

MaxQ High Performance Console

Incubated Orbital Shaker*

Operating and Maintenance Manual 7030435 Rev. 5

Visit us online to register your warranty www.thermoscientific.com/labwarranty



from cover:

- * Triple counter-balanced, single eccentric drive mechanism (U.S. Patent #5,558,437)
- * Horizontal, HEPA-filtered airflow design (U.S. Patent #5,577,837)
- * Test tube rack (U.S. Patent #5,632,388)

Models covered by this manual:			
Model	Number	Voltage, Frequency	Temperature Control
SHKE435HP	435	120VAC, 60Hz	Incubated
SHKE436HP	436	230VAC, 50Hz	Incubated

MANUAL NUMBER 7030435

5	40722	7/10/17	Added gas springs statement to Maitenance	bpg
4	41343	6/13/17	Moved F-Gas statement	bpg
3	41343	5/02/17	Added F-Gas statement, removed declaration of conformity	bpg
2	40230	8/1/16	Changed platform part number from 238054 to 238083 - pg 1-7	CCS
1	40139	4/15/15	Updated warranty information	CCS
0	28036/0S-751	9/8/14	Release 13 (435), Release 8 (436) - new control board	CCS



Important Read this instruction manual. Failure to read, understand and follow the instructions in this manual may 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. ▲

Warning Use MaxQ SHKE435HP/SHKE436HP Orbital Shaker to process non-flammable materials only!

Warning Grounding circuit continuity is vital for safe operation of this shaker. Never operate this unit with the grounding circuit disconnected. ▲

Material in this manual is for information 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 out of or related to the use of this manual.

When translated into other languages, the US English version of this manual is binding.

©2014 Thermo Fisher Scientific. All rights reserved.



Important operating and/or maintenance instructions. Read the accompanying text carefully.



Potential electrical hazards. Only qualified persons should perform procedures associated with this symbol.



Equipment being maintained or serviced must be turned off and locked off to prevent possible injury.



Hot surface(s) present which may cause burns to unprotected skin, or to materials which may be damaged by elevated temperatures.



WEEE Compliance: Thermo Fisher Scientific has contracted with companies for recycling/disposal in each EU Member State. For further information, send an email to weee.recycle@thermofisher.com.

- ✓ Always use the proper protective equipment (clothing, gloves, goggles, etc.)
- ✓ Always dissipate extreme cold or heat and wear protective clothing.
- ✔ Always follow good hygiene practices.
- ✓ Each individual is responsible for his or her own safety.

Do You Need Information or Assistance on Thermo Scientific Products?

If you do, please contact us 8:00 a.m. to 6:00 p.m. (Eastern Time) at:

1-740-373-4763 1-800-438-4851 1-877-213-8051 http://www.thermofisher.com service.led.marietta@thermofisher.com www.unitylabservices.com Direct Toll Free, U.S. and Canada FAX Internet Worldwide Web Home Page Tech Support Email Address Certified Service Web Page

Our **Sales Support** staff can provide information on pricing and give you quotations. We can take your order and provide delivery information on major equipment items or make arrangements to have your local sales representative contact you. Our products are listed on the Internet and we can be contacted through our Internet home page.

Our **Service Support** staff can supply technical information about proper setup, operation or troubleshooting of your equipment. We can fill your needs for spare or replacement parts or provide you with on-site service. We can also provide you with a quotation on our Extended Warranty for your Thermo Scientific products.

Whatever Thermo Scientific products you need or use, we will be happy to discuss your applications. If you are experiencing technical problems, working together, we will help you locate the problem and, chances are, correct it yourself...over the telephone without a service call.

When more extensive service is necessary, we will assist you with direct factory trained technicians or a qualified service organization for on-the-spot repair. If your service need is covered by the warranty, we will arrange for the unit to be repaired at our expense and to your satisfaction.

Regardless of your needs, our professional telephone technicians are available to assist you Monday through Friday from 8:00 a.m. to 6:00 p.m. Eastern Time. Please contact us by telephone or fax. If you wish to write, our mailing address is:

> Thermo Fisher Scientific (Asheville) LLC 401 Millcreek Road, Box 649 Marietta, OH 45750

International customers, please contact your local Thermo Scientific distributor.

Table of Contents

Section 1	Installation	1-1
	Pallet Hold-down Shipping Brackets	1-2
	Location	1-2
	Install the Cabinet Stand-off Bolts	1-2
	Chamber Drain	1-3
	Condensate Drain	1-3
	Foot Pedal	1-3
	Installing the Platform	1-4
	Assemble Flask Clips	1-6
	Install Flask Clips	1-7
	Install Test Tube Holders	1-8
	RS-232 Interface Connector	1-9
	Connect the Remote Alarm	1-11
	Lid Security Lock	1-11
	Connect to Electrical Power	1-12
Section 2	Operation	2-1
	Control Panel Operation	2-2
	Quick Start-Up	2-3
	Factory Default Settings	2-3
	Change Temperature, Speed & Time Settings	2-4
	Change Temperature	2-4
	Change Speed	2-5
	Change Time	2-5
	Change from Hold to Countdown	2-5
	Shaker Alarms	2-6
	Overtemp Shutdown	2-7
	Cycle Complete	2-7
	Power Failure	2-8
	RPM Tracking	2-8
	Check Belt	2-8
	Sensor Fault	2-9
	I emperature High or Low	2-9
	Platform Stalled	2-9
		2-10
	Turn the Audible Alarm On and Off	2-10
	Set Alarm Limits	· · · 2-11
	Jul 1 Mallil Lillillo	

