

Thermo Integration Agent Quick Start Guide

This Quick Start Guide describes the steps to install and configure the Thermo Integration agent data service and other supplementary components. Refer to the Thermo Integration Agent Data Service Installation Guide (IG) sections noted for more detailed instructions.

The installation of Thermo Integration agent data service involves the following three tasks:

- Pre-installation tasks [IG: Section 1.1]
- Installation tasks [IG: Section 1.2]
- Post-installation tasks [IG: Section 1.3]

Task 1: Pre-installation [IG 1.1]

Before you begin the Thermo Integration agent data service installation, you must first install the prerequisite hardware and software:

- Integration HPIC System Instrument PC
- Chromeleon CM7.2.SR4 (build 8179) or higher

If the user mode is enabled in Chromeleon, then

the exclude-user settings in the Chromeleon Administration Console must first be configured. See IG Sections 1.1.3 – 1.1.5.

Use the Thermo Password utility after installation to store the excluded-user credentials for Thermo Integration agent data service use. See IG Section 1.4 for details.

Task 2: Installation [IG 1.2]

The installation involves the following tasks:

- Thermo Integration agent data service installation [IG 1.2.1]
- Thermo Remote desktop server silent installation [IG 1.2.2]

Thermo Integration agent data service installation [IG 1.2.1]

Before installing Thermo Integration agent data service, ensure that the minimum required version of Chromeleon is installed on the machine to avoid exiting with an error message.

For installation, use the ThermoInstaller CD and perform the following steps in the installation wizard:

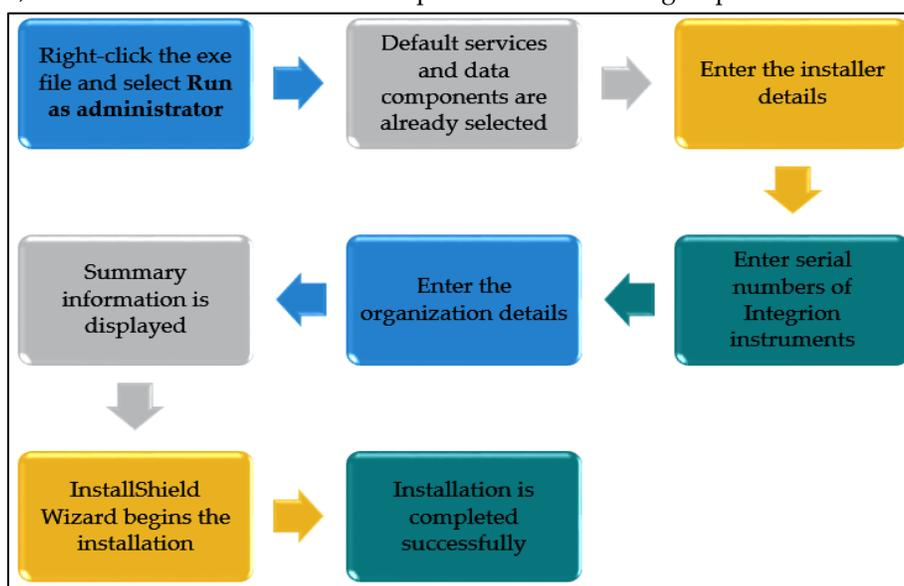


Figure 1: Installation Wizard step

Thermo Remote Desktop Server silent installation [IG 1.2.2]

Thermo Integrion agent data service installer performs a silent installation of the remote desktop server. No other action is required.

Task 3: Post-installation [IG 1.3]

Complete the following post-installation tasks:

- Verify the installation [IG 1.3.1]
- Verify the remote desktop server [IG 1.3.2]

Verify the installation [IG 1.3.1]

To verify if the Thermo Integrion agent data service is successfully installed, perform the following actions:

- Verify that the installer received an email
- Verify creation of the **ThermoGateway** folder at the following location:
C:\Program Files (x86)\
- Verify creation of the **xGate.log** file in the following folder:
C:\Program Files (x86)\
ThermoGateway\

- Verify creation of the **ContactInfo.txt** file in the following folder:
C:\Program Files (x86)\
ThermoGateway\
- Verify that the organization details entered for Thermo Chromeleon Gateway and assets in the installation are reflected in the **Provisioning.xml** file found in this folder:
C:\Program Files (x86)\
ThermoGateway\
- Verify that **ThermoDesktopServer**, **ThermoChromeleonGateway**, and **ThermoIntegrionDataService** services are visible with Started status and Automatic startup type in the **Services** program.

Verifying the remote desktop server installation [IG 1.3.2]

To confirm successful installation of the remote desktop server:

- Verify creation of the **Server** folder at the following location:
C:\Program Files (x86)\
ThermoGateway\Desktop\Server\
- You have already verified that the **ThermoDesktopServer** service is visible in the **Services** window.

For detailed information about installation and configuration of Thermo Integrion agent data service, see *Installation Guide*. If you have any queries or need support related to this guide, please send an email ULS.RemoteSupport@thermofisher.com.