click **Configuration** (⚙️) → **Users**.
2. Click on the user profile for which you want to reset the password. The user details appear at the right of the screen.
3. Click **Options** (⋮) at the top right of the form and select **Reset password**.
4. In the password reset window, click **Yes** to proceed with the change or **No** to return to the previous screen:

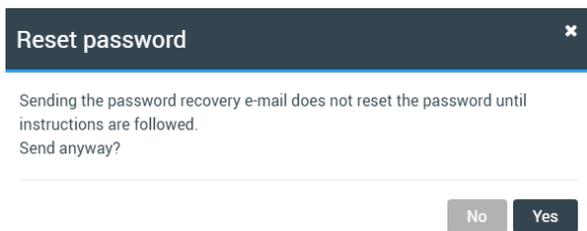


Figure 35. Password reset confirmation

5. The password recovery email is sent to the user with instructions to reset the user password as shown below:

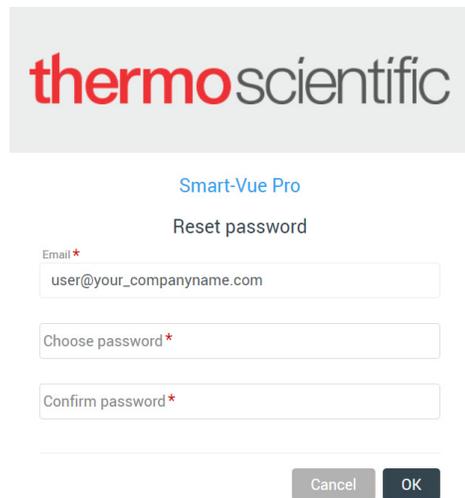


Figure 36. Resetting a user's password

6. The change is taken into account when the user clicks **OK**.
7. The regular login screen is displayed the next time the user logs in and user is prompted to enter the new password. The user cannot login without changing the password.

Note: If the user's account is not locked or expired, it is still possible to connect even if the password has not been reinitialized.

Adding an Account Profile Picture

Smart-Vue Pro can personalize your display with a profile picture when you are logged in to the system. The picture is shown in the upper right-hand corner of the screen.

Picture recommendations:

- Maximum size: < 1 MB
- JPG or PNG format

To add a profile picture:

1. Click on your user profile image (👤) in the upper right-hand corner of the screen, then click **Profile** to open your profile directly (or click **Users** in the left-hand menu and select your name in the user list).
2. Click **Options** (⋮) in the top right-hand corner of the window and select **Edit photo**.
3. Click **Choose file**:

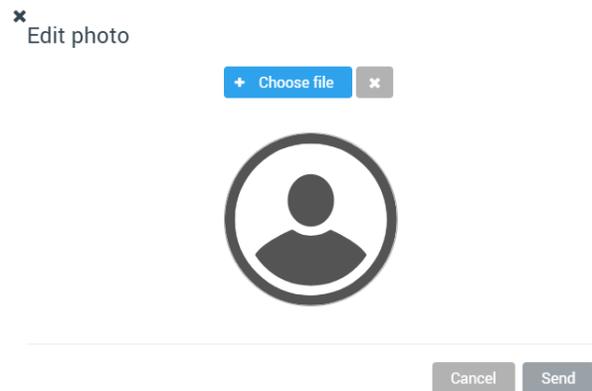


Figure 37. Selecting a profile picture

4. Browse your computer to locate the image you want to use and then click **Open**.

5. The image is uploaded and displayed in Smart-Vue Pro.

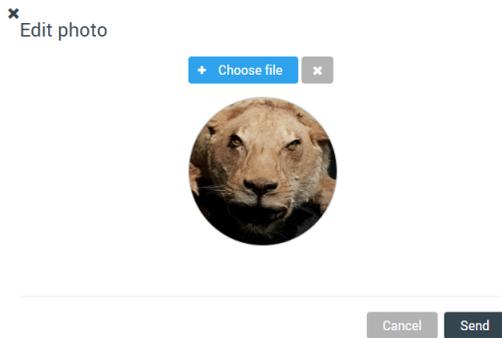


Figure 38. Deleting a profile picture

6. Click **Send** to save your changes and update your profile.
7. To delete the picture linked to the user profile, click () in the screen shown above.

Account Profile Overview

To view/edit your general profile information:

1. Click on your user profile image () in the upper right-hand corner of the screen, then click **Profile**. The screen contains two main sections:

Information: This section indicates the status of your profile (Active / Inactive), its creation date and expiry date (if entered), your user level, the language selected and the PIN code to be typed on a Smart-Vue Pro data logger screen to acknowledge alarms. You may edit this information by clicking on the **Edit** icon ().

Localization settings: Display format for time zone, time format, date format, measurement unit, and decimal separator. You may edit this information by clicking **Options** () → **Edit localization settings**.

2. Click **Save** to record your changes, or **Cancel** to return to the previous screen without saving changes.

User Access Profiles (Roles)

Initially created with a “User” profile, Smart-Vue Pro users can have a variety of different access rights to match tasks, responsibilities and the organization to which they belong.

The concept of profiles described in this section is closely tied to that of your “organization”.



CAUTION: Users with a user profile must be given rights within the organization in order to view and manage configured equipment or data logging information.

An Application Manager profile can manage all aspects of the system, whereas a user has limited access to application features and menus. User access is further refined by assignment to the organizational as a whole or one or more sites and departments.

Table 5. Access differences between Application Managers and Users

Main menu items	Application Manager profile	User profile	
		Access	Visibility based on Organization-Site Department
Home (Dashboards)	All	Limited	Yes
Equipment	All	Limited	Yes
Alarms	All	Limited	Yes
Reports	All	Yes	Yes
Data loggers	All	No	No
Sensors	All	Yes	No
Infrastructure	All	No	No
Configuration	Company Users Profiles Organizations Time slots Days off Alerts Data logging templates Plans Authentication modes	Organizations	Yes
Help	Yes	Yes	No

Adding a New Profile

Profiles are comprised of different access rights or roles. A user may have different roles in different parts of an organization. For example, an user could be limited to viewing information in one department and only be allowed to acknowledge alarms in another.

A profile is assigned to a user based on their roles and responsibilities in your organization. Profiles enable you to set specific permissions and authorizations for each type of user.

Smart-Vue Pro starts with three default profiles: **Manager, User and Acknowledger**. Only Application managers can add or customize user profiles.

To add a new user profile:

1. In the main menu, click on **Configuration**  **Profiles**.
2. Then click on  **(Add profile)**:

Profiles (3)  

Name ↕	Description ↕	
Manager	Authorized to manage equipment and data logging	 
Acknowledger	Authorized to acknowledge alarms	 
User	View only	 

Figure 39. Available user profiles

3. The **Create profile** window opens as shown here:

x

Create profile

Name *

Description

Rights

Manage organization

Manage data logging

Acknowledge alarms

Manage metrology parameters

Figure 40. Adding a new user profile

Fill in the fields as required:

Name: Enter a name for the new user profile.

Description: Enter a description for the user profile (optional).

Rights: The rights listed here define what a user is and is not allowed to do with Smart-Vue Pro.

To grant permissions, check the right(s) you want to assign to the profile:

Manage organization: The user can set up and configure the system's organizational structure and include other users as allowed.

Manage data logging: The user can set up data logging for equipment.

Acknowledge alarms: The user can acknowledge alarms.

Manage metrology parameters: The user can update calibration settings.

To assign all these permissions at once, click on the check box above the rights list, which selects all check boxes.

Your Organization

Company Information

You may update all company information at any time, except for the main contact e-mail address.

General Information

To update your company account:

1. Click **Configuration**  → **Company**.

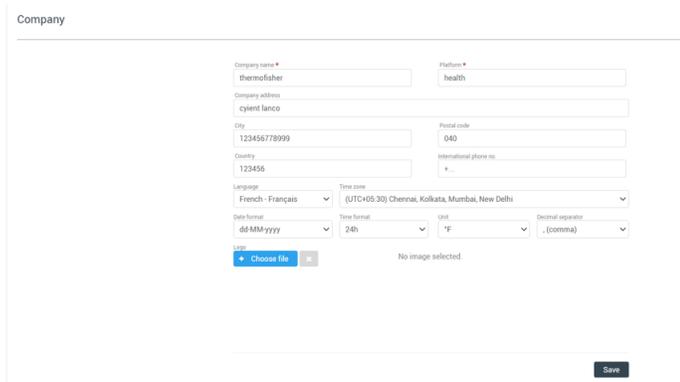


Figure 41. Company account information

2. Fill in the required fields (marked with *) as appropriate.

Company name: Enter the name of the company.

Platform: Unique company identifier, filled in automatically based on the company name entered.

Company address: Enter City, Postal code, Country, International phone number relevant to the company.



CAUTION: Make sure you enter the telephone number in international format, with the prefix “[country code]”, such as: +14153817894.

Do not include any extra digits or leading zeros.

The following settings apply as a default for your system. They can be overridden by each user’s individual settings for local operation and display purposes, but this information is used as a common reference for the company in case users have conflicting profile settings for shared actions. For example, if you have e-mail alerts sent to user in different regions, English could be the default company language for a same alert message.

Language: Choose the desired default language for your company. Users are free to choose their personal preference at any time.

Time zone: Select the time zone based on the primary geographical area for the company.

Date format: Select how dates are displayed in the application.

Time format: Select how time is displayed in the application. Readings can be displayed in the **12-hour (AM/PM)** or **24-hour** format.

Unit: Temperature can be displayed in degrees **Celsius (°C)** or **Fahrenheit (°F)**.

Decimal separator: Select the character to use as a decimal separator in numerical values.

Logo: If you want to add a logo image for your company, click on **Choose file** (JPG or PNG format, maximum size 1 MB). This image will be used in a future release to personalize application features

3. Make the desired changes directly in the form and then click on **Save** to save the information.

Managing License Keys

Access to Smart-Vue Pro is based on one or more service subscriptions linked with the number of sensors in your system. The following services require a subscription:

- **Smart-Vue Pro web application** and all integrated features (including 24/7 alerts by e-mail only) for a specified number of measurement points.
- **Smart-Vue Pro Alert notification system**, an additional service that provides 24/7 alert notification by telephone voice call and/or SMS-text message (to a cellular telephone).



CAUTION: Contact an authorized Thermo Scientific representative for more information.

When you sign up for one (or both) of these subscriptions, you will receive the associated voucher number(s) to enter in Smart-Vue Pro.

To view information about your Smart-Vue Pro solution or add a new license key:

1. Click **Configuration** (⚙️) → **License keys**.

This page shows information related to your licenses, such as activation key codes and expiration dates. The Thermo Scientific technical support team may ask you to provide this information if you contact them for assistance.

2. To enter a new voucher number, click (+) **(Add license key)**.

License keys (1) (+)

License key	Product	Description	Expiration
SJAAA-AAASB-AEAM-AQQLS-GICAA-YLVA4	Smart-Vue Pro	3 measuring points	25/04/2020

Figure 42. Adding a new license key

3. Enter your new voucher number in the pop-up window:

Add licence key
✕

Cancel
Save

Figure 43. Entering the new license key number

4. Click **Save** to register your license or **Cancel** to discard changes and return to the previous screen.
5. Once the license key is registered, the activation key number is shown in the table along with the license type and expiration date.

License keys (2) (+)

License key	Product	Description	Expiration
SJAAA-AAASJ-A2GAQ-CPQBD-AQAAA-A3AP4	Smart-Vue Pro	20 measuring points	2021-04-14
CJAAA-AAACA-AGKAA-AQYRZ-MGDAA-YOKDU	SmartVueAlert		2020-09-27

Figure 44. Monitoring system’s licenses

License Expiration (90-day Reminder)

As the expiration date for any of your licenses draws closer, a reminder is displayed in the Notifications menu (🔔) as shown here:

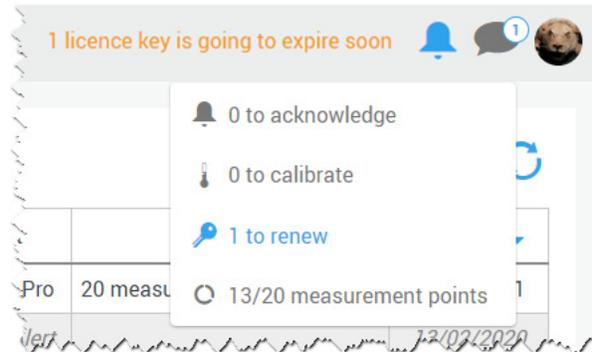


Figure 45. Reminder to renew license

This reminder is displayed starting 90 days before the license expiration date and remains present until the issue is resolved.

Lines containing license keys that have reached their expiration date are shown in Grey with italic text:

Licence keys (2) (+)

License key	Product	Description	Expiration
SJAAA-AAASJ-A2GAQ-CNIFD	Smart-Vue Pro	20 measuring points	10/21/2020
<i>CJAAA-AAACA-AKAAA-ANIIS</i>	<i>SmartVueAlert</i>		<i>10/29/2019</i>

Figure 46. Expired licenses indicated in Grey

Renewing or Upgrading your License

To renew the license within the expiration date or upgrade your current license (such as to expand the number of authorized measurement points), enter the new license number as described earlier in **Managing License Keys**.

Organization with Sites and Departments

The Smart-Vue Pro monitoring system is comprised of one **Organization**, which may be sub-divided into **Sites** and **Departments**. This flexibility enables you to create a logical hierarchy for managing your system based on geographical, operational or any other criteria that may be useful to you.

For example, you could create a hierarchy as shown in the following figure

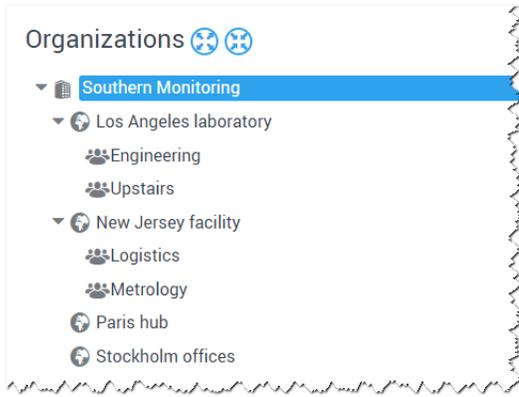


Figure 47. Creating a hierarchy that is appropriate for your organization

Users, equipment and infrastructure devices are assigned to the various entities in the overall organization. Users may even have different profiles in different parts of the organization such as someone who is a manager in the Los Angeles laboratory but has view-only rights for the facility in New Jersey.



CAUTION: Equipment and infrastructure devices (such as receivers) can only be assigned to one entity in the hierarchy.

An organization can have from one to three levels (Organization, Site and Department). There is no requirement to create sub-divisions with this feature, though it tends to make management much more intuitive than a “flat” structure. You may associate users at whichever level of the organization works best for you. Here are three examples:



Figure 47a. Sample organizational structures with users assigned at level 1

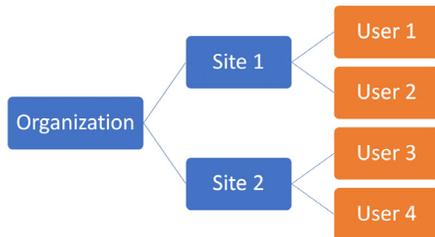


Figure 47b. Sample organizational structures with users assigned at level 2

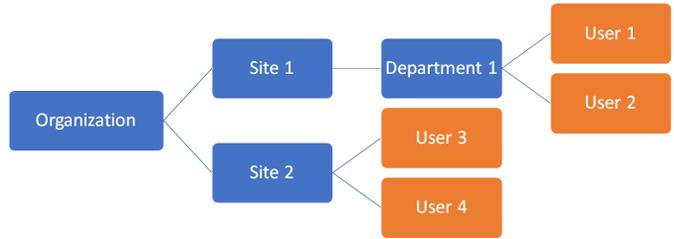


Figure 47c. Sample organizational structures with users assigned at level 3

Assigning an User to a Department

When you first start, you have one organization with the Application Manager – the user who created the account – listed as Manager (profiles are discussed in the previous chapter). Application managers are associated automatically with all the nodes in the organizational structure. Those users cannot be removed.

You must first create users (see **Users**) in order to add them to your sites and departments.

To associate a user with a department:

1. In the **Configuration** → **Organizations** menu, click **Associate users (+)** (1).

Name	Email	Profile	
Anystown Jane	j.anystown@southernmonitor...	Acknowledger	
Parsons Alfred	a.parsons@southernmonitor...	Manager	

Figure 48. Adding a user to a department

The **Associate user** window opens as shown here:

Associate users

Selected users will be associated with **Los Angeles laboratory** with selected profile.

Profile *

Available	Selected
<input type="text" value="Modjo Michael (com@wanadoo.com)"/>	<input type="text" value="Richards Robert (r.richards@southernmonitor...)"/>

Figure 49. Associate user window

2. Depending on the department to which the user is assigned, select the profile you want to assign to them

using the **Profile** drop-down list. A profile determines the actions that the user can perform and the objects that they can manage in Smart-Vue Pro.

There are three user profiles:

- **Acknowledger:** This user profile allows you to acknowledge alarms issued by the system for the sensors they are monitoring. You may also disable those sensors, if necessary.
- **Manager:** A Manager can access all application features and manage other users.
- **User:** User rights are limited to viewing only. These users do not have access to management and configuration windows.

3. Add users to your organization:

Available: Select a user from the list of Available users. Then click **>** to move the user to the list of Selected users on the right-hand side.

To select multiple users at a time, press and hold the Ctrl key and click on each of the users you want to select. Click **>** to add them to the list of Selected users. To select all available users, click **>>**.

Selected: To remove an user from your organization, click on that user in the list of Selected users and then click **<** to move the user to the list of Available users on the left-hand side.