Section 2 (continued)	Remote Alarm System	-14
	Cycle Complete	-14
	Power Failure	-14
	RPM Tracking 2	-15
	Check Belt	-15
	Sensor Fault	-15
	Temperature High or Low 2	-15
	View Total Operating Hours	-16
	Heat %	-16
	Software Version	-17
	Overtemp Sensor Readings 2	-17
	Menu Map	,-18
Section 3	Maintenance	3-1
	Gas Springs	3-1
	Platform and Cabinet Cleaning	3-1
	Control Panel	3-1
	Cleaning or Replacing Air Filter	3-1
	Preventive Maintenance	3-2
Section 4	Service	.4-1
	Alarms and Alarm Conditions	4-1
	Change HEPA Filter	4-2
	If the Shaker Will Not Operate	4-2
	Spare Fuses	4-2
	Circuit Boards	4-4
	Temperature Sensors	4-5
	Blower Fan Motors and Heating Elements	4-5
	Heater Element Circuit Breaker	4-6
	Tune the Cabinet	4-6
	Service the Drive Belt	4-6
	Calibrate Speed (RPM)	4-7
	Calibrate the Temperature	4-8
	Alarm Messages	4-9
Section 5	Specifications	5-1
Section 6	Parts List	6-1
Section 7	Electrical Schematics	7-1
Section 8	Warranty Information	8-1

Section 1 Installation



Figure 1-1. MaxQ SHKE435/SHKE436HP Console Incubated Orbital Shaker

The shipping carton should be inspected upon delivery. When received, carefully examine for any shipping damage before unpacking. If damage is discovered, the delivering carrier should specify and sign for the damage on your copy of the delivery receipt.

Open the carton carefully making certain that all parts are accounted for before packaging materials are discarded. After unpacking, if damage to any of the contents is found, promptly report it to the carrier and request a formal damage inspection.

Important Failure to request an inspection of damage within a few days after receipt of shipment absolves the carrier from any liability for damage. Call for a damage inspection promptly.

	 Model SHKE435HP and SHKE436HP Console Orbital Shakers are shipped with the following materials: 2 - Keys for the lid lock (packaged and attached to outside of unit) 1 - T-handle 5/32" hex wrench 2 - Platform alignment studs ¼-20 1 - Shaker platform 6 - Grade 8, 5/32" hex socket flat head screws (provided with platform) 2 - 3/4" Open end wrench 2 - ¼-20 x 7" Stand-off bolts with rubber caps 1 - Phillips screwdriver for flask clip installation and removal 1 - 8¾" Phillips screwdriver for flask clip installation and removal 1 - Line cord (country of destination)
Pallet Hold-down Shipping Brackets	To secure the console shaker to the shipping pallet, hold-down brackets are attached to slots in both sides of the cabinet. The brackets are fastened to the wood pallet with lag screws (Figure 1-2). Figure 1-2. Bracket
Location	Install the shaker on a firm, level surface in an area free of dust and dirt. To allow for lid opening, the back of the shaker must be at least 4 ½ inches from the wall. As the electrical plug is the "mains disconnect" for the unit, the electrical wall outlet must remain accessible at all times.
Install the Cabinet Stand-off Bolts	To maintain the 4½ inch minimum distance between the rear of the unit and the wall to allow for unobstructed lid opening, two ¼-20x7" bolts are included in the parts bag.
	Screw the bolts into the threaded holes on the back of the shaker cabinet (Figure 1-3). The bolts should be screwed in by hand to the limit of the threads. Further tightening is unnecessary. Put the protective rubber caps, also supplied in the parts bag, over the heads of the bolts. $Figure 1-3$. Rear View



Figure 1-4. Chamber Drain Hose Location

Chamber Drain A drain is provided in the bottom of the chamber for convenience when cleaning or removing spills (Figure 1-4). A clear vinyl hose and plastic valve is connected to the drain and accessed by removing the front grille assembly and the lower front panel. The grille is removed by gently pulling it off. It is held in place by six push-in type retainers.

To remove the cabinet panel located in back of the grille, remove the six Phillips screws; three on the top and three on the bottom. It may also be necessary to loosen the two left side Phillips screws which hold the foot pedal assembly to the shaker frame.

Condensate Drain A 1/4" stainless steel condensate drain is located on the back of the shaker to remove any water that may collect in the air ductwork. Refer to Figure 1-5.



- Figure 1-5. Drain Location
- **Foot Pedal** The lid of the console shaker is counterbalanced for ease of opening and closing. A foot pedal on the lower right of the front of the cabinet is also provided for operator convenience and ready access to the chamber.

Installing the Platform

Note If your platform is already installed, skip this section and continue to next section.

Caution Remove the shipping bracket and install the shaker platform before plugging in or attempting to operate the unit. ▲

After removing the orbital mechanism shipping bracket and installing the platform, remove this protective decal from the control panel to begin shaker operation.









To protect the shaker's orbital mechanism during shipment, a sheet metal shipping bracket (Figure 1-7) is installed and must be removed before the unit can be operated. Using a $7/16^{\circ}$ and $9/16^{\circ}$ hex wrench, remove the three $1/4^{\circ}$ and the single $3/8^{\circ}$ screws. Retain this hardware for future shipping.

This shaker accommodates either a 5/16" nominal heavy-duty, 29.5" wide x 18" front-to-back platform.

All shaker platforms are attached to their orbital mechanisms with six 1/4-20 hex socket flathead screws Grade 8. These screws are hardened and should not be exchanged with any other screw type. The 5/32" hex socket wrench, included with the shaker, must

be used when attaching the platform. Refer to Figure 1-8.

> Figure 1-8. T-Handle Wrench and Hex Socket Head Screw

Caution Do not attempt to use a Phillips head screwdriver. ▲



Installing the Platform (continued)

1. Insert the two 1/4-20 alignment pins into the two mounting holes identified in Figures 1-9 and 1-10.



Figure 1-9. Hole Locations (Universal Shaker Platform Shown)



Figure 1-10. Hole Locations

- 2. Rotate the drive mechanism until the four mounting holes generally match the holes in the platform.
- 3. Place the platform onto the shaker and over the alignment pins.

Installing the Platform (continued)

- 4. Move the platform in an orbital motion until one or more of the center mounting holes are located.
- 5. Insert the hex socket head screws as the four holes are located. Do not tighten the screws.
- 6. <u>Remove</u> the ¹/₄-20 alignment pins and replace them with the remaining two hex socket screws.
- 7. Tighten all screws using the T-handle wrench.

Caution Use only the hex socket flat screws to fasten the platform, and only the T-handle wrench to tighten the screws. Torque these screws to 10 ft-lbs. Check these screws monthly if the unit is operated at or near maximum speed (525 RPM). ▲

Assemble Flask Clips

Each Flask Clip up to 6.0 liters in size comes with a metal spring that must be installed onto the clip. For flask clips through 500 ml, insert the end of each spring into the holes at the top of the clip leg (Figure 1-11).



Figure 1-11. Clip Mounting Screw



The 2 liter, 2.8 liter, 4 liter, 5 liter, and 6 liter Flask Clips use two sets of metal springs and rubber spring tubes. On these larger clips, the springs are installed by hooking their ends together as illustrated in Figure 1-12. The upper spring and spring tubes should be installed prior to mounting the clip to the platform. The lower spring and spring tubes, however, are placed around the bottom of the clip legs after the flask clip is fastened to the platform.

Note that the rubber spring tubes are placed between the clip legs.

Install Flask Clips

Model SHKE435HP and SHKE436HP shakers accommodate glassware in numbers and sizes from ninety-one 25 ml flasks to four 6 liter flasks. All platforms have mounting holes for flask clips and test tube racks made by other manufacturers. Listed below are the dedicated platform kits available for these shakers.

Dedicated Platform Number	No. of Clips	Flask Size (ml)	Springs per Clip	Screws per Clip
238017	91	25	1	1
238018	91	50	1	1
238019	39	125	1	1
238051	30	250/300	1 (w/ 1 lg pad)	1
238021	24	500	1	1
238022	15	1 L	1	5
238023	12	2 L	2 (w/ 10 tubes)	5
238024	6	4 L	2 (w/ 10 tubes)	5
238083	5	5L	2 (w/ 12 tubes)	5
238025	4	6 L	2 (w/ 12 tubes)	5
238026	6	2.8 L	2 (w/ 10 tubes)	5
238020	40	250/300	1 (w/ lg pad)	1

Table 1-1. Dedicated Platform Kits

Flask clips can be attached anywhere on the platform and flasks can be inserted into any flask clip as the counter-balanced design of these shakers compensates for unbalanced loads.

The flask clips are supplied with the proper screws and can be attached to the platform with a standard Phillips screwdriver or the screwdriver provided with the unit.

Figures 1-11 and 1-12 illustrate the installation of the flask clips. Note that clips for 1, 2, 2.8, 4, 5, and 6 liter flasks use five screws. The 250/300ml flask clip has an adhesive-backed flask cushion pad that is installed on the flat base of the clip body. A hole is provided in the pad for the mounting screw.

Install Test Tube Holders

The Accessory Test Tube Racks and Test Tube Rack Holders are available in four sizes and are listed in Table 1-2.

Table 1-2. Test Tube Racks and Holders

Part No.	Description
950040	Test Tube Rack, 10-13 mm size
950060	Test Tube Rack, 16-20 mm size
600074	Test Tube Rack, 21-25 mm size
600075	Test Tube Rack, 26-30 mm size
600076	Adjustable-Angle Test Tube Holder w/ Rack, 10-13mm
600077	Adjustable-Angle Test Tube Holder w/ Rack, 16-20mm
600078	Adjustable-Angle Test Tube Holder w/ Rack, 21-25mm
600079	Adjustable-Angle Test Tube Holder w/ Rack, 26-30mm
600088	Universal Adjustable-Angle Test Tube Holder, 10-25mm
600089	2 Tier Micro-Plate Rack
600090	3 Tier Micro-Plate Rack
194024	#10-24 pan head Phillips screws for mounting holders to platforms

All the Test Tube Rack Holders are adjustable into seven positions, swinging and locking at 15°, 30° and 45° in either direction. Figure 1-13 illustrates the Test Tube Rack Holder with rack in place. To remove the rack, spread the metal tabs on either end of the holder and lift out the plastic Test Tube Rack.



Figure 1-13. Test Tube Rack Holder

Install Test Tube Holders (continued)

To install the Test Tube Rack Holder onto the shaker platform, remove the rack and rotate the swing-bed of the holder 90° by pulling the knobs of the locking pins on either end of the holder outward. The pins are locked outward by turning the knob 1/4-turn (Figure 1-14). Attach the tray to the platform with the screws provided.



Figure 1-14. Rotate Rack

RS-232 Interface Connector

The MaxQ Console Orbital Shaker is equipped with an RS-232 Serial Communication Interface for the remote transmission of data. An RJ-11 telephone style connector is located on the left side of the incubator. A cable with RJ-11 plugs and an RJ-11 to DB-25 adapter are required. Refer to Figure 1-18 for connector locations on the shaker back panel. Figure 1-15 identifies the RS-232 and Remote Alarm pin contacts.

The data is "dumb terminal" formatted, which permits interfacing with either a computer or a serial printer.



Figure 1-16. Remote Alarm Output to Screw Terminal Connection

RS-232 Interface Connector (cont.)



Figure 1-17. Pin Connections

Three wires are used for the RS-232 interface:

- 1. Transmit data (/TXD) pin 2 DB-25 connections
- 2. Receive data (/RXD) pin 3 DB-25 connections
- 3. Signal ground (GND) pin 7 DB-25 connections

The data format is:

The data transfer sequence is transmitted in the following format. X refers to the numerical time, speed and temperature.

(NUL)XXX:XX(H)(SP)(SP)XXXRPM(SP)(SP)XX.XC(SP)(LF)(CR)(EOT)

NUL	Null character (0)
SP	Space (32)
LF	Line feed (10)
CR	
EOT	End of transmission (4)
H for tin	

RS-232 Interface Connector (cont.)

Model SHKE435HP and SHKE436HP shakers transmits= time, RPM and temperature information one minute after power is first applied to the unit, then every 60 minutes thereafter unless the shaker receives either a <Ctrl><Q> or a <Ctrl><S>.

The shaker's microprocessor responds to two commands from the remote:

<Ctrl><Q> (XON)

The shaker will immediately transmit time, speed, and temperature data upon receiving a <CTRL><Q> and will reset the 60 minute data transmission interval timer.

<Ctrl><S> (XOFF)

The shaker will stop serial data transmission upon receiving a <Ctrl><S> until a <Ctrl><Q> is received or power is cycled.