To remove multiple users, press and hold the Ctrl key and, while holding down the key, click on each of the other users you want to remove. Click **<** to move them to the list of Available users on the left. To remove all users, click **<<**.

4. Click **Save** to save your changes or **Cancel** to exit this screen without saving changes.

Adding a Site

To add a new site:

1. Click **Configuration (⚙️) → Organizations**.
2. Click on the name of your company account and then click **Add site** (1).

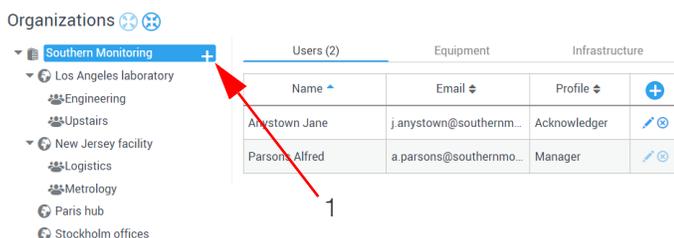


Figure 50. Adding a new site

3. Enter a name for the new site in the pop-up window and click **Save**.

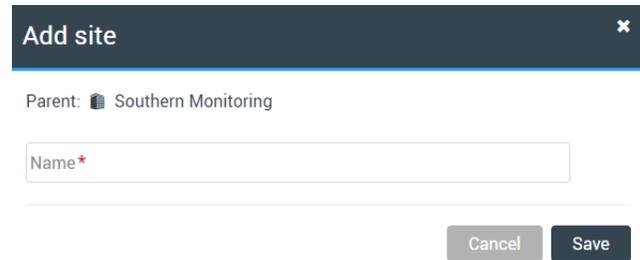


Figure 51. Entering the new site name

The new site appears in your organization's tree structure. You may click **⊕** **⊖** or the triangles (**▼**) in the tree to expand and contract the view.

Adding a Department

1. Click **Configuration (⚙️) → Organizations**.
2. In the organization's tree structure, click on the site and then on Add department (2):

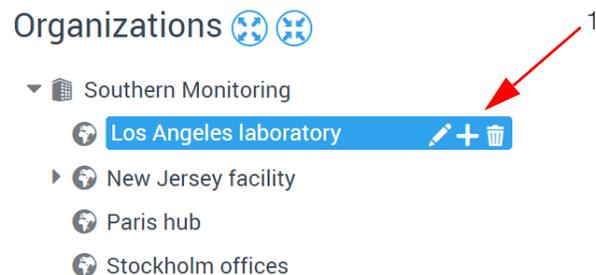


Figure 52. Adding a new department

3. Enter a name in the pop-up window for the new department and click **Save**:

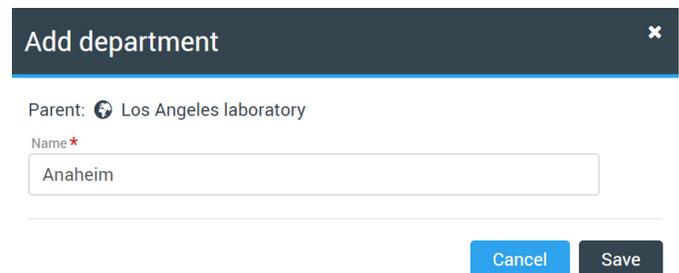


Figure 53. Entering the new department name

The new site is listed in your organization's tree structure:

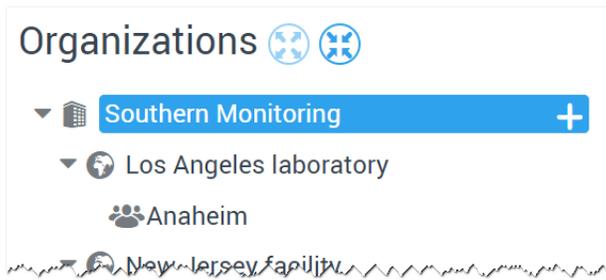


Figure 54. Organizational structure

Editing / Deleting a Site or Department

To manage different sites and departments, click on the desired node in the organizational structure and use the available options:

New To add a new site or department, click .

Delete To delete a site or department, click .

Edit To edit a site name, click . Modify or enter a new name and click  to confirm your changes.

Adding Plan Images

Smart-View Pro allows you to visually place your equipment on floor plan representing the layout of your organization. Plans are useful for identifying the locations of the sensors associated with the equipment you are monitoring. You may add as many plans as required.

To add a plan:

1. In the main menu, click **Configuration**  → **Plans**.
2. Click **Add plan** () to import a plan image.

Plans (1) 

Name	Equipment	Description	
Lab 2nd floor	0		

Figure 55. Managing plans

The create plan window opens as shown here:

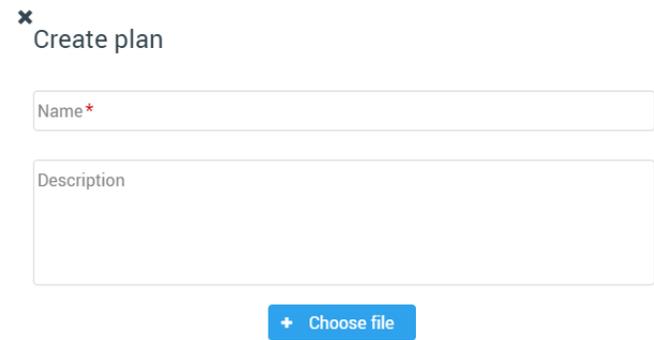


Figure 56. Adding a new floor plan

3. Enter a name for the floor plan in the **Name** field.
4. You may add a description in the **Description** field (optional).
5. Click **Choose file** and select the desired image of your premises (JPG, PNG or GIF format up to 1 MB maximum size). A preview image is displayed after you load the floor plan:

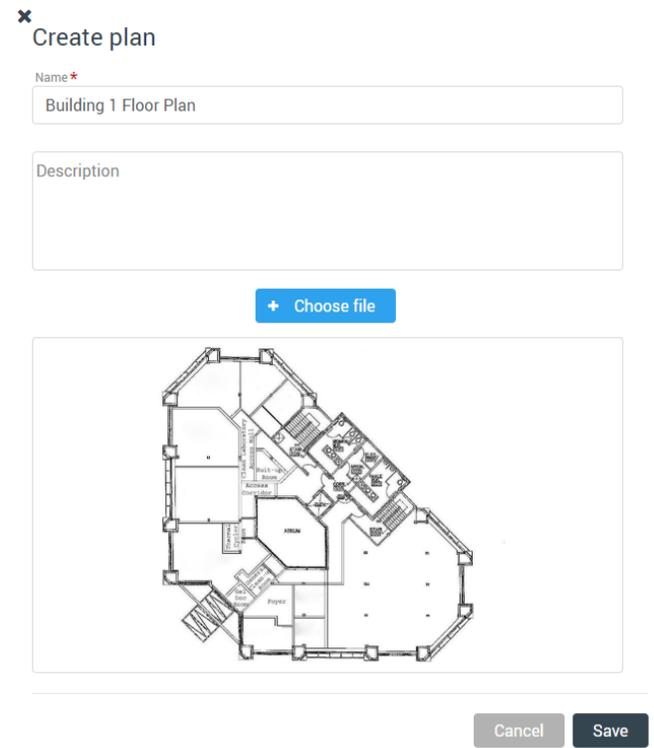


Figure 57. Adding a new floor plan

6. You may click on the thumbnail image to open the enlarged floor plan.

Placing Equipment on the Plan

Once you have loaded one or more floor plan images, place the equipment on the image according to its physical location.

To place equipment on the plan:

1. In the main menu, click **Equipment** (🏠).
2. Select the equipment you want to place on the plan. The equipment details are shown on the right-hand side of the screen.
3. Click **Options** (⋮) in the upper right-hand corner of the equipment window and then click **Edit location**.
4. Select the desired floor plan using the **Plan** drop-down list to display a larger image.
5. Place your equipment by clicking the desired location on the image. The blue location pin (📍) moves to show the new position (1). The grey pin represents the previous equipment position, which disappears after you save the new equipment location.

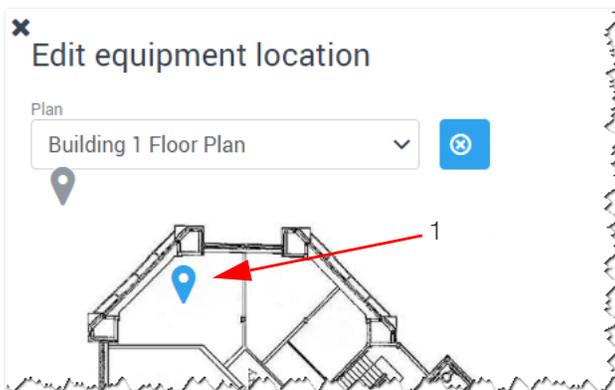


Figure 58. Placing an equipment on the plan

If you want to remove the equipment from the image, you may click on the Remove association button (⊗).

6. Click **Save** to save the equipment location (or removal) or click **Cancel** to return to the previous screen without saving the changes.

Infrastructure Devices

This section allows you to add different infrastructure components that comprise your monitoring system, such as a LoRaWAN receiver, an alarm device, etc.,.

Your LoRaWAN connection must be available to interact with LoRaWAN enabled devices, that is, a Thermo Scientific LoRaWAN-enabled receiver must be installed and configured to communicate with your Smart-Vue Pro system.

Adding Infrastructure

Follow these steps to add a new component to the system manually:

1. Click **Infrastructure** in the main menu and then click :

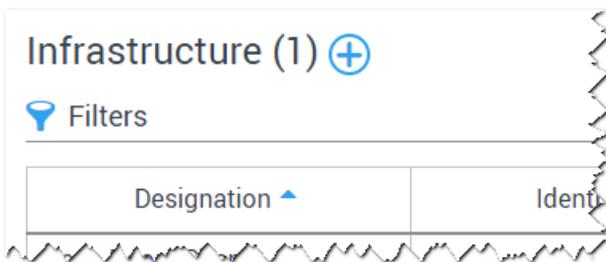


Figure 59. Click to add infrastructure

The **Create infrastructure** window opens as shown here:

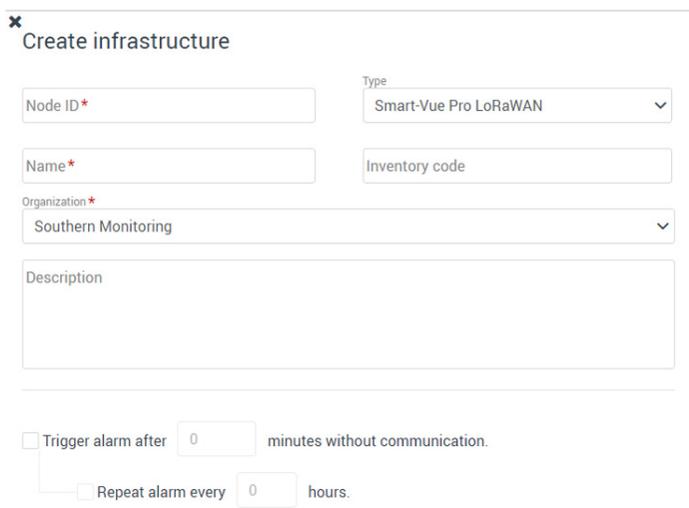
The image shows a "Create infrastructure" form. It has a close button (X) in the top left. The form contains several fields: "Node ID*" (required), "Name*" (required), "Organization*" (required, with a dropdown menu showing "Southern Monitoring"), "Type" (dropdown menu showing "Smart-Vue Pro LoRaWAN"), "Inventory code", and "Description" (text area). At the bottom, there are two checkboxes for alarm settings: "Trigger alarm after" (with a numeric input field set to 0 and the text "minutes without communication.") and "Repeat alarm every" (with a numeric input field set to 0 and the text "hours.").

Figure 60. Adding infrastructure details

2. Enter the details of the infrastructure element. Fields marked with a red asterisk (*) are required:

Serial number/Node ID: Enter the serial number on the back of your infrastructure element (device).

The system identifies the element based on its serial number. If you enter an incorrect serial number, a red exclamation mark is shown in front of the corresponding field. Check the serial number and try again. If the problem persists, contact technical support.

Type: For infrastructure elements, the Type is selected manually. It is therefore essential to enter the correct identifier (e.g. serial number, node ID). If the type does not correspond to the entered identifier, the application cannot communicate with the device.

Name: Enter a name for the infrastructure component. This name is used as the reference in all lists.

Inventory code: (Optional) Enter the inventory code of the infrastructure element for information purposes.

Organization: Use the drop-down list to select the department to which the infrastructure element is associated. It may be helpful to associate topology with an infrastructure element to target precise alert management and alarm acknowledgment rights.

Note: To add other departments, follow the procedure described in **Organization with Sites and Departments**.

Description: (Optional) You may enter a description of the component for information purposes.

Trigger alarm after: Check this option and enter the delay (in minutes) after which the system will trigger an alarm if the infrastructure component is not responding. This mechanism ensures early notification in case of problems or a communication error.

Repeat alarm every: Check this option to set the repeat frequency for this alarm and specify the desired time interval in the entry field.

3. Click **Save** to save your infrastructure component, or **Cancel** to close this window without saving changes.

Editing Infrastructure

To manage existing infrastructure, use the following options:

1. In the main menu, click **Infrastructure**.
2. Click on a line in the table to view or edit the desired device.

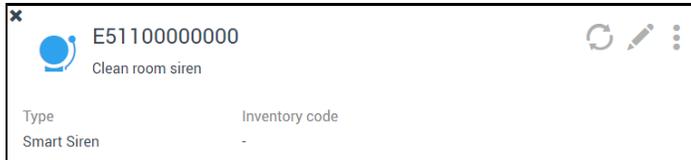


Figure 61. Editing infrastructure details

3. Click **Edit** (✎) on the right-hand side of the screen and edit information directly as necessary.
4. Several other functions are available under **Options** (⋮):

Options	Category	Description
Options (⋮)	Alarm history	Click to disable the siren's functionality (in the Infrastructure list, a filter enables you to display or hide disabled infrastructure).
	Simulate an alarm	You may use this function to simulate an alarm as described in the previous section
	Disable	Opens the Alarm list to see any alarms involving this device (a filter is applied to show "Infrastructure" with the siren's serial number)
	Remove	Click to remove the siren from your system. The siren is marked as "removed" in the Infrastructure list. If you wish to restore the same siren, you must create it again by entering the serial number.
	Refresh	Refreshes the infrastructure information window.

5. Click **Save** to keep your changes or **Cancel** to return to the previous screen without saving changes.

Adding a LoRaWAN Receiver as Infrastructure

LoRaWAN wireless technology is a key feature of the Smart-Vue Pro system.

In terms of configuration, the LoRaWAN receiver is entirely configured on the receiver itself, as described in the Smart-Vue Pro LoRaWAN Receiver User Guide, which you may access online via the **Help** menu of the Smart-Vue Pro web application. That is, the receiver contains its own configuration interface, which includes information on sending data to your Smart-Vue Pro system.

You are not required to configure the receiver via the Smart-Vue Pro web application. However, if you do not add the receiver to your Smart-Vue Pro infrastructure via the web application, then you will not receive alerts in case a technical problem arises.



CAUTION: Even if it is not technically required to add your LoRaWAN receiver as infrastructure via the Smart-Vue Pro web application, we recommend that you include it along with your other infrastructure devices in order to receive technical alerts in case of any problems.

Data Loggers

The features described in this section are reserved for application managers.

Adding Data Loggers

Adding a Data Logger Manually

Follow these steps to add a Smart-Vue Pro data logger to the system manually:

1. In the main menu, click Data Loggers (📁) and click (+) to enter all the details for a new data logger, or (📄) to use the selected data logger as a template for the new one (1):

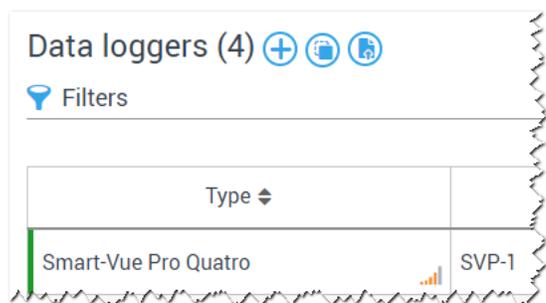


Figure 62. Data logger management window

The **Create data logger** window opens as shown here:

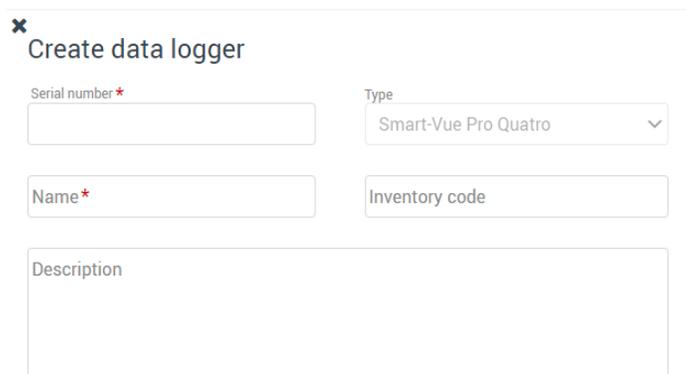


Figure 63. Data logger management window before entering the serial number

2. As you type in the data logger serial number, additional fields are displayed:

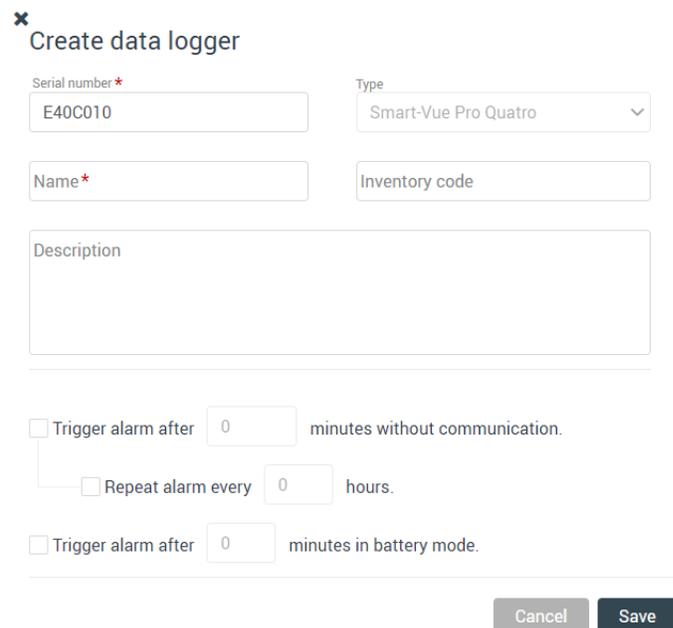


Figure 64. Adding a new data logger

3. Fill the data logger details as required. Options are displayed according to the type of data logger. The fields marked with a red asterisk (*) are required.