Connect the Remote Alarm

IMPORTANT USER INFORMATION

CAUTION! Stored product should be protected by a redundant 24 hour/day monitoring system with alarm capability. An interconnect jack and thermocouple are installed for centralized monitoring, should on-board system fail.

Lid Security Lock

An internal SPDT relay is provided to monitor alarms and is connected by a RJ-11 (telephone style) jack on the rear of the cabinet. The remote alarm provides NO (normally open) and NC (normally closed) output. Figure 1-16 identifies the pin contacts and may be wired to a central remote alarm location or to an independent alarm system. Figure 1-15 identifies the pin contacts. Figure 1-18 shows the location of the Remote Alarm Connector.

A modular to modular cable (Stock No. 190388) and an RJ-11 telephone style terminal converter (Stock No. 190392) or equivalent may be used to convert the remote alarm output to a screw terminal connection. Refer to Figures 1-16 and 1-17.

To protect the contents of the shaker or prevent tampering or unauthorized access, a security lock is located on the right side of the lid (Figure 1-18). Two keys for this lock are in the parts package attached to the outside of this unit when shipped.



Connect to Electrical Power

Connect the line cord to the power inlet in the back of the unit.

See the serial tag on the side of the unit for electrical specifications or refer to the electrical schematics at the end of this manual.

Caution Connect the orbital shaker to a grounded dedicated circuit. The On/Off switch is the mains disconnect device for the orbital shaker. Position the unit so the switch is easily accessible. ▲

Section 2 **Operation**



Figure 2-1. Front View

Model SHKE435HP and SHKE436HP shakers are microprocessorcontrolled incubated orbital console shakers designed to accommodate a wide variety of flasks, test tubes and other glassware. The control system is easily programmed and stores the user-defined time, temperature and speed settings in battery-supported memory that remain even when the shaker is turned off and unplugged.

The platform speed controller continuously adjusts for line voltage fluctuations and provides smooth transitions with consistent control. The circuitry is designed to slowly bring the platform up to speed and down to a stop to prevent liquid splashing from flasks or test tubes.

An insulated lid with viewing port is counter-balanced for easy opening by hand or foot pedal. A safety interlock requires that the lid be closed for the drive motor, circulating fans and heating elements to operate.

Caution It may take up to one minute to bring the platform up to full speed. Never leave the shaker unattended when starting it.

Caution Make sure all flasks and test tube racks are firmly seated in the clips and check the security of the flask clip and platform attachment screws monthly.

Caution Do not operate the shaker at maximum speed without a load. \blacktriangle

Control Panel Operation

The control panel on these units has a liquid crystal display and eight operating buttons which are identified by word and symbol. During programming, the up and down arrows increase and decrease the numerical values of time, platform speed, or temperature. Press and hold either arrow to cause the values to scroll in that direction; hold for more than five seconds to increase the scrolling speed.

When changing the system configuration, the down arrow advances the display to the next screen while the up arrow returns the display to the previous screen. Pressing the Time, RPM or Temperature button selects the parameter above it to be changed, while the up and down arrows increase and decrease the numerical values, respectively, or toggle between two different options. Pressing and holding either arrow will cause the values to scroll in that direction; holding for more than five seconds will increase the scrolling speed.

The START button begins platform operation as defined by the Time and Speed setpoints, while the STOP button halts the platform. Chamber temperature control begins upon power-up as defined by the Temp setpoint.

The alarm indicator and alarm silence button complete the shaker control panel. When in alarm, the unit sounds an audible warning and flashes the three red indicators. Depending upon the error detected, pressing the Silence button turns off the audible alarm. However, the three red indicators continue to flash until the alarm condition is corrected. For most alarms, the audible warning will sound again in about fifteen minutes if the condition persists.

The alarm features are discussed in more detail in the "Shaker Alarms" section of this manual.



Figure 2-2. Control Panel

Quick Start-Up Caution If the unit is shipped or stored in very cold conditions, allow the unit to warm to ambient temperatures before using. ▲

Caution This unit should be operated by trained personnel only, as described in this manual. All appropriate personal protective equipment should be worn as required. ▲

At power-up, one of the the screens at right will appear on the display for 10 seconds (where X.XX is the current software revision) before it

Software Version # 435/436 REL X.XX

shows the Actual and Setpoint times, speeds and temperatures similar to those illustrated in Figure 2-2. For convenience, this is called the Operating Screen throughout this manual.

Initially, the Actual values along the top of the liquid crystal display will differ from the Setpoint values shown along the bottom. The Actual numbers will change as the unit continues to operate.

- Time With the time set at Hold, the time showing in the upper left portion of the LCD will begin to count upward, showing the total operating hours and minutes. The system will reset to 00:00 whenever the unit is stopped and restarted, using the Stop and START buttons. The system will not reset if the unit is turned off and on using the power switch, or if the shaker door is repeatedly opened and closed.
- **Speed** The speed shown in the upper center portion of the LCD will indicate the present platform speed. It will display zero RPM at rest and will gradually rise to the setpoint speed after the START button is pressed and the platform begins to rotate.
- **Temperature** The temperature shown in the upper right portion of the LCD will indicate the ambient temperature inside the shaker and will gradually move toward the setpoint value.

Factory Default Settings

The values shown in Figure 2-2 are factory default settings. Other factory settings are shown in the table below.

Function	Default
Audible Alarm	ON
RPM Tracking Limit (fixed)	5 RPM
Temperature Tracking Limit	10°C
Over Temperature Shutdown	83°C - 85°C
All Remote Alarms	ON

Table 2-1. Factory Settings

Factory Default Settings (continued)	The Console Shaker has been shipped from the factory with the following default settings:	
	Time: When the shaker is turned on for the first time, the liquid crystal display will show 00:00H. (Hold time) This means the unit is set to record accumulated operating time. Any programming changes in the Time settings are made in increments of five minutes.	
	Speed: The display shows the unit ready to operate at 25 RPM. Programming changes in speed are made in increments of 1 RPM. However, if the up or down arrows are held for about two seconds, t display will scroll in that direction.	
	Temperature: The display shows the operating temperature set at 37°C. Changes to the Temperature program settings are made in increments of 0.1°C. However, if the up or down arrows are held for about two seconds, the display will scroll in that direction.	
	The Console Shaker can be easily programmed to meet the most demanding laboratory requirements using its microprocessor-based technology. The following sections outline the procedures for changing the settings and programming the control system.	
Change Temperature, Speed & Time Settings	All programming or setting changes start from the Operating Screen as typically illustrated in Figure 2-2.	
	The instructions to program the Model SHKE435HP and SHKE436HP shakers are written in a step-by-step format. For convenience, the instructions begin and end at the Operating Screen.	
	Note At any time during programming or changing configuration settings, if no control panel buttons are pressed for about fifteen seconds, the display automatically returns to the Operating Screen, storing and acting upon any changes made. New settings are also stored and acted upon immediately when an arrow button is pressed. ▲	
Change Temperature	 Press the button beneath the temperature setpoint (Temp °C). The Run temperature value will begin to flash. 	
	2. Press the up or down arrows to set the new Run temperature in 0.1°C increments. Hold either arrow button to scroll. However, if the up or down arrows are held for about two seconds, the display will scroll in that direction.	
	3. Press the temperature button again to return to the Operating Screen. The temperature can be set over a range of 5.0 to 60.0C. However, Model SHKE435HP and SHKE436HP shakers may not control temperature properly if the temperature is set less than 10.0C above ambient termperature.	



- **Change Speed** 1. Press the button beneath the speed setpoint. The speed value will begin to flash.
 - 2. Press the up or down arrows to set the new speed in 1 RPM increments. Hold either arrow button to scroll. However, if the up or down arrows are held for about two seconds, the display will scroll in that direction.
 - 3. Press the speed button again to return to the Operating Screen. The speed can be set over a range of 25 to 550 RPM.
 - **Change Time** Model SHKE435HP and SHKE436HP shakers manage operating time in two ways:



- **Hold** When time is set to Hold, the value shown in the Actual portion of the display represents total operating time and may be reset at the operator's convenience. The shaker will continue to count upwards even if the console lid has been repeatedly opened and closed, or turned off and on with the power switch. The time will, however, reset to 00:00 when the STOP button is pressed and the unit then restarted by pressing the START button.
- **Countdown** When the Hold setpoint is changed to Countdown entering a time value in hours and minutes, the platform will operate for that period and automatically stop. The display will show the total time in the Setpoint segment and the operating time remaining in the Actual part of the display, as the microprocessor counts down to zero.

Change from Hold to Countdown

1. Press the button beneath the time setpoint. Hold will begin to flash.





- 2. Press either arrow to access the countdown time setpoint. The last preset time setpoint will begin to flash.
- 3. Press the up or down arrows to set the desired operating time in five minute increments. Hold either arrow button to scroll in that direction. However, if the up or down arrows are held for about two seconds, the display will scroll in that direction.

Change from Hold to Countdown (continued)

4. When the desired elapsed time is set (8 hours, 30 minutes in this example), press the time button to return to the Operating Screen. Pressing the START button will start the platform and begin the countdown sequence. When 00:00 is reached, the platform will automatically stop and the Cycle Complete alarm will sound.



Figure 2-4. Time Set

Shaker Alarms The MaxQ SHKE435HP/SHKE436HP control system monitors and provides alarms for nine operating parameters.

Table 2-2. Alarm Parameters			
Parameter	Alarm Message	Remote Alarm	
Overtemp Setpoint Status	Overtemp Shutdown	No	
Cycle Status	Cycle Complete	Yes	
Loss of Input Power	Power Failure	Yes	
RPM versus Setpoint	RPM is High, RPM is Low	Yes	
Drive Belt Integrity	Check Belt	Yes	
Temp Sensor Integrity	Main Temp Sensor, Over Temp Sensor	Yes	
Temp Control Status	Temperature is High, Temperature is Low	Yes	
Platform Movement Status	Platform Stalled	No	
Motor Drive Board Input Power Integrity	Check Fuse	No	

Both audible and visual alarm warnings for these nine parameters are provided by the shaker. Visual flashing of the three diagonal indicator lights on the control panel, a progression of alarm messages on the display, and an audible tone alerts the operator that an alarm condition has occurred, or currently exists.

Shaker Alarms (continued)	For convenience, the audible tone is muted by pressing the Silence button, but rings back in about 15 minutes, for most alarms, if the alarm condition is still present. However, the alarm indicator lights and alarm messages continue until the alarm condition is corrected by the operator. After the root cause of the fault has been corrected, pressing the Silence button will clear the alarm message from the display and stop the alarm indicator lights from flashing. The audible tone will ring back in about 30 minutes for the Check Belt and Check Fuse alarms if the alarm condition is still present. These alarm messages clear from the display when the unit is turned back on after correcting the alarm condition. The audible alarm feature may be turned off to suit operator or laboratory needs. As discussed in the Configuration section of this manual, the audible alarm feature may be turned off to suit the operator and laboratory needs. Refer also to the Alarm Message/Corrective Action chart in the Service section of this manual.
Overtemp Shutdown	Overtemp Shutdown alerts the operator that the overtemp setpoint has been exceeded by a few tenths of a degree. Actual 08:41 250 37.0 Setpoints Overtemp Shutdown
	The Overtemp Shutdown message displays and the heaters are turned off, but the platform and the blowers continue to operate.
	In the alarm state, the audible alarm is muted by pressing the Silence button, but rings back in about 15 minutes, if the alarm condition is still present. However, the message and indicator lights persist until the fault is corrected. Afterward, the remaining relevant alarm components are cleared by pressing the Silence button.
Cycle Complete	Cycle Complete alerts the operator that Actual 00:00 00 37.0 the end of the count-down running time Setpoints Cycle Complete has been reached.
	The Cycle Complete message shown displays and the platform stops.
	Press the Silence button to clear the alarm message from the display screen and mutes the audible alarm.