Serial number: Enter the serial number from the front of data logger / from the information screen of data logger. If you enter an incorrect serial number, a red exclamation mark is shown in front of the corresponding field. Check the serial number and try again. Contact technical support if the problem persists.

Type: This information is assigned automatically by the system based on the serial number you enter.

Name: Enter a name for the data logger. This name is used as the reference in all lists.

Inventory code: (Optional) You may enter an inventory code for the data logger for information purposes.

Description: (Optional) You may enter a description of the data logger for information purposes.



CAUTION: Alarms are always sent in real-time, without waiting for the programmed transfer interval.



CAUTION: It is a general practice to allow 3 communication failures before triggering an alarm.

4. Click **Save** to save the data logger in the database.

At first, the data logger is shown as follows:

Figure 65. Freshly added data logger

5. On your Smart-Vue Pro data logger: Tap the screen to open the menu and click **Sensors** → **Refresh** sensors to detect physically connected sensors and ensure that the data logger is up-to-date internally before continuing.

6. In the Smart-Vue Pro web application: Click **Refresh sensors on device**  icon (1) (above) to communicate with the data logger and update the screen with the latest information.

Designation	Ranges	Latest value	
0334120000001041 PT100	50.00 °C -180.00 °C	-	
DF34120000001441_I 4-20mA/0-5V/0-10V	20.00 %RH 4.00 %RH		
DF34120000001441_U 4-20mA/0-5V/0-10V	10.00 °C 0.00 °C		
E40C0104576F_T Sensor	85.00 °C -40.00 °C		

Figure 66. Example of data logger information after refresh

Adding a Data Logger with the Same Settings as an Existing One

If you are setting up a data logger with settings that are already configured, you can save time by clicking the  icon to use the selected device as a template for the new one.

Importing a Batch of Data Logger

If you would like to set up several data loggers at once, you may import a TXT or .CSV text file with values separated by a semi-colon character (;). You must include at least the device serial number and name, as shown in the following figure:

```

SerialNumber;Name;
E40C0104###1;Incubator 1
E40C0104###2;Incubator 2
E40C0104###3;Incubator 3
E40C0104###4;Incubator 4
E40C0104###5;Incubator 5
E40C0104###6;Incubator 6
E40C0104###7;Incubator 7
E40C0104###8;Incubator 8
E40C0104###9;Incubator 9
E40C0104###10;Incubator 10
E40C0104###11;Incubator 11

```

Figure 67. Sample .CSV import file with data logger serial numbers and names

You may also include the inventory code and data logger description as shown here:

```

SerialNumber;Name;InventoryCode;Description;
E40C0104576D;Incubator 1;123451;Lab data logger 1;
E40C01045731;Incubator 2;123452;Lab data logger 2;
E40C0104576F;Incubator 3;123453;Lab data logger 3;
E40C0104577B;Incubator 4;123454;Lab data logger 4;
E40C0104577A;Incubator 5;123455;Lab data logger 5;

```

Figure 68. Sample .CSV file with additional information

1. Click  to import a batch file.
2. Browse to locate the.TXT or .CSV file on your computer.
3. Click **Open** to import the list.

Viewing the Data Logger List

All the data loggers you add to the system are shown in the **Data loggers** screen:

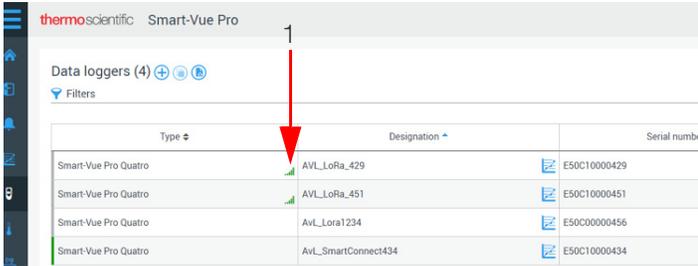


Figure 69. List of data loggers stored in Smart-Vue Pro

This screen displays all data loggers in the system, with columns containing the information below. The data logger's status is indicated by red-green color coding on left-hand edge of the table (1) (this does not indicate alarms related to readings, just the data logger itself). You may click on the up/down arrows (2) in the column headers to change the list for ascending/descending order.

Type: Type of data logger

Designation: A name you assign to the data logger. You may click on the graph icon (3) to see a data graph and other sensor information.

Serial number: Data logger serial number.

Version: Version of Firmware running on data logger.

Sensors: Number of sensors connected to the data logger.

Last activity: Date and time of last communication with the data logger.

Power status: Plugged into AC adapter (4) or battery level (5).



CAUTION: The number of sensors that you may use in your system is determined by your license. You can only use the allowed number of sensors simultaneously.

Viewing Data Logger Details

Regardless of your Smart-Vue Pro user level, you may always check the data for the data loggers you are authorized to view.

To view the graph for a given data logger:

1. Click **Data loggers** in the main menu.
2. Select a data logger.

The data logger details are shown on the right-hand side of the screen, with several sections:

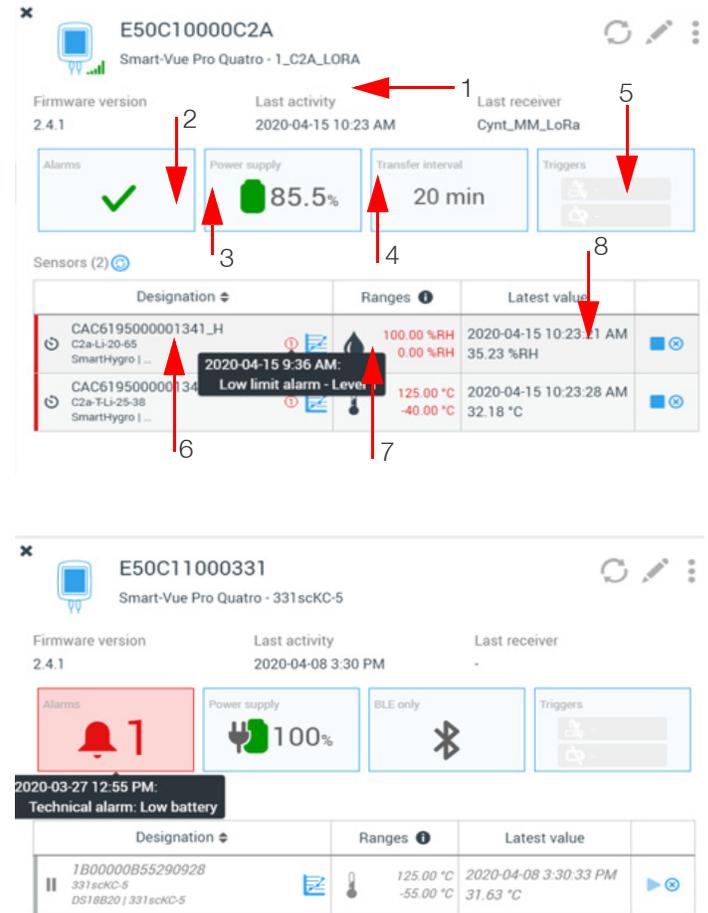


Figure 70. Data logger details

Table 6. Data Logger Details

SNo.	Description
1	The serial number, name and inventory code (if present) of the data logger
2	<p>Alarm counter indicating data logger status (sensor related alarm and unexpected stop indication are not shown in data logger alarm status). In case of alarm, the icon changes color and shows the number of alarms, for example:</p>  <p>Figure 71. Alarm counter in data logger details</p> <p>Hover over the alarm counter with your mouse to view alarm details. If there are no alarms, the indicator  is shown.</p> <p>Click on the alarm counter to get more information, acknowledge one or more alarms, or generate an alarm report. You will be directed to the alarm management screen (see Alarms & Alerts).</p>
3	Indicator showing the power source being used and estimated remaining battery level.
4	The Transfer interval indicates the frequency with which data is transmitted by the data logger to the Smart-Vue Pro system.
5	Alarm triggering and repeat rules, as defined when adding the data logger. Hover over the alarm counter with your mouse to view alarm details.
6	<p>Number of sensors associated with the data logger. The active icon  indicates sensors with data logging in progress.</p> <p>Sensors not yet programmed are greyed out, as indicated by the inactive icon .</p>
7	Click on the graph icon to view readings of sensor over time (measurements and associated events.) or download the report in PDF, Excel or Word format (see Viewing Sensor Data).
8	Sensor operating range.

Data Logger Actions

You may use the options available in the data logger details screen to manage your data loggers.

1. Click **Data loggers**  in the main menu and select a data logger in the list.
2. Click **Options**  (1) in the upper right-hand corner to access available options:

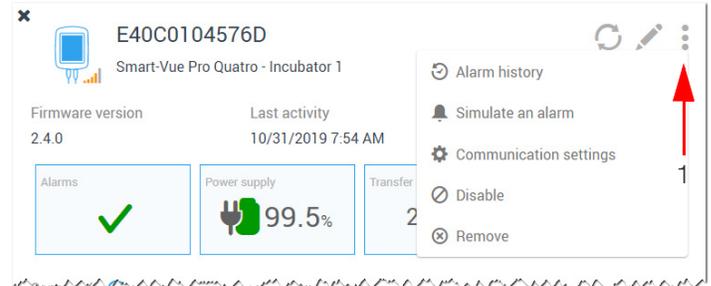


Figure 72. Data logger options menu

Alarm history: Opens the alarm page for the selected data logger (see **Viewing Alarm Details**).

Simulate an alarm: Allows you to simulate the occurrence of an alarm on this data logger for testing purposes. The simulation generates a Technical alarm (Communication lost, Device mismatch, Lost power supply, Low Battery) that is listed in the Audit Trail as an “Alarm simulation”.

Communication settings: This menu provides technical information regarding data logger communications. Reserved for qualified technicians. Incorrect use of data logger settings could cause your system to malfunction. Contact Thermo Scientific technical support for more information.

Disable: Enables you to effectively turn this device off and stop data logging.

Remove: Enables you to completely remove this data logger from your system and stops data logging. You may need to do this in the event of a data logger swap or hardware problem. If you choose this option, you must enter your password to confirm the operation.

Editing Data Logger Details

Follow these steps if you need to modify the information or default values for your data logger:

1. Click **Data loggers** () in the main menu and select a data logger in the list.
2. Click **Edit** () icon. This opens a window identical to that described in **Adding Data Loggers**.

Note: You may not edit the serial number or data logger type fields.

3. The following information is non-editable:
 - Firmware version: the data logger's firmware version number.
 - Last activity: the last time the data logger communicated with the server.
4. Click **Save** to keep your changes or, **Cancel** to return to the previous screen without saving changes.

Sensors

Sensors that are connected physically or wirelessly to your Smart-Vue Pro data loggers are added to the Smart-Vue Pro web application automatically.

For Smart-Vue Pro to recognize the sensor(s) that are physically connected to a Smart-Vue Pro data logger, click **Refresh Sensor** and **Synchronize** in the data logger's **Sensors** menu. Once detected, these sensors are shown in the sensor management window. From the sensor management window, you may:

- Access detailed information about the sensor.
- Display readings in a graph, along with various other information about the selected sensor.
- Manage calibration parameters.

Viewing the Sensor List

Sensors connected to the Smart-Vue Pro data loggers are listed in the Sensors screen (🌡️):

Type	Designation	Data logging	Equipment	Calibration	Latest value
⚡	DF34120000001441_J 4-20mA/0-5V/0-10V Lab 1st...	🔄 Control room	Control room	🕒	11/21/2019 8:12 AM 15.91 %RH
⚡	DF34120000001441_U 4-20mA/0-5V/0-10V Lab 1st...	🔄 Control room	Control room	🕒	11/21/2019 8:12 AM 5.31 °C
🌡️	0334120000001041 PT100 Lab 1st Floor	🔄 Samples	Downstairs Freezer	🕒	11/21/2019 8:14 AM -20.15 °C
🌡️	E40C0104576F_T Internal Sensor Lab 1st Floor	⏸	-	🕒	-
🌡️	A655443322110028 DS18B20 Room 1	🔄 Samples	Samples Cold storage	🕒	11/21/2019 8:11 AM 5.00 °C
🌡️	E40C0104576D_T Internal Sensor Room 1	⏸	-	🕒	-
CO ₂	B635120000001541 CO2 Room 1	🔄 Incubator	Cold room 1 Samples	🕒	11/21/2019 8:11 AM 5.31 %

Figure 73. Sensor list in Smart-Vue Pro

This screen displays all sensors with the following information:

Type: Visual indication of sensor type.

Description: Sensor serial number, sensor type and associated data logger.

Data logging: Data logging status: running (🔄) or idle (⏸). Move your mouse over the icon to see data logging settings.

Equipment: The equipment associated with the sensor.

Calibration: Hover your mouse over the field to see calibration parameters or click to open the calibration management window.

Latest value: Date and value of the most recently recorded reading.

Viewing Sensor Details

Regardless of your user level in Smart-Vue Pro, you may always check the data for sensors you are authorized to view.

To view detailed information about a sensor:

1. In the main menu, click **Sensors**.
2. Select the desired sensor.
3. The sensor details are shown on the right-hand side of the screen, with several sections:

Equipment Samples **Device Room 1**

Limits alarms **Technical alarms** **Data logging** **Latest value**

Data logging

Name
Samples

Start mode
Immediate

Start
11/21/2019 7:35 AM

Reading interval
1 min

Latest value date
11/21/2019 8:52 AM

End
-

Limits
1: 2.00 °C | 8.00 °C

Figure 74. Sensor details

Laboratory: Using the drop-down list, select the calibration laboratory that issued the calibration certificate.

Identifier: Enter the certificate reference number.

Date: Click in the Date entry field and choose the date on which the sensor was calibrated from the calendar. The current date is displayed by default.

Note: If calibration certificate information had been entered previously for this sensor, the date you enter here must not be earlier than that of the previous certificate.

Calibration and Reminder Frequency

Once you have added calibration information, you may choose to receive notification as a reminder to recalibrate the sensor by a given date.

In the **calibration** screen, click **Edit** (✎) and enter the number of months after the certificate date at which the system will remind you to recalibrate the sensor:

x Calibration parameters ✎ +

Sensor		
Serial number	Unit	
033412000001041	°C	
Parameters		
Coefficient A	Coefficient B	Coefficient C
1.0099	0.009	0.000100
Drift	Uncertainty	
0.00	0.00	
Certificate		
Laboratory	Identifier	Date
2	BABA	02/01/2018
Reminder notification		
No reminder notification. ✎ 🗑		
Show calibration notification <input type="text" value="6"/> months after certificate date. ✓ ✕		

Figure 76. Sensor calibration parameters

Click **Save** (✎) to update sensor information, **Cancel** (✕) to undo any changes you may have made, or **Delete** (🗑) to remove notification.

4-20 mA, 0-5 V, 0-10 V Loop Sensors

Description

4-20 mA current and 0-10 V voltage Smart-Sensors can be connected to industry-standard devices that generate appropriate output signal. The signal can be measured to provide status information for a variety of applications, including wind monitors, particle counters, ultra-low-temperature (ULT) freezers and temperature, gas, VOC, and relative humidity devices, and more. Read values are converted into useful information by the Smart-Vue Pro web application, then displayed both in the web application and on the Smart-Vue Pro data logger display.



Figure 77. 4-20 mA / 0-10 V Smart-Sensor

Table 8. Sensor wiring

Black	Common ground
Red	0-10 V input
White	4-20 mA input

Connect the wires according the type of device you want to monitor.



CAUTION: Do not connect all three wires

A sensor may only monitor one of the two input parameters (red or white wire, but not both).

The maximum voltage on the 4-20 mA input wire is 2 V. Exceeding this value will damage your sensor.

Mapping Values for Meaningful Readings

For sensor readings to be relevant in your context, you must map the low values (0 V or 4 mA) and high values (10 V or 20 mA) to match the limit values produced by your connected device. For example, on a 4-20 mA device, the low 4 value could correspond to a CO₂ level of 0% and the high 20 mA value could correspond to 10% CO₂. In that case, you would map 4 to 0 and 20 to 10 in Smart-Vue Pro. You may use other units to meet your needs.

To do this:

1. Click **Sensors** in the left-hand menu.
2. Select the sensor you wish to modify from the sensor list.
3. Click **Calibration** (CAL).
4. Edit the values in the **Range** zone (1) (below) to meet your needs.
Low: The lower value output by your device.
High: The higher value output by your device.
Unit: The measurement unit you would like to display in Smart-Vue Pro.
5. Click **Save** to save your changes or **Cancel** to close this window without saving changes.