Power Failure	Power Failure alerts the operator that Actual 00:00 00 37.0 electrical power to the unit was Setpoints Power Failure interrupted, under specific operating conditions.
	While the system returns to normal operation when power is restored, the alarm message remains and the audible tone continues to sound to alert the operator. Both the display message and the audible tone are cleared by pressing the Silence button.
	Note The alarm will not occur if the power failure is less than 15 seconds in duration, while the unit is shaking. ▲
	If power is interrupted for more than 1½ hours while the unit is turned on but not shaking, a Power Failure alarm will occur. The purpose of the alarm in this case is to alert the user that an extended duration power failure occurred during the Hold interval after a timed shaking operation, or during a period of incubation only. This alarm will also occur any time the unitis turned on after being turned off for more than 1½ hours (such as when the unit is shipped from the factory, or when it is returned to use after a period of storage).
RPM Tracking	RPM Tracking alerts the operator by either alarm message shown below that the platform speed has varied ±5 RPM. In the alarm state, the audible alarm is muted by pressing the Silence button, but rings back in about 15 minutes, if the alarm condition is still present. However, the message and indicator lights persist until the fault is corrected. Afterward, the remaining relevant alarm components are cleared by pressing the Silence button.
Check Belt	Check Belt alerts the operator that the Actual O8:41 00 37.0 drive belt may have broken, is slipping Setpoints Check Belt because it needs tightened, or something is slowing or preventing platform movement.
	In the alarm state, the audible alarm is muted by pressing the Silence button, but rings back in about 30 minutes, if the alarm condition is still present. However, the message and indicator lights persist until the fault is corrected. Afterward, the remaining relevant alarm components are cleared by pressing the Silence button.

Sensor Fault	Sensor Fault alerts the operator that either of the shaker's two temperature sensors have failed, by an appropriate alarm message as shown.	Actual 08:41 250 37.0 Setpoints Main Temp Sensor Actual 08:41 250 37.0 Setpoints Over Temp Sensor
	In the alarm state, the audible alarm is m button, but rings back in about 15 minu present. However, the message and indic corrected. Afterward, the remaining relev by pressing the Silence button.	uted by pressing the Silence tes, if the alarm condition is still ator lights persist until the fault is rant alarm components are cleared
Temperature High or Low	Temperature High or Temperature Low alerts the operator that the operating temperature of the shaker has risen above or fallen below the programmed temperature tracking limit control point, by an appropriate alarm message as show	Actual 08:41 250 47.0 Setpoints Temperature is High Actual 08:41 250 27.0 Setpoints Temperature is Low
	In the alarm state, the audible alarm is m button, but rings back in about 15 minu present. However, the message and indic corrected. Afterward, the remaining relev by pressing the Silence button.	nuted by pressing the Silence tes, if the alarm condition is still ator lights persist until the fault is rant alarm components are cleared
Platform Stalled	Platform Stalled alerts the operator that free platform movement is inhibited. Th motor will automatically shut off and attempt to restart after approximately 15 audible alarm is muted by pressing the S about 15 minutes, if the alarm condition message and indicator lights persist until will continue to cycle on and off until th unit is turned off. On motor restart, the are automatically cleared, but the alarm to the Silence button	Actual 08:41 0 23.7 Platform Stalled seconds. In the alarm state, the ilence button, but rings back in is still present. However, the the fault is corrected. The motor e obstruction is removed, or the audible alarm and indicator lights nessage will remain until pressing
	Caution Turn off unit power when remo	ving any platform obstruction to

prevent possible injury.

Check Fuse Check Fuse alerts the operator that there Actual O8:41 0 23.7 is no communication with the motor Setpoints Check Fuse drive circuit board (historically the primary drive motor fuse has blown). In the alarm state, the audible alarm is muted by pressing the Silence button, but rings back in about 30 minutes, if the alarm condition is still present. However, the message and indicator lights persist until the fault is corrected. When the unit is turned on after fuse replacement, all alarm indicators are automatically cleared.

Warning Fuse replacement must be performed by qualified service personnel. See Service section. ▲

Change Configuration

To access the system Configuration menu, press the down arrow, the up arrow and the Silence button in that sequence.



Pressing the up arrow returns to the Operating Screen.

During the following configuration procedures, menu options are given to either modify a setting as it appears in sequence, or scroll past to the next item. If no selection is made by pressing a button or arrow, the display will revert to the Operating Screen in about 15 seconds. The complete configuration menu is shown in the chart at the end of this section.

Note In these procedures, values and settings for time, temperature, speeds, alarms, and so forth are shown on the display screens. These numbers are for example only and may not be the values encountered when programming your unit. ▲



Set Alarm Limits (continued)	The following screen appears and the present over-temperature alarm setting flashes. Actual Overtemp Alarm 64.4
	Change the temperature setting by pressing the up or down arrow. When set, press any of the three buttons (Time, Speed, Temp) to save the new setpoint and return to the previous screen or press nothing for about 15 seconds to save the new setpoint and return the display to the Operating Screen.
	Note The Overtemp Alarm setpoint values are calculated from the hardware and will not include every numerical value between the upper and lower limits. ▲
	When the overtemperature setpoint is exceeded by a few tenths of a degree, the control system will shut the shaker down by turning off the heaters.
	The Overtemp Shutdown warning shown above will be displayed, the indicator lights will flash and the audible tone (if not turned off) will sound.
	In the alarm state, the audible alarm is muted by pressing the Silence button, but rings back in about 15 minutes, if the alarm condition is still present. However, the message and indicator lights persist until the fault is corrected. Afterward, the remaining relevant alarm components are cleared by pressing the Silence button.
Set the Temperature Alarm Tracking Limit	The Temperature Tracking alarm activates whenever the operating temperature goes above or below the setpoint temperature by a user selectable value in the range of 1°C to 20°C. The limit is set at the factory as 10°C above and below the temperature setpoint.
	Note The above and below limits will always be the same value. \blacktriangle
	To change this limit, open the Configuration menu by pressing the down arrow, the up arrow, and the Silence button, then the down arrow again, in the sequence shown at the right.
	The screen shown below appears on the display: Actual Actual For more Setpoints Audible Alarms
	Then press the Temperature button beneath Alarms.