✕ Edit calibration parameters

Sensor

Serial number: DF3412000001441 Unit: mA

Parameters

Coefficient A: 1.0499	Coefficient B: 0.050	Coefficient C: 0.000500
Drift: 0.00	Uncertainty: 0.00	

Range **1**

Low (4 mA): 0.00	High (20 mA): 100.00	Unit: %RH
------------------	----------------------	-----------

Certificate

Laboratory: Oceansoft Lab	Identifier: BBA	Date: 12/01/2019
---------------------------	-----------------	------------------

Figure 78. Mapping 4-20 mA / 0-10 V values

The unit you specify here is shown in the Smart-Vue Pro web application as well as on the data logger.



CAUTION: The mapping from 4 to 20 mA or 0 to 10 V is strictly linear.

The 0-10 V sensor is also used for 0-5 V output devices. Simply adjust your Low and High values accordingly.

Simulating a Sensor Alarm for Testing

You may trigger a limit alarm on a sensor for testing purposes to ensure that related alert rules and user contact work properly. To test a sensor alarm:

1. Login to Smart-Vue Pro as an Application manager.
2. In the main menu, click **Sensors** for a list of the sensors in your system.
3. Click on the sensor you want to test.
4. Click **Options** (⋮) and select **Simulate an alarm** from the drop-down menu. You will then be able to see the alarm in the Alarm list and Audit trail (indicated as "Simulation" in both places). Any related alert rules for the sensor are triggered, such as, e-mail or telephone alerts.

Removing a Sensor

You may remove sensors from the system, for example if they are no longer used or if they have been replaced on the Smart-Vue Pro data logger. To remove a sensor:

1. Login to Smart-Vue Pro as an Application manager.
2. In the main menu, click **Sensors** for a list of the sensors in your system.
3. Click on the sensor you want to remove.
4. Click **Options** (⋮) and select **Remove** from the drop-down menu. To remove a sensor, data logging must not be running.



CAUTION: Smart-Vue Pro data loggers update sensor information in Smart-Vue Pro automatically based on the physically connected sensor(s). When you plug in a new sensor (or even plug the same one back in if it had been removed accidentally), the sensor(s) will appear again in the Smart-Vue Pro web application.

Disabling / Enabling a Sensor

You may disable a sensor so that it cannot be used for data logging or you can enable a sensor that has been disabled. To do this:

1. Login to Smart-Vue Pro as an Application manager.
2. In the main menu, click **Sensors** from a list of the sensors in your system.
3. Click on the sensor you want to disable or enable.
4. Click **Options** (⋮) and select **Disable** (or **Enable**) from the drop-down menu. To disable a sensor, data logging must not be running.



CAUTION: Disabled sensors are no longer shown in the sensor list. To enable a disabled sensor, use the Filter option in the sensor list and select Display disabled sensors. You may then choose to enable it as described above.

Equipment

This chapter describes managing equipment in the Smart-Vue Pro system and assigning data loggers and sensors to that equipment.

The “equipment” in Smart-Vue Pro refers to the temperature-controlled chambers, cabinets, refrigerators, tanks, enclosures, and other spaces you may choose to monitor with Smart-Vue Pro data loggers.

Adding Equipment

To add equipment:

1. Click **Equipment** (📦) in the left-hand menu. There are two options for creating new equipment:
 - Click (+) to enter details manuallyOr
 - Click to select an existing piece of equipment and click (📦) to use that equipment as a template to create new equipment:

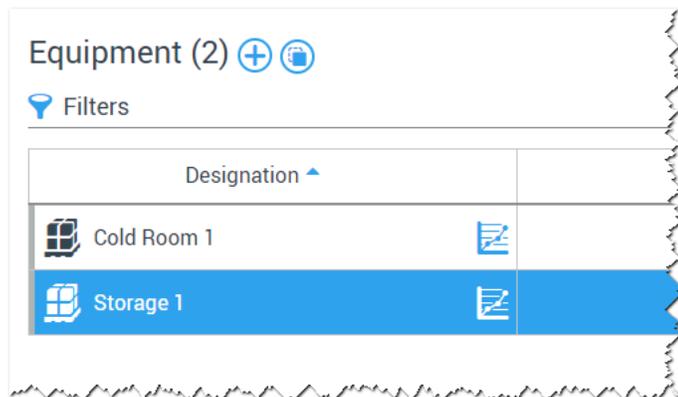


Figure 79. Equipment management window

The **Create equipment** window opens as shown here:

Create equipment

Name *

Inventory code

Icon Type

Organization * Southern Monitoring

Critical Not critical

Mobile Static

Figure 80. Adding new equipment

2. Fill the equipment details. The fields marked with a red asterisk (*) are required:

Name: Enter a name for your equipment. This name is used in all lists.

Inventory code: (Optional) You may enter an inventory code for the equipment for information purposes.

Type: Enter a type of equipment in the field.

Icon: Click on the icon and select the most appropriate symbol to visually represent the equipment.

Organization: Equipment is assigned to sites and departments.

You must assign an organization to your equipment using the pull-down menu. To add a department to this list, click **Configuration** (⚙️) **Organization** and follow the instructions provided earlier in **Organization with Sites and Departments**.

Critical / Not critical: Determine the criticality level for the equipment by clicking on the appropriate option.

This information is used to fine-tune the data logging mode and alert rules based on equipment criticality.

Mobile / Static: Indicate whether the equipment is dedicated to a static environment (refrigerators, freezers, water-baths) or a mobile environment (coolers, packages).

3. Click **Save** to save the equipment or **Cancel** to close this window without saving changes.

The equipment is listed in the equipment management screen:

Designation	Type	Sensors	Organization	Plan
Cold Room 1		0	Southern Monitoring	
Storage 1		0	Southern Monitoring	

Figure 81. Equipment list in Smart-Vue Pro

Associating Sensors with Equipment

This section describes assigning one or more sensors to monitor equipment.

To associate a sensor with equipment:

1. Click **Equipment** (🔧) on the left-hand side of the menu to see the list of equipments you are allowed to manage.
2. Click on the equipment to which you want to add a sensor.
3. Click on **Add** (+) (1):

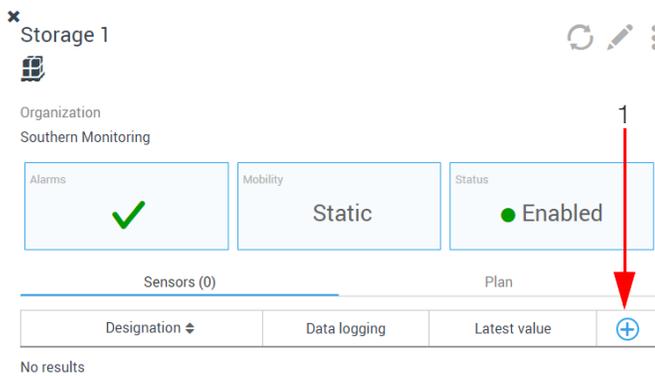


Figure 82. Equipment details

4. Tick one or more check boxes to select the corresponding sensor(s):

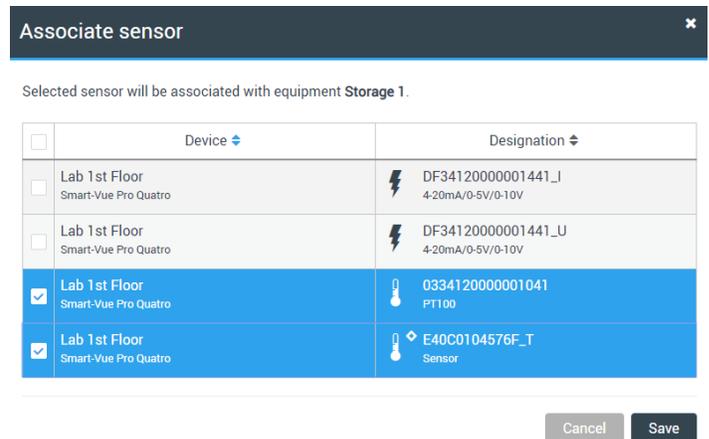


Figure 83. List of available sensors

5. Click **Save** to complete the association or **Cancel** to exit this screen without saving changes.

Dissociating Sensors from Equipment



CAUTION: You may only dissociate sensors from equipment if they are not being used for data logging. If the sensor is being used for data logging, you must first stop data logging as described in **Stopping Data Logging**, and then dissociate the sensor from the equipment.

To dissociate a sensor from equipment:

1. Click **Equipment** (🔧) in the main menu to see the list of equipment you are authorized to manage.
2. Click on the equipment from which you want to remove a sensor.

- Click **Remove association** (⊗) (1) next to the sensor you want to remove:

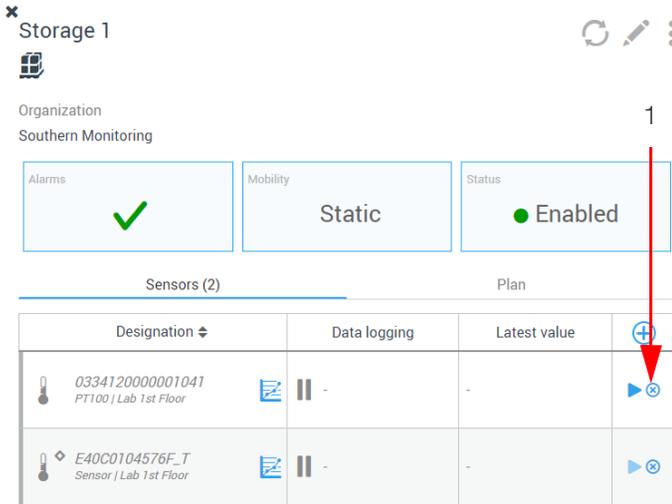


Figure 84. Dissociating a sensor from equipment

- Click **Yes** and confirm the operation when prompted, or click **No** to return to the previous screen:

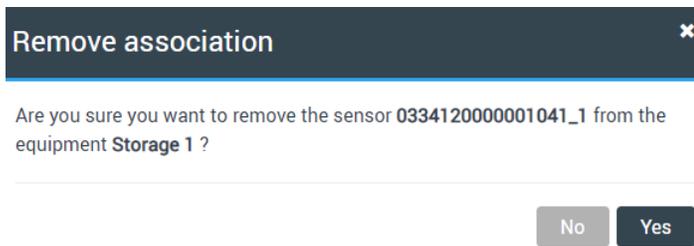


Figure 85. Confirming sensor dissociation

Viewing Equipment Details

Regardless of your user level in Smart-Vue Pro, you may always check the data for the equipment you are authorized to view.

For a complete view of all equipment details, including sensor readings:

- Click **Equipment** (📦) in the main menu.
- Select the desired equipment.

- The equipment details are displayed as shown below:

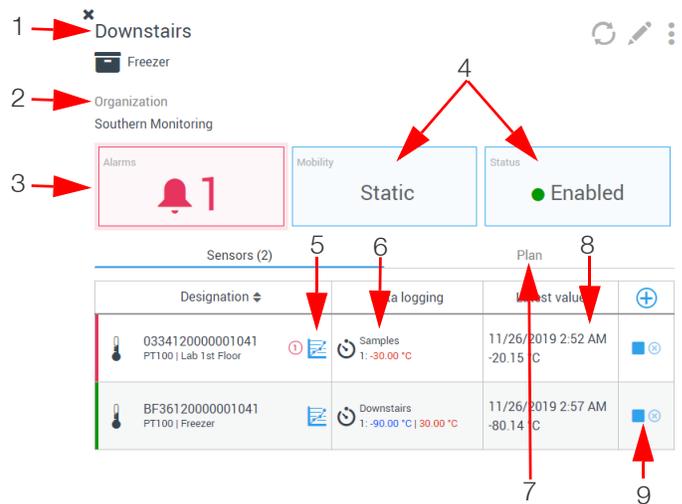


Figure 86. Equipment details

Table 9. Equipment Details

SNo.	Description
1	Equipment name and type.
2	Organization and department (if configured) to which the equipment is associated.
3	Current alarm indicator. An alarm is an “out-of-bounds” state or other error detected on a data logger or sensor. Hover your mouse over the indicator or click on it for details or to acknowledge an alarm.
4	Mobility indicator (Mobile or Static) and equipment status (Enabled or Disabled).
5	Opens a graph with all sensor readings.
6	Programmed upper and/or lower limits for the sensor and indication of data logging status.
7	If you have placed the equipment on a plan, click Plan to view the exact location of the equipment on the map.
8	Shows most recent values reading for this sensor. A dash (-) is displayed if no readings are available.
9	<ul style="list-style-type: none"> Click to associate other sensors with the selected equipment. Click to remove the sensor from its current equipment. Click to set sensor settings and start data logging (password required). Click to stop data logging (password required).

Modifying / Managing Equipment

You may use the options available in the equipment details screen to manage the equipment in your Smart-Vue Pro system.

1. Click on **Equipment** (📦) in the main menu.
2. Select the desired equipment.
3. Click on **Options** (⋮) in the upper right-hand corner of the screen and then click on one of the available options:

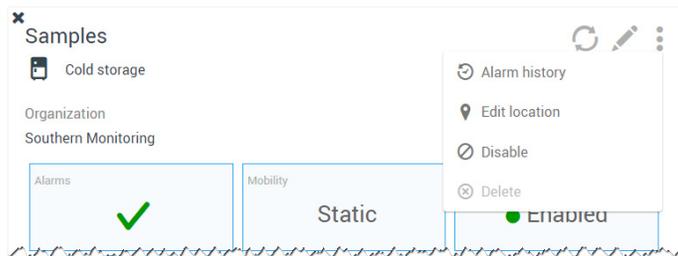


Figure 87. Equipment details

Alarm history: Opens the alarm page for the selected equipment (described in **Viewing Alarm Details**).

Edit location: This option is only available for "Static" equipment, allowing you to locate the equipment on a floor plan or change its position on the plan (see **Placing Equipment on the Plan**).

Note: Disable / Enable: Select this option to disable equipment or enable equipment that has been disabled. Disabled equipment is no longer shown in the equipment list or as a tile in the Watch Mode display.

Note: Equipment cannot be disabled if data logging is currently running on one of its sensors. To enable disabled equipment, use the Filter option in the equipment list and select **Display disabled equipment** to enable it.

Delete: Removes the selected equipment from the system.

Edit: To modify the information for existing equipment, click on **Edit** (✎). Click **Save** to keep your changes, or **Cancel** to return to the previous screen without saving changes.

Refresh: Click **Refresh** (🔄) to update the information in this window.

Deleting Equipment

You may delete equipment only if the following conditions are met:

- The equipment must not be associated with a sensor.
- If one or more sensors are associated with the equipment, you must first unpair the sensor(s) before deleting the equipment. See **Dissociating Sensors from Equipment** for more details on this feature.

To remove equipment:

1. Click **Equipment** (📦) in the main menu.
2. Select the desired equipment.
3. Click **Delete** (⊗) (1) to remove the selected equipment:

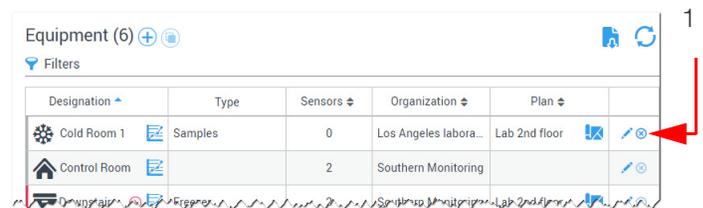


Figure 88. Equipment management window

4. Click **Yes** to confirm the operation when prompted, or click **No** to return to the previous screen:

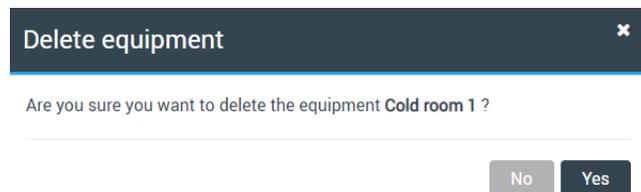


Figure 89. Confirming equipment removal

Data Logging

To use your Smart-Vue Pro data logger, you must first program data logging.



CAUTION: Data logging is configured at the equipment level. Once you have assigned a sensor to equipment (described in **Associating Sensors with Equipment**), you may set up data logging.

Data logging consists of recording physical parameters (such as temperature and humidity) over time. Sensor readings are recorded based on a set of configuration options that are used to manage your data logger while data logging.

Data logging parameters include:

- The sensor reading frequency.
- High and low alarm limits.

How does it work?

Here is a quick overview of the data logging flow with Smart-Vue Pro:

1. The Smart-Vue Pro data logger equipped with sensors is installed on the equipment to be monitored.
2. Sensor readings are collected by the Smart-Vue Pro data logger, which then sends the data to the Cloud or your server via LoRaWAN wireless connectivity.
3. Smart-Vue Pro enables you to manage every aspect of your system, equipment, users, and data. When limits are exceeded or other anomalies are detected, an alert is displayed on the Smart-Vue Pro data logger. You can receive alerts 24/7 from the Smart-Vue Pro system.

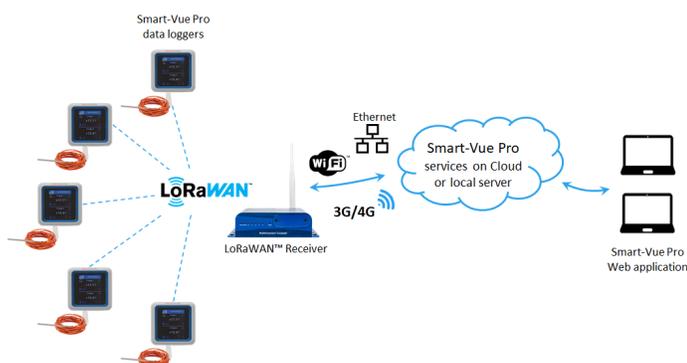


Figure 90. The Smart-Vue Pro data logger communicates with Smart-Vue Pro via an on-premises LoRaWAN receiver

Starting Data Logging



CAUTION: A sensor must be associated with equipment in order to start data logging.

To configure data logging:

1. You can start data logging either via **Equipment** (🔧) or **Data Loggers** (📄) in the left-hand menu as long as the sensor is associated with equipment. The instructions below are identical for both cases.
2. Click on the name of the equipment (1) (or data logger) for which you want to start data logging:

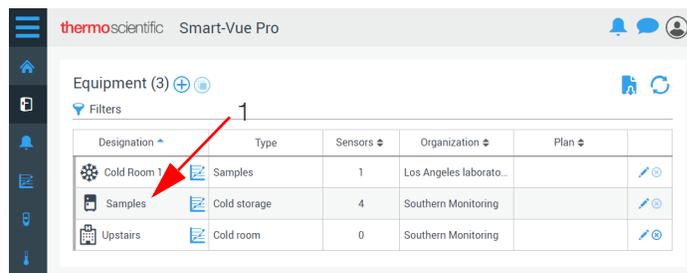


Figure 91. Click on a piece of equipment to set up data logging

3. Details and available sensors are shown on the right-hand side of the screen. Click **Start data logging** button (▶) (2) to access sensor settings:

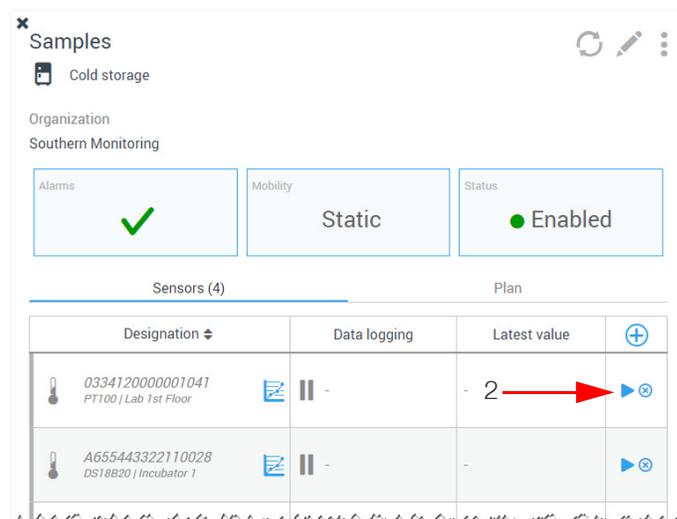


Figure 92. Choosing a sensor for data logging

- Enter your system password and click **Send** to continue.

Figure 93. Entering password to enter data logging settings screen

The **Start data logging** window opens, with several zones (in the image below):

- General sensor properties (1).
- Choosing to create a new data logging configuration or selecting an existing configuration (2).
- Entering specific parameters (3).

Figure 94. Defining data logging parameters

- Define the settings for a new data logging session as described in the next section or select an existing configuration.

Data Logging Details

If you are an Application Manager or have manager rights in your branch of the organization, you may create configuration templates with data logging settings that you can use later when programming similar sensors.

If there are no templates, the **New data logging configuration** field is selected by default. Previous settings are displayed if the sensor has already been used.

- Set the following information:

New configuration: This option is selected by default (configuration template feature not yet implemented).

Reading interval: The frequency with which the sensor records a reading. Use the drop-down list to select the value that meets needs.

Use drift/uncertainty: Select this check box if you want to take into account the deviation and uncertainty values provided for the sensor (see **Updating Parameters for Calibrated Sensors**).

- You may set low and high limits to ensure that the physical parameter(s) in the monitored environment remain(s) within a specific range of values.



CAUTION: You may enter up to three high limits and three low limits to benefit from multiple warnings if limit excursions persist. You may enter the limits in any order that you choose. The first one you enter is displayed on the Smart-View Pro data logger display.

Low limits: The lowest acceptable sensor reading (lowest temperature, lowest humidity level, etc). Any readings below the limit(s) will trigger an alarm. Click **+** to set up additional limits or **×** to remove limits you may have added.

Low delay: The duration of the reading that can remain lower than the low limit before an alarm is triggered (must be a multiple of the reading interval, in minutes).

High limits: The highest acceptable sensor reading. Any readings above the limit(s) will trigger an alarm. Click **+** to set up additional limits or **×** to remove limits you may have added.

High delay: The duration of the reading can remain higher than the high limit before an alarm is triggered (must be a multiple of the reading interval, in minutes).

If a reading exceeds these limits when data logging is running (and the delay is reached, if configured), you will be notified of the alarm by a color change on the data logger screen and outer ring, and via the system's alert mechanism.

While the data logger is in an alarm state, the Smart-View Pro data logger buzzer will also emit an audio alarm when plugged into AC power.

3. Click **Start** to begin recording sensor readings.

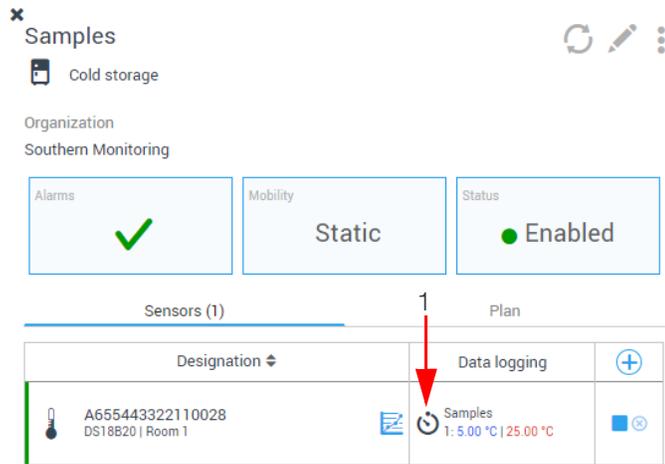


Figure 95. Equipment details with data logging in the process of starting

Data logging in the process of being started is indicated by a timer icon (🕒) in equipment details screen (1). The timer changes to (🕒) when data logging is running.

Stopping Data Logging

You may stop data logging that is currently running:

1. In the main menu, click **Equipment** (📁).
2. Click on the name of the equipment for which you want to stop data logging
3. Sensors for which data logging is in progress have a **Stop data logging** button (🛑). Click on the button (2) to stop data logging:

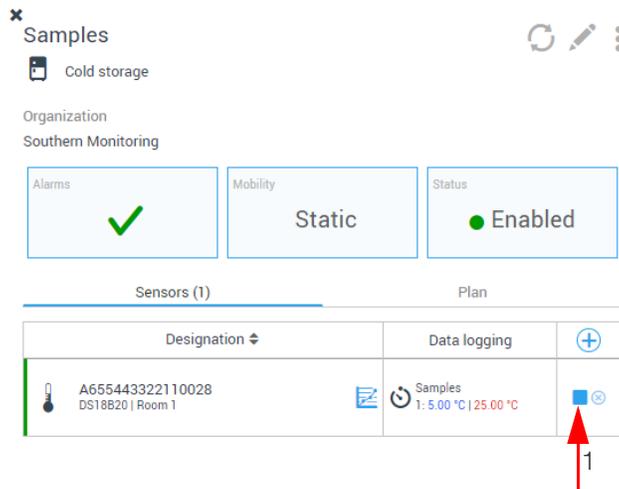


Figure 96. Stopping data logging

4. Confirm your action by entering your password in the pop-up window:

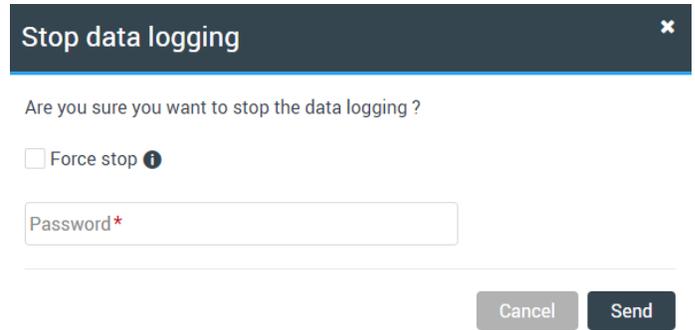


Figure 97. Entering your password to stop data logging

If you select **Force stop**, data logging is stopped at the next programmed transfer interval without retrieving the latest information from the data logger (readings or any other information). Any unretrieved information is lost.

If you do not select **Force stop**, data logging is stopped at the next programmed transfer interval, but all the information is retrieved from the data logger. The operation may take longer with this option.



CAUTION: Do not confuse the data logger's wireless data "transfer" interval, with the sensor "reading" interval configured for each sensor.

Viewing Sensor Data

Sensor data is first stored in the data logger's internal memory and then transferred wireless to the Smart-Vue Pro platform at programmed intervals. To preserve data logger battery life under normal operation, transmission to the server occurs periodically, but not at every reading. If you modify settings in Smart-Vue Pro, the values on the Smart-Vue Pro data loggers screen will be updated when the next reading interval occurs.

There are several ways to access data collected by a sensor:

Equipment details: Provides equipment details, including graphs.

Sensor list: Sensor readings displayed in a table with data export options.

To access the list of the readings recorded by a sensor:

1. Click **Readings** (📄) in the Equipment, Data Loggers or Sensors screens.
2. A color graph is displayed showing readings recorded by the selected sensor:



Figure 98. Example of a sensor graph

The **Readings** window contains two tabs (1):

Graph: The graph shows readings in a visual format over time. You can display information from several sensors running data logging in a given piece of equipment at the same time:

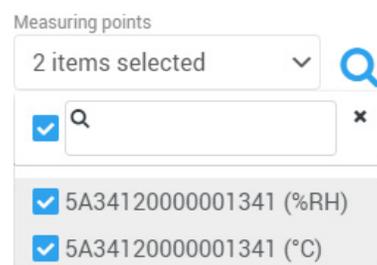


Figure 99. Readings window

Data: A list is displayed showing all data and events that occurred throughout the data logging session (readings, time-stamps, start, stop, limits exceeded, technical alarm, etc.,).

Displaying Sensor Graph

All sensor readings are shown in graph to track variations over the entire data logging cycle:

- The vertical axis shows the reading value.
- The horizontal axis shows the time scale.
- High limits are represented by red horizontal lines.
- Low limits are represented by red horizontal lines.
- Alarms (high and low limits) and events (technical alarms) are displayed in color. A key describing different colors is shown at the bottom of the graph.
- Highest, lowest, and average readings calculated during the data logging cycle are shown at the bottom right-hand side of the graph.

Selecting the Display Period

By default, the graph shows readings for the past 7 days.

To adjust the dates for the graph display:

1. Click **Start date** and/or **End date** fields (1) and choose a month, day, and year from the date selection calendars.



Figure 100. Selecting the display period

2. Click (Q) to confirm your choice and refresh the graph.
3. The two scroll buttons next to the **Start** and **End** dates (2) are used to change the period for which you would like to display graphs. Click the arrows to move back or forward by one week.

Including Events in Graph Display

You may choose to include events in the graph displays, such as “Data logging started”, “Data logging stopped”, “Low battery alert”, etc.,

To include events, click **Display events** toggle switch at the top of the window to adjust the display:

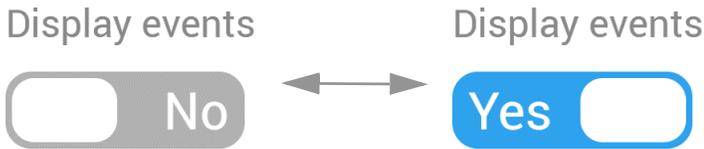


Figure 101. Event display toggle switch

Zoom and Control Tools

When you move the mouse cursor (1) over the graph, a tool-tip displays reading values and timestamps:

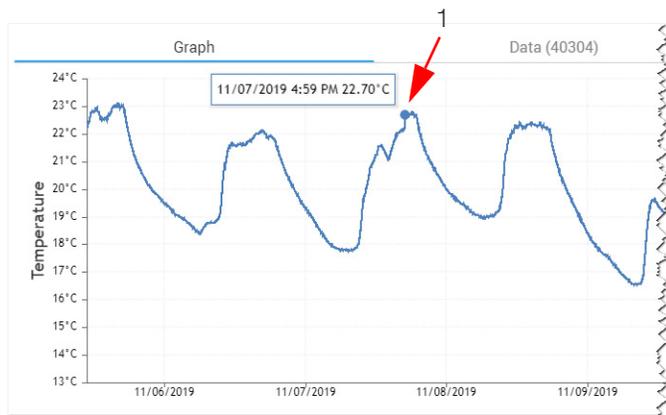


Figure 102. Reading graph with high and low limits

If **Display events** is selected, events are indicated with the appropriate background color on the graph. Move the mouse cursor over the graph to see event details (2):

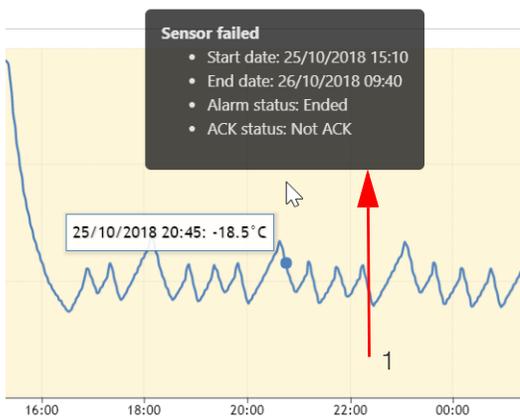


Figure 103. Graph with event details

You may zoom into the graph to examine details more closely:

1. Left-click and hold the mouse button and drag the cursor over the desired location in the graph.
2. Release the mouse button to zoom in on the selected area (2).

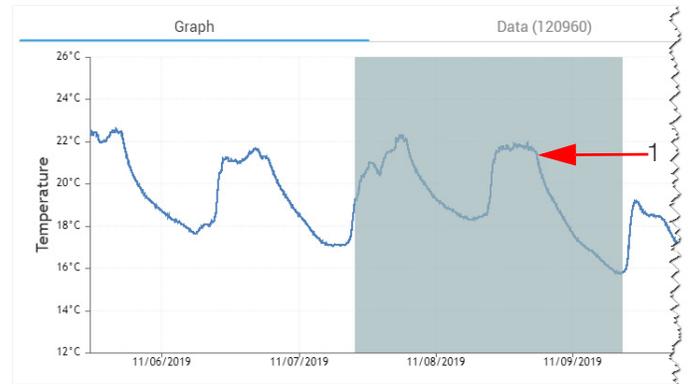


Figure 104. Selecting the zoom area

The selected area is enlarged.

3. Repeat these steps to get a higher zoom level if needed.

As you zoom in on the graph, additional tools are shown in the upper right-hand corner of the graph (3):



Figure 105. Zoom and Control tools

Icon	Description
	Enables you to use your mouse to scroll right and left through the graph. When you click on this button, the mouse cursor turns into a multi-directional arrow. Then click anywhere on the graph and drag right or left while holding the mouse button to focus on particular areas of the graph.
	Turns off panning mode.

Icon	Description
	Clears zoom or pan effects and restores the graph to its original scale.

Exporting Data

Click the **Download** button () at the top of the screen to generate a report based on the filters applied to the graph display. Select the format (PDF, Word or Excel).

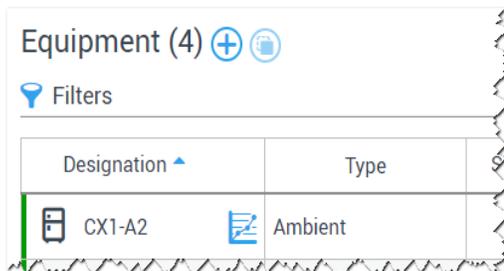
List of Sensor Readings

To display detailed sensor readings:

1. Click **Readings** icon () via the Equipment, Data Loggers or Sensors menu:



OR



OR

Data loggers (36)   

Filters

Type	Designation	S num
Smart-Vue ...	1Cyient...	E50C

Figure 106. Accessing sensor data

2. Click on the Data tab (1):

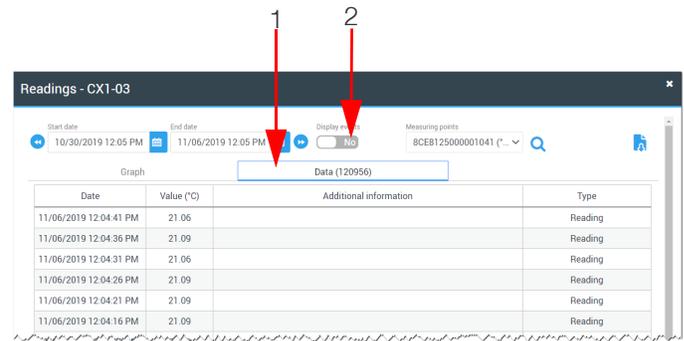


Figure 107. List of readings/events

3. You may include the list of events for the selected sensor using the **Display events (Yes/No)** toggle button (2).

The table includes the following information:

- Event date
 - Reading value
 - Event details
 - Event type
4. You may use the date range selector to adjust the dates for displaying events.

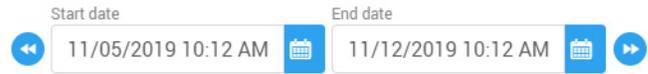


Figure 108. Selecting the date range for displayed data

Use the right and left arrow buttons ( and ) to move forwards or backwards through the calendar.

Creating Reports

Smart-Vue Pro enables you to export several different reports concerning your equipment.