Section 2 Operation

Set Temperature Alarm Tracking Limit (continued)	From the screen below, press the Temperature button beneath Tracking (Trckng).
	The following screen will appear and the present Temperature Tracking alarm limit Setpoints 10.0
	Change the Temperature Tracking limit by pressing the up or down arrow. When set, press any of the three buttons (Time, Speed, Temp) to save the new setpoint and return to the previous screen or press nothing for about 15 seconds to save the new setpoint and return the display to the Operating Screen.
	When the chamber temperature rises above or falls below the temperature tracking limit, the appropriate message is displayed (at right), the indicator lights flash and the audible tone sounds (if not turned off).
	In the alarm state, the audible alarm is muted by pressing the Silence button, but rings back in about 15 minutes, if the alarm condition is still

button, but rings back in about 15 minutes, if the alarm condition is still present. However, the message and indicator lights persist until the fault is corrected. Afterward, the remaining relevant alarm components are cleared by pressing the Silence button.

Remote Alarm System

Most of the alarm states described previously (see Table 2-2) can alert a remote alarm monitoring system through an internal SPDT relay connected to an RJ-11 jack on the rear of the shaker cabinet. Refer also to 'Connect the Remote Alarm' in Section 1. For the convenience of the laboratory, these remote alarms can be individually turned on or off. Any of the remote alarms set to On will activate the internal relay.

Note The remote Overtemp Shutdown, Platform Stalled and Check Fuse alarms cannot be deactivated. ▲

To set the remote alarms to On or Off, open the Configuration menu by pressing the down arrow, up arrow, and the Silence button, and then the down arrow three times, in the sequence shown at right. The screen shown at right will appear on the display.



Press the Speed button beneath Remote (Rmte). The alarms will be shown in the following sequence.

Cycle Complete Toggle the Cycle Complete alarm with either the up (On) arrow or the down (Off) arrow. Pressing Temp

button beneath Next advances the display to the next alarm, saving the shown Cycle Complete setting to memory.



If no buttons are pressed, the display automatically returns to the Operating Screen after about 15 seconds, saving the selection to memory.

 Power Failure
 Toggle the Power Failure alarm with either the up (On) arrow or the down (Off) arrow. Pressing the Temp button beneath Next advances the display to the next alarm, saving the Power Failure setting to memory.
 Power Failure:
 On Next

If no buttons are pressed, the display automatically returns to the Operating Screen after about fifteen seconds, saving the selection to memory.

Section 2 Operation

RPM Tracking	Toggle the RPM Tracking alarm with either the up (On) arrow or the down (Off) arrow. Pressing the Temp button beneath Next advances the display to the next alarm, saving the RPM Tracking setting to memory.
	If no buttons are pressed, the display automatically returns to the Operating Screen after about 15 seconds, saving the selection to memory.
Check Belt	Toggle the Check Belt alarm with either the up (on) arrow or the down (off) arrow. Pressing the Temp button beneath Next advances the display to the next alarm, saving the Check Belt setting to memory.
	If no buttons are pressed, the display automatically returns to the Operating Screen after about 15 seconds, saving the selection to memory.
Sensor Fault	Toggle the Sensor Fault alarm with either the up (On) arrow or the down Setpoints Sensor Fault: On Next (Off) arrow. Pressing the Temp button beneath Next advances the display to the next alarm, saving the on/off setting to memory.
	If no buttons are pressed, the display automatically returns to the Operating Screen after about 15 seconds, saving the selection to memory.
Temperature High or Low	Toggle the Temperature High or Low alarm with either the up (On) arrow or the down (Off) arrow. Pressing the Temp button beneath Return returns the display to the previous screen, saving the on/off setting to memory.
	If no buttons are pressed, the display automatically returns to the Operating Screen after about 15 seconds, saving the selection to memory.

View Total Operating Hours

Whether the unit has been operated in Hold or Countdown modes, and/or has been turned off and unplugged many times, the microprocessor control system maintains a running total platform operating hours.

To view this information, open the Configuration menu by pressing the down arrow, up arrow and Silence button, then the down arrow three times, in the sequence shown at right.

The screen shown at right will appear on the display:

Pressing the Time button beneath RunHrs shows total accumulated run hours as displayed in the illustration at the right. When finished, press any of the three buttons (Time, Speed, or Temp) to save the new setpoint value and



return to the previous screen, or press nothing for about 15 seconds to save the new setpoint value and return to the Operating Screen.

Heat % Heat percent is intended for factory use only, but can be helpful in troubleshooting the heat control system.

To view this information, open the Configuration menu by pressing the down arrow, up arrow and Silence button then the down arrow three times, in the sequence shown at right. The screen shown at right will appear on the display:

Press the Temp button beneath Heat %.

Main Heat % is the percentage of time that the chamber heater is turned on during a five second period.



Example: If the heater is being cycled on for two seconds and off for three seconds, the Heat % value is 40 percent.

When finished, press any of the three buttons (Time, Speed, or Temp) to save the new setpoint value and return to the previous screen, or press nothing for about 15 seconds to save the new setpoint value and return to the Operating Screen.

Software Version

Software Version is for factory use only and will be important if troubleshooting the microprocessor programming is ever necessary.

To view this information, open the Configuration menu by pressing the down arrow, up arrow and Silence button, then the down arrow four times, in



the sequence shown at right. The screen below will appear on the display:



Software Version *#* XXXXXXX

Press the Time button beneath SwVers and the screen at above right will appear, showing the Model SHKE435HP/SHKE436HP software version in the control system memory.

To return to the previous screen, press the Time button. To return to the Operating Screen, wait about 15 seconds.

Overtemp Sensor Readings

Overtemp Sensor Readings is for factory use only and will be important if troubleshooting the microprocessor programming is ever necessary.

To access this screen, press the down arrow, up arrow, Silence button, then the down arrow button four more times.

The screen shown at the right will appear on the display.

Press the Speed button beneath O-Temp and the screen at the right will appear, showing the temperatures being read by the Overtemperature sensor.