Lists

Many screens in the application offer a **Download** button (), notably those presenting lists of equipment, sensors, alarms, etc. Several format options are displayed when you click on the download button (1).

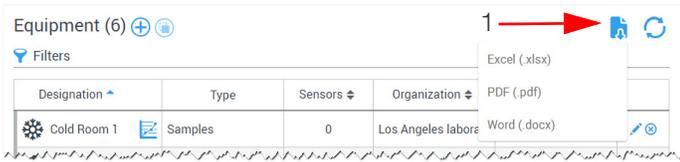


Figure 109. Data export tools

The available options are:

Excel (.xlsx): Creates XLSX format file that you can open with Microsoft Excel.

PDF (.pdf): Creates PDF printout of this table.

Word (.docx): Creates DOCX format file that you can open with Microsoft Word.

Audit Trail

The audit trail is a key feature for traceability in equipment monitoring solutions, notably with respect to the FDA’s CFR Part 11 guidelines. Smart-Vue Pro tracks every system action in a non-modifiable list that provides complete details about system events, such as:

- Alarms
- User login/logout
- Configuration changes
- New equipment/devices/sensors
- Profile updates
- Alert rule creation/modification/testing
- Data logging start/stop
- System messages
- Password changes

To view your system’s audit trail:

1. Click **Reports**  **Audit trail.**
2. Here you may click on:
 - (1) to open filter options and refine the display by choosing the types of information you want to see.
 - (2) to download an XLSX, PDF or DOCX version of the displayed list.



Figure 110. Audit trail list

3. The header fields show the event date, user, type, and description. You may sort the list by date or user by clicking on the headers using your mouse.
4. You may expand any event item by clicking on (3) for details on a particular event, as shown here:

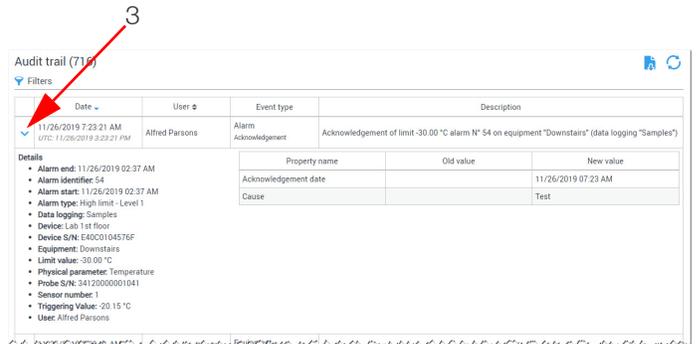


Figure 111. Complete details for a specific event

Exporting the Audit Trail

As described above in section **Lists**, you may also download audit trail information in XLSX, PDF or DOCX format. The following figure shows a sample first page from a PDF download:



Figure 112. Downloading an audit trail report

Alert Devices

The Smart-Vue Pro system supports several “electronic” alert notification mechanisms, notably:

- E-mail (native to all Smart-Vue Pro systems).
- SMS/Text message (optional via Smart-Vue Pro Alert).
- Automated voice call (optional via Smart-Vue Pro Alert).

In addition, you may use hardware alert devices via LoRaWAN® wireless technology:



Figure 113. Smart Siren (audio & visual alerts)



Figure 114. Smart Remote Contact (triggers an external smart remote contact device connect)

This chapter describes adding the hardware devices to your system.

Smart Siren

To add a Smart Siren to your system:

1. Click **Infrastructure** (📶) in the left-hand menu, then **Add** (+).
2. In the screen shown below, fill in the fields to create the new Smart Siren:

✕ Create infrastructure

Serial number*	Type Smart Siren
Name*	Inventory code
Organization Southern Monitoring	Time zone (UTC-08:00/-07:00) Pacific Time (US & Ca)
Description	

Technical alarm: None

Limits alarm: None

Dry contact alarm: None

Allow snooze

Trigger alarm after 0 minutes without communication.

Repeat alarm every 0 hours.

Trigger alarm after 0 minutes on battery mode.

Cancel Save

Figure 115. Adding a Smart Siren to your system

Serial number: Enter the serial number provided on your product.

Type: Select Smart Siren. If you select the incorrect type of device here, it will be adjusted automatically upon the system’s first connection.



CAUTION: We recommend that you test new alert devices by simulating alerts. You may test the siren directly. **Options** (⚙️) → **Simulate an alarm** once the device has been created, and/or in the full context of testing an alert rule by simulating a sensor alarm (see **Simulating a Sensor Alarm for Testing**).

Name: Enter a name to use as a reference in your system.

Inventory code: Enter an inventory code for this device (optional).

Organization: You must assign the device to a branch of your organization. Choose the level that is appropriate for your hierarchy.

Time zone: Enter the time zone in which the device is located in order to make sure that it is activated at the right moment (as defined in Alert rules).

Description: You may enter information here as desired (optional).

Technical alarm: Choose how the siren should react when receiving a technical alarm, such as a low data logger battery:

None = no reaction

Buzzer and light = Audio sound + red siren light

Light only = Red siren light only

Limit alarm: Choose the siren type when receiving a limit alarm, such as a temperature excursion (None, Buzzer and light, Light only).

Smart Remote contact alarm: Choose the siren type when receiving a limit alarm, such as a temperature excursion (None, Buzzer and light, Light only).

Allow snooze: Enables you to press the snooze button on the siren to turn off the light and/or buzzer until the next alarm occurs.

Trigger alarm after X minutes without communication:

The system checks communication with all devices periodically. If the system fails to communicate with the siren, and then fails again after “X” minutes that you specify, an alarm is triggered (such as an e-mail or text message).

Repeat alarm every X hours:

Enter the frequency with which the system should trigger an alarm again if the siren is still not communicating.

Trigger alarm after X minutes on battery mode:

If the siren switches from AC power to battery mode, an alarm is triggered after the number of minutes you specify here.

The new siren is listed in the Infrastructure list shown here:

Infrastructure (1) 

 Filters

Designation 	Identifier 	Organization	Version	Last activity 
Control Room Remote Contact Smart Remote Contact	E51100000001	Southern Monitoring		

Figure 116. Siren shown in Infrastructure list

Testing your Smart Siren

We recommend that you test any newly installed alert devices to ensure smooth operation in your context. To test the siren:

1. Click on the siren name in the list.
2. Click **Options**   **Simulate an alarm**.

Note: This alarm is useful for testing the siren device itself, outside the context of any alert rules.

3. Confirm details in the window that opens:

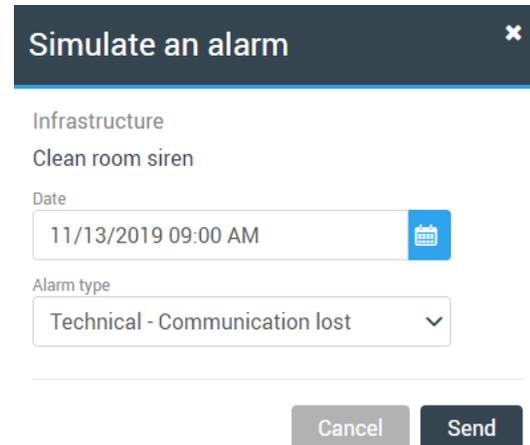


Figure 117. Details for simulating an alarm

Use the drop-down list to choose one of the available alarm types for the siren: “Technical – Communication lost”, “Technical – Lost power supply”, or “Technical – Low Battery”.

4. Click **Send** to launch the test, or **Cancel** to return to the previous window without taking any action.

Other Smart Siren Functions

Once you have added the siren to your system, you may include it in alert rules to notify users in case of alarms.

Alarms & Alerts provides details on how to configure alerts. Other functions (1) are available by clicking on the siren in the list, which opens the siren details as shown here:

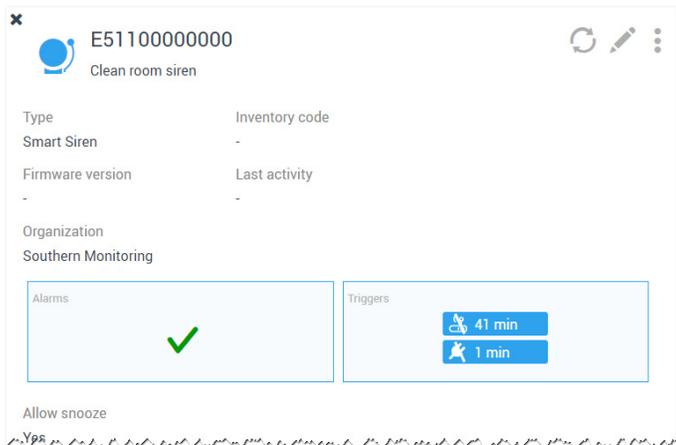


Figure 118. Additional options for Smart Siren management

Options	Category	Description
Options (⋮)	Alarm history	Opens the Alarm list so you can see any alarms involving this device (a filter is applied to show “Infrastructure” with the siren’s serial number) .
	Simulate an alarm	You may use this function to simulate an alarm as described in the previous section.
	Disable	Click to disable the siren’s functionality (in the Infrastructure list, a filter enables you to display or hide disabled infrastructure).
	Remove	Click to remove the siren from your system. The siren is marked as “removed” in the Infrastructure list. If you wish to restore the same siren, you must create it again by entering the serial number.
	Refresh	Refreshes the siren information window.
	Edit	Opens the Edit infrastructure window in which you may modify all the same details you entered when creating the siren.

Smart Remote Contact

To add a Smart Remote Contact device to your system:

1. Click **Infrastructure** (🔌) in the left-hand menu, then **Add** (+).
2. In the screen shown below, enter the fields to create the new Smart Remote Contact:

Figure 119. Adding a Smart Remote Contact device to your system

Serial number: Enter the serial number provided on your product.

Type: Select Smart Remote Contact. If you select the incorrect type of device here, it will be adjusted automatically upon the system’s first connection.



CAUTION: We recommend that you test new alert devices by simulating alerts. You may test the siren directly via **Options** (⋮) → **Simulate an alarm** once the device has been created, and/or in the full context of testing an alert rule by simulating a sensor alarm (see **Simulating a Sensor Alarm for Testing**).

Name: Enter a name to use as a reference in your system.

Inventory code: If you have an inventory code for this device, you may enter it here (optional).

Organization: You must assign the device to a branch of your organization. Choose the level that is appropriate for your hierarchy.

Time zone: Enter the time zone in which the device is located in order to make sure that it is activated at the right moment (as defined in Alert rules).

Description: You may enter information here as desired (optional).

Technical alarm: Choose how the remote contact device should react when receiving a technical alarm, such as a low data logger battery:

None = no reaction

Relay 1 = Triggers state change on relay 1 (terminals 1-2-3)

Relay 2 = Triggers state change on relay 2 (terminals 4-5-6)

Relay 1 & relay 2 = Triggers state change on both relays

Limit alarm: Choose how device reacts when receiving a limit alarm, such as a temperature excursion (None, Relay 1, Relay 2 or Relay 1 & relay 2)

Smart remote contact alarm: Choose how device reacts when receiving a limit alarm, such as a temperature excursion (None, Relay 1, Relay 2 or Relay 1 & relay 2)

Trigger relays when power supply is unplugged:

Determines relay behavior if AC power is lost. You may click on **Show connections** to see images of the different wiring options (details on relay open and closed configuration are provided below in section **Terminal Wiring**).

Trigger alarm after X minutes without communication:

The system checks communication with all devices periodically. If the system fails to communicate with the device, and then fails again after “X” minutes that you specify, an alarm is triggered (such as an e-mail or text message).

Repeat alarm every X hours:

Enter the frequency with which the system should trigger an alarm again if the device is still not communicating.

Trigger alarm after X minutes on battery mode:

If the remote contact device switches from AC power to battery mode, an alarm is triggered after the number of minutes you specify here.

The new remote contact device is listed in the Infrastructure list shown here:

Infrastructure (1) 

 Filters

Designation 	Identifier 	Organization	Version	Last activity 
Control Room Remote Contact Smart Remote Contact	E51100000001	Southern Monitoring		

Figure 120. Siren shown in Infrastructure list

Testing your Smart Remote Contact Device

We recommend that you test any newly installed alert devices to ensure proper operation in your context. To test the remote contact device:

1. Click on the siren name in the list.
2. Click **Options**   **Simulate an alarm.**

Note: This alarm is useful for testing the remote contact device itself, outside the context of any alert rules.

3. Confirm details in the window that opens:

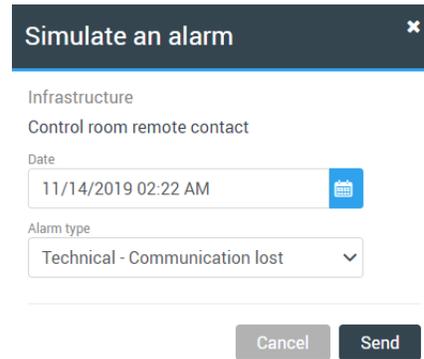


Figure 121. Details for simulating an alarm

Use the pull-down list to choose one of the available alarm types for the siren: “Technical – Communication lost”, “Technical – Lost power supply” or “Technical – Low Battery”.

4. Click **Send** to launch the test, or **Cancel** to return to the previous window without taking any action.

Other Smart Remote Contact Device Functions

Once you have added the remote contact device to your system, you may include it in alert rules to notify users in case of alarms. **Alert Devices** provides details on how to configure alerts. Other functions (1) are available by clicking on the device in the list, which opens details as shown here:

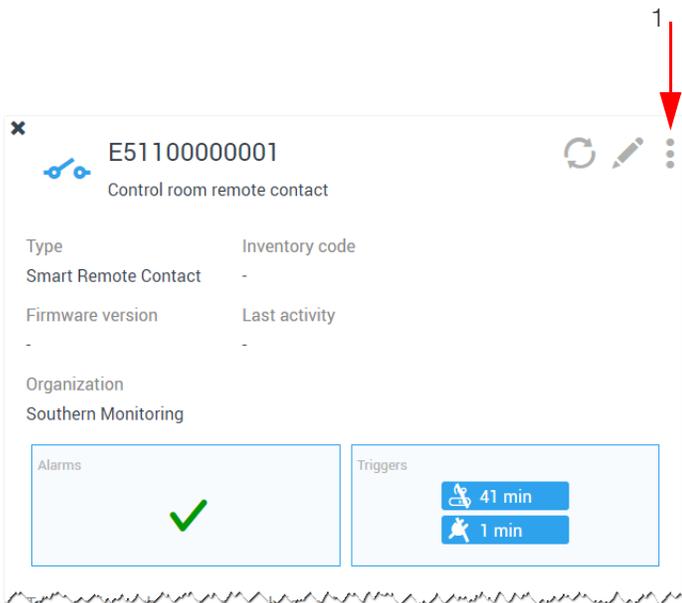


Figure 122. Additional options for Smart Remote Contact management

Table 10. Other functions for Smart Remote Contact

Options	Category	Description
Options (⋮)	Alarm history	Opens the Alarm list so you can see any alarms involving this device (a filter is applied to show “Infrastructure” with the device’s serial number)
	Simulate an alarm	You may use this function to simulate an alarm as described in the previous section.
	Disable	Click to disable the device’s functionality (in the Infrastructure list, a filter enables you to display or hide disabled infrastructure).
	Remove	Click to remove the remote contact device from your system. The device is marked as “removed” in the Infrastructure list. If you wish to restore the same siren, you must create it again by entering the serial number.
	Refresh	Refreshes the Smart Remote Contact information window.
	Edit	Opens the Edit infrastructure window in which you may modify all the same details you entered when creating the device.

Terminal Wiring

The terminal is wired as shown in the following figure (the images are shown in the Smart-Vue Pro Web application).

When "Trigger relays..." checkbox is not selected.

Default state: Pins 2&3 of relay1 and Pins 5&6 of relay2 are short circuited.

Alarm State: Pins 1&2 of relay1 and Pins 4&5 of relay2 are short circuited.

When "Trigger relays..." checkbox is selected.

Default state: Pins 1&2 of relay1 and Pins 4&5 of relay2 are short circuited.

Alarm State: Pins 2&3 of relay1 and Pins 5&6 of relay2 are short circuited.

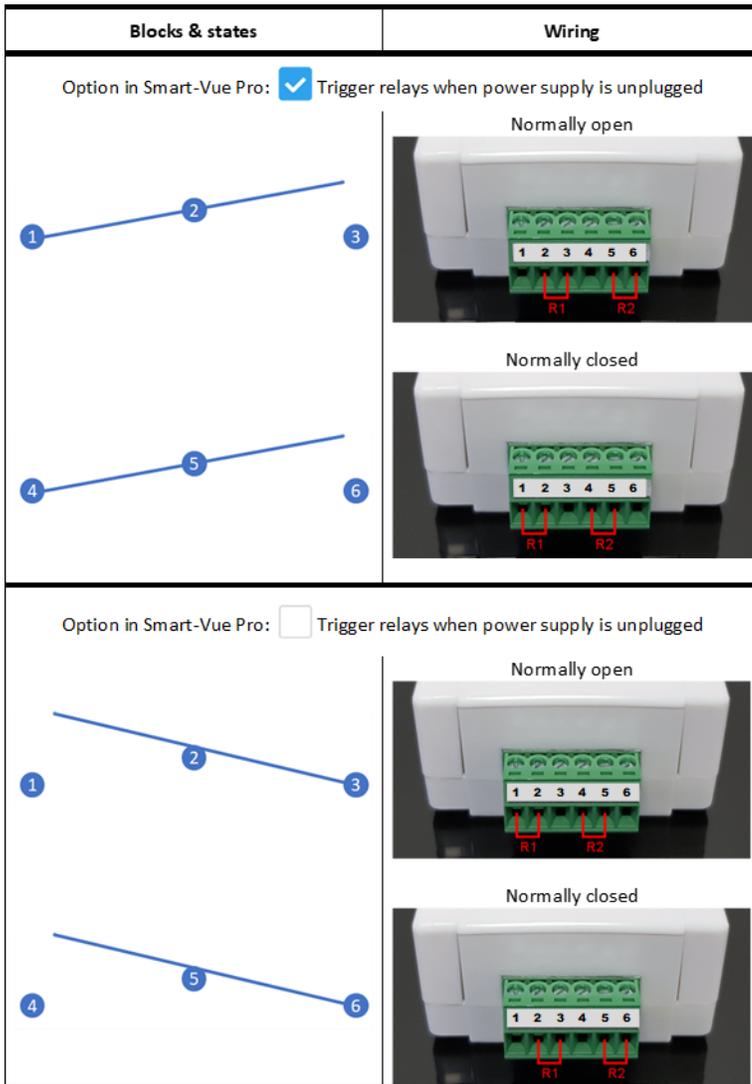


Figure 123. Terminals for open and closed relays

Connect the wires of your external devices to the terminal poles as appropriate for your needs.



CAUTION: The two blocks (1-2-3 and 4-5-6) may be cabled differently from each other.

For example, with **Trigger relays when power supply is unplugged** selected, 1-2-3 may be cabled for "Normally open" and 4-5-6 may be cabled for "Normally closed".

Alarms & Alerts

The Smart-View Pro monitoring system features an alert mechanism that notifies you if the sensors go outside their expected range or if some type of technical error occurs. For example, if a temperature sensor is configured to monitor a cold-storage unit between 2°C and 8°C, and the sensor reads a temperature of 9°C, then an alarm is shown in Smart-View Pro. The system can also send an alert to one or more users if an alarm is detected. It is therefore important to understand the distinction between alarms and alerts.

Handling Alarms

Alarms Highlighted in the Dashboard

The first place you will see an alarm visually is in the dashboard, in the **Watch Mode** screen.

When there are no errors, the equipment thumbnail is shown in green and the alarm indicator at the top right of the equipment details page displays a green OK icon (✓).

Here is a “healthy” system dashboard and equipment details window:

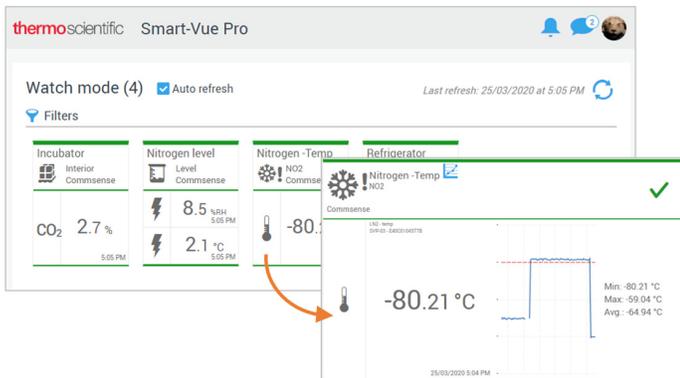


Figure 124. Dashboard and equipment details with no alarms

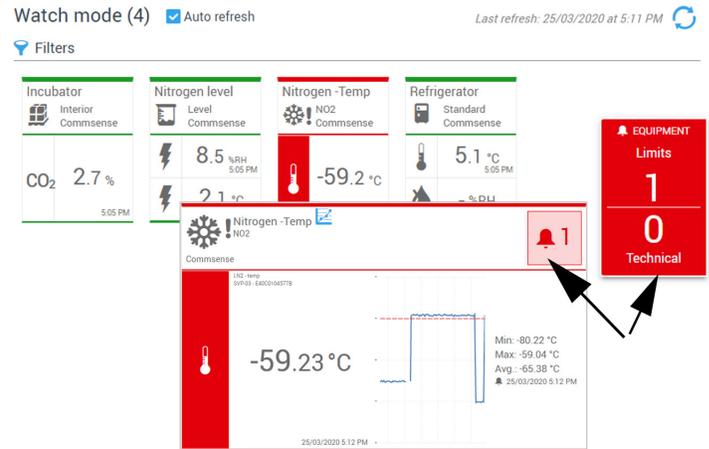


Figure 125. Dashboard and equipment details with an alarm

The alarm counters (1) increment to show:

1. On the main screen, the number of alarms currently on an equipment.
2. On the equipment details screen, the number of alarms currently on the equipment (i.e. on different sensors at the same time).

Viewing Alarm Details

When you see an equipment with errors:

1. Click on one of the alarm counters mentioned above (1) or click **Alarm** icon to open the alarm list in the left-hand menu (🔔).
 - If you click on the counter in the dashboard window or on 🔔 in the menu, you will see all current system alarms.
 - If you click on the equipment thumbnail, you will access the alarm directly (that is, with filters applied in the Alarm list to show alarm only).
2. The **Alarm** list opens and shows the alarm events for the selected equipment. Unfold alarm details by clicking on the blue arrow (↕) (2).

	Date	Designation	Equipment	Type	Status	
<input type="checkbox"/>	Start: 11/14/2019 6:40 AM	Sensor: 0334120000001041 ID: 9 - Temperature	Downstairs	High Limit	Ongoing	Acknowledge
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:21 AM	Sensor: A655443322110028 ID: 8 - Temperature	Samples	High Limit	Ongoing	Acknowledge
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:05 AM End: 11/14/2019 6:11 AM	Sensor: A655443322110028 ID: 7 - Temperature	Samples	High Limit	Ended	✓

2

3

Date	Designation	Equipment	Type	Status	
Start: 11/14/2019 6:40 AM	Sensor: 0334120000001041 ID: 9 - Temperature	Downstairs	High Limit	Ongoing	Acknowledge

Equipment	Plan	Organization
Downstairs		Southern Monitoring

Figure 126. Displaying alarm details

- Click the graph icon (📊) to open the sensor graph which shows the date/time when the alarm started.
- You may access the equipment directly by clicking on the link in the Equipment column (3).

Acknowledging Alarms

When alarms are displayed for your equipment, you must acknowledge the alarm in Smart-Vue Pro and troubleshoot the problem so that further alarms do not continue to occur.



CAUTION: It is important to note that the alarm status is still shown in red in the Smart-Vue Pro dashboard until the next reading transmitted by the data logger no longer indicates the alarm condition.

More importantly, acknowledging the alarm in the software does not actually fix any physical or technical problem encountered by the data logger and the equipment could still be in an alarm state.

Follow these instructions to acknowledge an alarm:

1. Click **Alarms** (🔔) in the main menu, or on the alarm counter in the dashboard or equipment details screens.

2. The alarm management screen displays equipment with alarms:

	Date	Designation	Equipment	Type	Status	
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:40 AM	Sensor: 0334120000001041 ID: 9 - Temperature	Downstairs	High Limit	Ongoing	Acknowledge
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:21 AM	Sensor: A655443322110028 ID: 8 - Temperature	Samples	High Limit	Ongoing	Acknowledge
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:05 AM End: 11/14/2019 6:11 AM	Sensor: A655443322110028 ID: 7 - Temperature	Samples	High Limit	Ended	✓

Figure 127. Alarm management and acknowledgment screen



To organize the display of unacknowledged alarms, click **Filters** and select **Not ACK** from the **Acknowledgment** drop-down list, or select any other criteria depending on your needs.

3. To acknowledge a single alarm, click [Acknowledge](#) next to the alarm you want to acknowledge.
4. You may acknowledge several alarms at once by clicking on their check boxes in the first column:

	Date	Designation
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:40 AM	Sensor: 0334120000001041 ID: 9 - Temperature
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:21 AM	Sensor: A655443322110028 ID: 8 - Temperature
<input checked="" type="checkbox"/>	Start: 11/14/2019 6:05 AM	Sensor: A655443322110028

Figure 128. Acknowledging several alarms at the same time

5. Click [Acknowledge](#) at the top of the screen. If one of the alarms in your selection has already been acknowledged, it will be deselected automatically and thus not affected by the current acknowledgment.
6. The **Alarm acknowledgment** pop-up is shown on the right-hand side of the screen:

✕ Acknowledge alarm

Alarm type: High limit Alarm status: Ongoing

Cause: ▼

Corrective action:

Preventive action:

[Cancel](#) [Save](#)

Figure 129. Enter the details for acknowledging an alarm

- To acknowledge the alarm, you must enter an incident cause in the **Cause** field. Any text you enter in this field is stored and may be reused in the future by clicking on the drop-down list (1).
- The **Corrective action** and **Preventive action** fields are optional and can be used to include further information with the acknowledgment.
- Click **Save** to save your changes (and confirm by entering your password) or **Cancel** to discard changes and return to the previous screen.
- Once the alarm has been acknowledged, the green check mark (✓) is shown on the alarm line.



Remember, acknowledging an alarm does not necessarily eliminate any physical problems at the sensor, data logger or equipment levels.

Alarm Report Subscriptions

Smart-Vue Pro can generate alarm reports automatically and send them at programmed intervals.



Users with Application Manager rights can assign reports to other users, whereas regular users can only create reports for themselves, without being able to assign them to other users.

Scheduling a Report

To schedule a report and assign users:

- Click **Alarms** (🔔) → Schedule a report (📄).

	Date	Designation	Equipment	Type	Status	
<input type="checkbox"/>	Start: 19/03/2020 11:31:49 End: 24/03/2020 10:37:19	Data logger: E40CO...	Refrigerator	⚠️	Ended	Acknowledge
<input checked="" type="checkbox"/>	Start: 19/03/2020 10:31:49 End: 19/03/2020 10:40:58	Sensor: 5A3412000...	Refrigerator	↓	Ended	✓
<input checked="" type="checkbox"/>	Start: 19/03/2020 10:17:14 End: 19/03/2020 10:26:14	Sensor: 5A3412000...	Refrigerator	↓	Ended	✓

Figure 130. Scheduling an alarm report to send to one or more users

The report is created with default parameters (not editable as of this writing):

Schedule a report ✕

The report will be scheduled with the following parameters:

Report type
Alarms
Organizations
All
Types
All
Acknowledgment
All
Status
All

Cancel Previous Next

Figure 131. Default report parameters

- Click **Next** and edit details for the report:

Schedule a report ✕

Name*

Periodicity
 Weekly Monthly

Monday Tuesday Wednesday Thursday
 Friday Saturday Sunday

At 00:00

Data range
Previous 7 days

Cancel Previous Next

Figure 132. Set up the report to meet your needs

Name: Enter a name to refer to this report.

Periodicity: Select whether the report should be produced on a weekly or monthly basis.

Weekly: Select one or more days to generate the report:

Weekly Monthly

Monday Tuesday Wednesday Thursday
 Friday Saturday Sunday

Figure 133. Report printed on selected days every week

Monthly: Select the day of the month to print the report:

Weekly Monthly

The

Figure 134. Report printed on the specified day once a month

At: Indicate the time of the report to be produced.

Data range: Specify the time period for which you want to include data (from “Previous 7 days” to “Previous month”).

- Users with “User” rights in the system can only create reports for themselves, so they must click **Save** at this point, or on **Previous** to return to the scheduling screen, or **Cancel** to close the window without saving changes.
- Users with “Application Manager” rights must click **Next** to select the users (including themselves) who will receive the report in PDF format via e-mail:

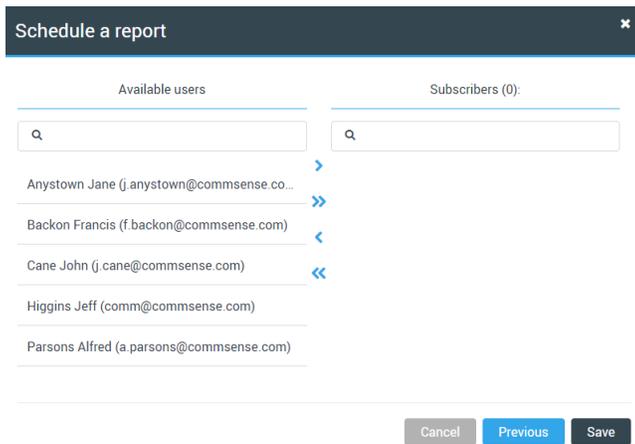


Figure 135. Include users in report distribution

To add users:

Double-click on a user-name or click to move the selected user from **Available** to **Subscribers**.

To select multiple users at a time, press and hold the **Ctrl** key and click on each of the users you want to select. Click on to add them to the list of **Subscribers**.

To select all available users, click on

To remove users:

To remove a user from a report, double-click on a user-name in the list of **Subscribers** (on the right-hand side) or click on a user and click on to move the user to the list of Available users on the left-hand side.

To remove multiple users, press and hold the **Ctrl** key and, while holding down the key, click on each of the other users

you want to remove. Then click to move them to the list of **Available** users on the left.

To remove all users, click

- Click **Save** when done, **Previous** to return to the scheduling screen, or **Cancel** to close the window without saving your changes.

Editing Scheduled Reports and Subscriptions

You may edit and update reports that you have created (as described in the previous section) and reassign users to those reports.

To modify a report:

- Click **Reports** **Report subscriptions**.

Type	Name	Scheduling parameters	Owner	Subscribers
Alarms	Weekly report	At 08:00, on the first Monday - Previous 7 days	Higgins Jeff	1

Figure 136. Include users in report distribution



CAUTION: The user who created the reported is indicated as the “Owner”. You may change the owner via Options as described below.

- To run the report right now, click .
- To delete the report, click .
- To access report configuration, click anywhere on the line with the report (1).

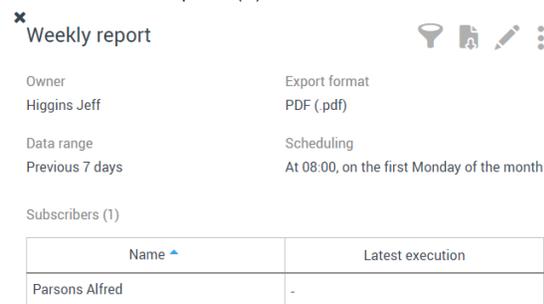


Figure 137. Report summary screen

This screen contains several options for managing the report:

Table 11. Options for Managing Report

Options	Category	Description
		Filter – Displays the parameters used to filter data and establish the report. As of this writing, the parameters may not be edited.
		Generate the report – You may click on this icon to generate the report at any time.
		Edit – Enables you to change scheduling options for the report. The options here are the same as those used to create the report (as described in the previous section).
	Options – Contains the following items:	
		Change owner – The report owner is allowed to manage report details. If the report is assigned to multiple subscribers, ownership may only be assigned to another Application Manager. A regular User cannot be the owner of a report that is sent to other people.
		Edit subscribers – Allows Application Managers to edit the list of users who receive the report via e-mail.
		Disable – Deactivates automatic generation of the report. You may Enable the report subsequently if you want (to show reports that have been disabled, open Filters at the top of the screen and select “Show inactive”)
		Remove – Completely removes the report from the system.

Alert Notifications

Smart-Vue Pro offers several alert notification mechanisms:

- By default, every system provides alert notification via e-mail and every user profile is required to include an e-mail address.
- **Smart-Vue Pro Alert** is an optional web-based solution that adds SMS/text messaging and automated voice call support. You must enter a valid license key in order to use Smart-Vue Pro Alert notification features, as described in **Managing License Keys**.

Smart-Vue Pro can send alerts to one or more users configured in Call Groups. Each user is configured with an e-mail address to be used for this purpose. To implement alerts, you must therefore create one or more Call Groups and create alert rules to determine the equipment and data loggers for which those groups will receive alert e-mails.

Configuring Alert Rules

An alert is a notification sent by the system to users when the system observes an alarm condition or potential problem in the monitored environment.

Alert rules can be configured and assigned to different call groups in order to notify, by telephone and/or e-mail, users designated to handle an alarm category. You may define different criteria for one or more nodes in your organization, or according to the geographical areas you cover or sensor categories. Alert rules are based on “Time slots” and “Weekends and holidays”, defined in the system. These entities enable you to handle a variety of different cases, including Day/Night and weekend modes.

Follow these steps to configure an alert rule:

1. Click **Configuration**  → **Alerts** → **Add alert rule** .
2. The **Create alert rule** screen is displayed:

Figure 138. Setting up an alert rule (Step 1)

- Fill the in fields according to your needs. The fields marked with a red asterisk (*) are required:

Name: Assign a friendly name to identify the alert rule.

Description: Enter a description of the alert, for your information.

Alarm type: Select one or more type of alerts (Limits, Dry Contact (for future) or Technical).

Limit alerts are the high and low limit values that you may configure for sensor readings.

Technical alerts concern technical issues with your sensors and receivers, such as low battery and communication errors. A given rule can only apply to one or the other of these types of alerts.

Items in organizations: Choose the relevant organization(s) and/or department(s).

Alerts for: Select the type of equipment for which alerts will be sent: Infrastructure, mobile equipment, Shipments, or Static equipment.

Criticality: Select the criticality level(s) for which the alert rule applies (critical and/or not critical) as configured for equipment.

Repeat alert if Check this option to set the repeat frequency for this alert if the alarm condition persists and specify the desired time interval.

- Click Next to set up the users who will be alerted and when, as described below.

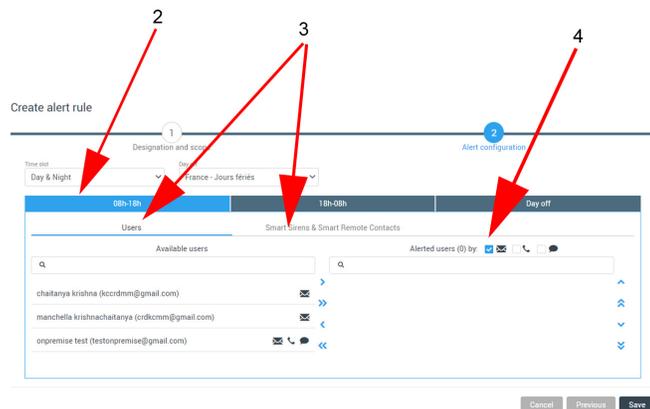


Figure 139. Assigning users to time periods in an alert rule (Step 2)

In the above image, time periods are shown on three tabs (2). Remember to assign users and/or alert devices to each period.

You may configure notifications to be sent to users or devices (3), notably Smart Sirens and/or Smart Remote Contact configured in your system.

- Select the **Time slot** and **Days off** for the alert:

Time slot: Use the drop-down menu to select the alarm period from the available options. Other time slots may be added as described in **Time Slots**.

Days off: Use the drop-down list to select Days off periods (such as weekends and holidays). Other days off may be added as described in **Setting Days off (Weekends and Non-working days)**.

Alert Rules with Users

Add or remove users in the **Users** tab (3) of the screen as shown in **Figure 138** and **Figure 139**:

- Determine how users will be alerted (4): via e-mail (✉) automated voice call (☎) or SMS / Text message (💬). Click the check box → to assign one or more alert types.
- To add users: Double-click on a user-name or click ➤ to move the selected user from Available to Alerted. To select multiple users at a time, press and hold the Ctrl key and click on each of the users you want to select. Click on ➤ to add them to the list of Selected users. To select all available users, click ➤➤.
- To remove users: To remove a user from a rule, double-click on a user-name in the list of Alerted users (on the right-hand side) or click on a user and click on ⏪ to move the user to the list of Available users on the left-hand side. To remove multiple users, press and hold the Ctrl key and while holding down the key, click on each of the other users you want to remove. Then click on ⏪ to move them to the list of Available users on the left. To remove all users, click ⏪⏪.
- To change the contact sequence: Users are contacted in the order in which they are listed in the alert rule. The top user is contacted first, etc. You may click ⏴ and ⏵ icons to move the selected user up or down one line at a time. You may click on the ⏴⏴ and ⏵⏵ icons to move the selected user to the very top or bottom, respectively.
- Click **Save** to save the rule (or add alert devices, as described in the next section), click **Previous** to return to step 1 or click **Cancel** to close this window without saving any changes.

Alert rules with Smart Siren or Smart Remote Contact Devices

To add or remove alert devices, click **Sirens & Smart remote contacts** tab (5) of the alert rule screen:



Figure 140. Assigning alert devices to time periods in an alert rule (Step 2)

1. Select one or more sirens (🔊) or smart remote contact devices (🔗) in the “Available” column.
2. To add devices:
Double-click on a device name or click ➤ to move the selected device from **Available** to **Alerted**.
To select multiple devices at a time, press and hold the **Ctrl** key and click on each of the devices you want to select. Click ➤ to add them to the list of “Triggered” devices. To select all available devices, click ➤➤.
3. To remove devices:
To remove a device from a rule, double-click on a device name in the list of **Triggered** devices (on the right-hand side), or single-click on a device name and click ⬅ to move the device to the list of **Available** devices on the left-hand side.
To remove multiple devices, press and hold the **Ctrl** key and, while holding down the key, click on each of the other devices you want to remove. Then click ⬅ to move them to the list of **Available** devices on the left.
To remove all devices, click ⬅⬅.
4. To change the trigger sequence:
Devices are triggered in the order in which they are listed in the alert rule. The top device is triggered first, etc.
You may click ⬆ and ⬇ icons to move the selected device up or down one line at a time.
You may click on the ⬆ and ⬇ icons to move the selected device to the very top or bottom, respectively.

5. Click on **Save** to save the rule (or add users as described in the **previous** section), on **Previous** to return to step 1, or on **Cancel** to close this window without saving any changes.

Modifying or Deleting Alert Rules

To disable or change the conditions or recipients of an alert rule, click on click **Configuration (⚙️) → Alerts**.

The alert rules you defined earlier are shown in the table with the configuration parameters and characteristics of the equipment involved.

1. To modify an alert rule, click Edit (✎) for the rule you wish to modify:

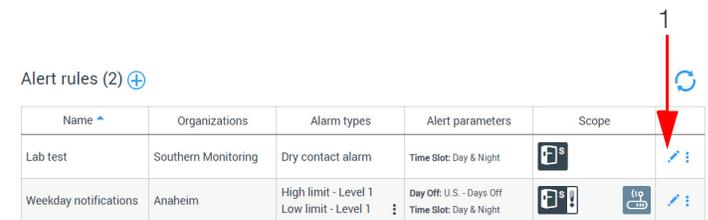


Figure 141. List of alert rules saved in the system

2. Make the desired changes (with all the same options as described in the previous sections when creating the rule).
3. Click **Save** to save your changes or **Previous** to return to the previous screen without saving changes.
4. For more options, click **Options (⋮)** for a drop-down menu with the following options:

Create using as template: This option enables you to create a new alert rule based on the currently selected rule.

Simulate an alarm: Enables you to validate the alarm scenario by sending a test alert using the selected rule. No alert is listed in the Alarm list, but people are contacted and/or devices are triggered as if an alarm occurred.

Disable: Deactivates the alert rule. You may reactivate it at any time selecting the alert rule and clicking on “Enable” in the drop-down menu.

Delete: Deletes the alert rule from the system.

Time Slots

As described earlier in **Configuring Alert Rules**, you may configure specific time slots during which your system will send alerts to users. These time slots may then be used in the context of one or more alert rules.

To add a time slot:

1. In the main menu, click **Configuration**  → **Time slots**.
2. Click **Add time slot** ()

Time slots (1) 

Name 	Description 	
Day & Night	Standard office hours	 

Figure 142. Adding a new time slot

3. The **Create time slot** window opens on the right-hand side of the screen:

x Create time slot

Name*

Description

Slots

Name* From  To  

Figure 143. Time slot details

4. Enter the details of the new time slot. The fields marked with a red asterisk (*) are required:

Name: Enter a name for the time slot.

Description: Enter a description of the time slot.

Slots: Enter a name for the time slot and then, define the start and end time. These parameters are used by the application to determine the start and end of the alarm period.



To avoid confusion, all times are indicated in local time of the user who will be receiving the alert or the device to be triggered.

1. Click **Add** () to add up to a total of three time slots, if needed.
2. Click **Save** to save your changes or **Cancel** to discard changes and return to the previous screen.

Modifying or Deleting Time Slots

To change the time slots of your alert rules:

1. In the main menu, click **Configuration**  → **Time slots**.

Time slots (2) 

Name 	Description 	
Day & Night	Standard office hours	 
Weekends	Not at the office	 

Figure 144. Time slot list

2. Click **Edit** () to modify a time slot.
3. Make the necessary changes.
4. Click **Save** to save your changes or **Cancel** to discard changes and return to the previous screen.
5. If you want to delete a time slot, select the time slot and click **Delete** ()

Setting Days off (Weekends and Non-working days)

Smart-Vue Pro enables you to manage non-working days such as weekends, public holidays and other days off. You may use this feature to fine-tune your system's alert rules.

To add days off:

1. In the main menu, click **Configuration**  → **Days off**.

Days off (2) 

Name 	Description 	
France - National holidays		 
U.S. - National holidays		 

Figure 145. Adding days off (non-working days)

- Click **Add days off (+)**, which opens the **Create days off** window:

Figure 146. Entering details for days off

- Fill the fields according as desired. The fields marked with a red asterisk (*) are required:

Name: Enter a name as a reference for the day(s) off. You will be able to use this name when you create alert rules.

Description: (Optional) You may enter a description of the day off for information purposes.

Weekend configuration: Set the start and end times that are considered as part of the weekend, for example, for on-call staff. These settings are used by the application to determine the start and end of the weekend alarm period.

- Click **Next** → **+ Add days off**

Figure 147. Adding days off to be used in alert rules

- Select dates in the **From** and **To** fields.
- If the day is to be considered as a day off every year, click the **Yearly** check box.
- You may remove a day off by clicking on **Remove (-)** next to the item to be deleted.
- Click **Save** to save the day(s) off in the system, **Previous** to return to the previous screen without saving changes or **Cancel** to discard your changes.

Modifying or Deleting Weekends and Days off

To edit a weekend or day off:

- In the main menu, click **Configuration** ⚙️ → **Days off**

Name	Description	
France - National holidays		
U.S. - National holidays		

Figure 148. List of days off in your system

- Select the line you want to change, click **Edit** (✎).
- Make the desired changes.
- Click **Save** to save your changes or **Cancel** to discard changes and return to the previous screen.
- If you want to delete a day off configuration, select the desired line and click **Delete** (✖).

Getting Help

User Manuals

User manuals for the Smart-Vue Pro web application and related products are available directly from the application.

Click **Help** → **Online help** to download the latest documents:

Help

Read the latest online user guides to for detailed information on using the Smart-Vue Pro solution and getting the most from its features.

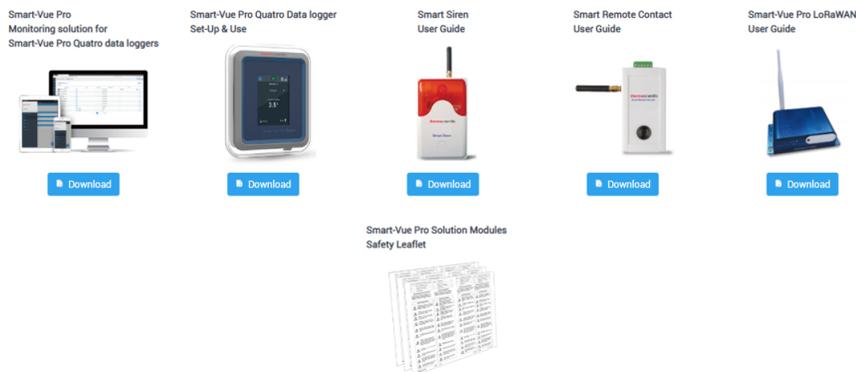


Figure 149. Access to Smart-Vue Pro and companion product user guides

Viewing Smart-Vue Pro Version Release Notes

If the Smart-Vue Pro web application has been updated since the last time you logged in, the release notes concerning the new version are displayed on your screen:



Figure 150. Latest release notes

This message is not displayed for new users.

You may click **Do not show these notes again** to stop the message from being displayed when you log in again. Click **Close** to close the window and continue with Smart-Vue Pro.



You may show all Smart-Vue Pro release notes at any time by clicking **Help** → **Release notes**.

Appendix

Alarms

Table 12. Alarm Details

Type	Meaning	Web App alarm color
High Limit Alarm	Sensor input reading is higher than high limit value set.	Red
Low Limit Alarm	Sensor input reading is lower than low limit value set.	Red
Sensor Fail	Input Sensor has failed.	Red
Invalid sensor	When a different sensor is replaced during data logging.	Red
Communication lost	When data logger or infrastructure (LoRaWAN) is not communicating as per the setting in web app i.e., web app triggers an alarm after X minutes without communication with Data logger or Infrastructure.	Red
Power lost	When data logger is not under AC power and communicating on battery power, as per the setting in web app i.e., web app triggers an alarm after X minutes without AC power supply.	Red
Unexpected stop	When you restore power, the data logger is operational but the data logging session is not restarted. Then SmartVue Pro web application triggers the Unexpected Stop alarm to notify you that your equipment is not being monitored.	Red
Low battery	The battery on the SVP Quatro is low.	Red

FAQ

1. I just installed a new data logger. It seems to be working fine, but no information is updated in the Smart-Vue Pro dashboard.

- Information is automatically sent to the system at the end of each programmed “transfer period”. The transfer period is generally longer than the sensor reading period, so you may simply need to wait for the cycle to begin in order to have data in the web application.
- On the Smart-Vue Pro data logger, you may use the **Synchronize** feature to trigger communication between the system and the data logger.
- You may also refresh the application screen using the **Refresh** button () that is present throughout Smart-Vue Pro.

2. I used the Smart-Vue Pro web application to activate data logging on a Smart-Vue Pro data logger. The data logger now displays the temperature and session name as expected, however, I still do not see any readings in the web application. What should I do?

If the data logger screen shows information other than “Data logger waiting for configuration”, it means that communication from the server was successful. The edges of the tile in the Watch Mode screen (in the web application) should no longer be gray. Just wait until sensor readings are uploaded at the next transfer interval.

3. My data logger was data logging and lost power (unplugged or batteries removed). However, when the power was restored, the web application showed an “Unexpected Stop alarm”. What could be the reason?

When you restore power, the data logger is operational but the data logging session is not restarted. The data logger therefore still needs your attention and hence, the Smart-Vue Pro web application triggers the “Unexpected Stop alarm” to notify that the equipment is not being monitored.

4. Can a data logger be removed which is currently running a data logging session?

Yes, you can remove a data logger from your Smart-Vue Pro system even if data logging is running. You should only do this in cases where there is a problem with the data logger, or if a data logger had to be swapped while

still running. In those cases, the Smart-Vue Pro web application can no longer stop the data logging session via software and you may need to remove the data logger from the system.

5. Where can I find information about using the Smart-Vue web application with my Smart-Vue Pro data loggers running in Bluetooth mode and the Smart Connect mobile application?

The Smart Connect User Guide provides details about using the mobile application in conjunction with the web application.

6. Are all features available when using the Smart-Vue Pro web application and the Smart Connect mobile application for my data loggers in Bluetooth-only mode?

Most viewing and display features are the same, regardless of whether data is collected via a LoRaWAN receiver or pushed by your Bluetooth-enabled device running Smart Connect. Some configuration features are specific for LoRaWAN operation and do not apply when using Smart Connect, such as “Trigger alarm after n minutes without communication.”

7. Why an unexpected stop alarm shows in ongoing state although it is acknowledged?

Unexpected stop event occurs when the data logging is in progress and battery/power supply is removed.

When the battery (or plug your power supply) is placed, the device reboots. However, all data loggings are stopped so the data logger sends this alarm to notify the customer that your product (ex: vaccine) is no longer monitored.

8. Why doesn't it show any error while creating a company even if the license key or last digit is changed?

Validation of license is done when one accepts and tries to create company. To ensure the security, it is a best practice to check the license key on server side and not on front side (javascript). The error is shown in the final page and user won't be able to create a company when last digit of the license key is changed.

9. Explain the following error “Data logger configuration mismatches server”.

If the physical data logger has a different state than the status on the web app, this alarm is created. For example, starting a data logging, performing a factory reset and communicating with the server.

10. What is the significance of system user and support user in Audit Trail records?

System user: When the action is performed by the application.

Note: It is not a real user (sending an email, etc.)

Support user: When the technical support uses the special account to perform some actions.

11. What is the significance of setting time interval for communication lost separately in Infrastructure and Data logger pages?

Infrastructure (gateway) communicates with server in every 10 seconds, however, it updates its activity in every 5 minutes.

A LoRa gateway takes 5 minutes to start. Communication lost alarm is triggered after two missing reading intervals +1 boot duration. Hence, a 15 minutes time interval is defined for communication lost in Infrastructure page.

12. Why does the time stamp show as 01/01/0001 before starting the data logging in web app watch mode screen?

Until the first data push to cloud we find the time stamp as 01/01/0001. Once the data is synced to cloud we observe the related time stamp.

WEEE Compliance

WEEE Compliance. This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. It is marked with the following symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on our compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive are available at www.thermofisher.com under Services & Support.

Great Britain



WEEE Konformität. Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2012/19/EU erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliederstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter www.thermofisher.com unter Services & Support.

Deutschland



Conformità WEEE. Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2012/19/EU in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Thermo Fisher Scientific ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Thermo Fisher Scientific con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Thermo Scientific che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito www.thermofisher.com in Servizi e Supporto.

Italia



Conformité WEEE. Ce produit doit être conforme à la directive euro-péenne (2012/19/EU) des Déchets d'Equipements Electriques et Electroniques (DEEE). Il est marqué par le symbole suivant. Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur www.thermofisher.com sous Services et Assistance.

France



Contact Information

Thermo Fisher Scientific products are backed by a global technical support team ready to support your applications. We offer cold storage accessories, including remote alarms, temperature recorders, and validation services.

Visit www.thermofisher.com or call:

Countries	Sales	Services
North America	+1 866 984 3766	(800) 438-4851
India	1800 22 8374, +91 22 6716 2200	+91 22 6716 2200
China	+800 810 5118, +400 650 5118	+8621 68654588
Japan	+81 3 5826 1616	+81 3 3816 3355
Australia	+61 39757 4300	1 300 735 292
Austria	+43 1 801 40 0	+43 1 801 40 0
Belgium	+32 53 73 42 41	+32 2 482 30 30
France	+33 2 2803 2180	+33 2 2803 2180
Germany	0800 1 536 376, +49 6184 90 6000	0800 1 536 376
Italy	+32 02 95059 552	+39 02 95059 552, 432 254 375
Netherlands	+31 76 579 55 55	+31 76 571 4440
Nordic/Baltic/CIS	+358 9 329 10200	+358 9 329 100
Russia	+7 812 703 4215	+7 812 703 4215
Spain/Portugal	+34 93 223 09 18	+34 93 223 09 18
Switzerland	+41 44 454 12 22	+41 44 454 12 12
UK/Ireland	+44 870 609 9203	+44 870 609 9203
New Zealand	+64 9 980 6700	+64 9 980 6700
Other Asian Countries	+852 2885 4613	+852 2885 4613
Countries not listed	+49 6184 90 6000	+49 6184 90 6000

Thermo Fisher Scientific Inc.
 275 Aiken Road
 Asheville, NC 28804
 United States

Find out more at thermofisher.com/cold