Actual Over Temp. 37.2 Setpoints

When finished, press any of the three buttons (Time, Speed, or Temp) to save the new setpoint value and return to the previous screen, or press nothing for about 15 seconds to save the new setpoint value and return to the Operating Screen.
Selecting Hold or Countdown Time Setting Operating Speed Setting Operating Temperature

Orbital Shaker Menu Map



Orbital Shaker Menu Map

Turning the Audible Alarm On and Off



Calibrating Speed Calibrating Temperature Viewing Total Unit's Running Time Viewing Percent Heat

Orbital Shaker Menu Map



Setting Overtemperature Alarm Value Setting Temperature Tracking Limit Value

Orbital Shaker Menu Map



Page Four

Turning the Individual Remote Alarms On and Off

Orbital Shaker Menu Map



Note: Numerical values and alarm settings shown here are for reference only and may not match any specific shaker

Viewing Software Version Viewing Overtemperature Sensor Reading

Orbital Shaker Menu Map



Section 3 Maintenance

	The Model SHKE435HP and SHKE436HP shakers use a brushless DC motor and oversized, permanently-lubricated bearings, requiring no maintenance.	
Gas Springs	The gas springs should be checked periodically, and ideally every six months. The opening force, as measured from the front lip from a closed position, should be below 100 N (22.5 lbf) maximum. If the force is above this value, the gas springs should be replaced. If a force measurement is not possible, the gas springs should be replaced every two years.	
Platform and Cabinet Cleaning	The anodized brushed aluminum platform and powder-coated steel cabinet surfaces can be cleaned with common laboratory materials. However, liquids should not be allowed to enter the shaker cabinet from under the platform. All spills should be cleaned up immediately. If necessary, remove the platform. Follow 'Installing the Platform' procedure in Section 1 when re- installing the platform.	
Control Panel	The control panel has sealed push buttons and a liquid crystal display. It may be cleaned with a mild detergent and dried with a soft cloth.	
Cleaning or Replacing Air Filter	The air filter is located behind the grille on the front of the cabinet. The grille is held in place by six press-in type retainers and is easily removed by grasping it by the edges and pulling it off.	
	The air filter is held in place by four retaining springs (Figure 3-1) and is easily removed. It may be washed in water with a mild detergent and dried between two lint-free towels.	
	Retaining Springs	



Figure 3-1. Air Filter

PREVENTIVE MAINTENANCE
Shakers
Your equipment has been thoroughly tested and calibrated before shipment. Regular preventive maintenance is important to keep your unit functioning properly. The operator should perform routine cleaning and maintenance on a regular basis. For maximum performance and efficiency, it is recommended the unit be checked and calibrated periodically by a qualified service technician.
The following is a condensed list of preventive maintenance requirements. See the specified section of the operating manual for further details.
We have qualified service technicians, using NIST traceable instruments, available in many areas. For more information on Preventive Maintenance or Extended Warranties, please contact us at the number below.
Cleaning and calibration adjustment intervals are dependent upon use, environmental conditions and accuracy required.
Tips for all shakers:Use only our standard flat-head screws for flask clips.Use only our standard round-head screws for test tube racks, holders and utility trays.

401 Millcreek Road, Box 649 • Marietta, Ohio 45750 USA • 740-373-4763 USA and Canada 800-438-4851 • Telefax: 740-373-4189 • http://www.service.led.marietta@thermofisher.com

efer to Manual Section	Action	Daily	Monthly	Yea
ł	Clean the unit with mild detergent and wipe dry as needed		~	

~

Refer to Manual Section	Action	Daily	Monthly	Yearl
-	Clean the unit with mild detergent and wipe dry as needed		>	
	Clean the window with a mild detergent and wipe dry		^	
	Check under the platform for broken glass or other debris.		^	
4	Inspect air filter. Clean as needed		^	
£	* Check and document calibration of temperature, alarms, speed and time, as applicable			>
	* Verify operation of circulation fan motor			>
2	Change the HEPA filter, as needed			>

* Qualified service technicians only

Preventive Maintenance for SHKE435HP Series Shakers

Section 4 Service

Caution The procedures outlined in this section must be performed by persons experienced in servicing and maintaining laboratory equipment. Lockout and tagout electrical power connections whenever removing cabinet panels or working on electrical or motor control components. To avoid damage to solid state electrical components, proper grounding techniques must be observed whenever working on this shaker. ▲

With the exception of the chamber HEPA filter, the Model SHKE435HP and SHKE436HP Orbital Shakers contain no user-serviceable components.

Alarms and Alarm Conditions

If the microprocessor control system senses a fault, malfunction or abnormal operating condition, alarm messages appear on the display. These messages will be helpful should service or repair assistance be necessary. Refer to the table below and the alarm matrix at the end of this section.

Table 4-1. Alarms

Alarm Message	Fault Condition
Overtemp Shutdown	System shutdown due to overtemperature condition
Main Temp Sensor	Temperature sensor has failed
Over Temp Sensor	Temperature sensor has failed
Temperature is High	Temperature tracking has sensed higher temperature than setting
Temperature is Low	Temperature tracking has sensed lower temperature than setting
RPM is High	RPM tracking has sensed shaker speed higher than setting
RPM is Low	RPM tracking has sensed shaker speed lower than setting
Power Failure	Power has failed during shaker operation or shaker power is off for more than 1.5 hours
Cycle Complete	Countdown to zero time has been reached. Unit stops.
Check Belt	Motor V-belt has broken or slipped
Audible is Disabled!	Continuously notifies operator that audible alarm has been disabled
Platform Stalled	Free movement of the platform has been obstructed
Check Fuse	Power loss to motor drive circuit board, most likely the primary drive motor fuse has blown

Change HEPA Filter

The HEPA filter is located on the left side of the chamber, and is accessed by pulling up the four black press-in fasteners and sliding the cover off. Refer to Figure 4-1.



Figure 4-1. HEPA Filter and Temperature Sensors Locations

If the Shaker Will Not Operate

If the shaker platform won't operate with the unit plugged in and the power switch turned on, the following conditions may be present:

- The lid is open. Lower the lid to its fully closed position.
- Time countdown has been reached. Reset the time, or change to continuous operation (Hold).

Spare Fuses

S Three spare fuses are provided with this shaker and are taped to the underside of the control panel plastic frame. The plastic frame is attached to the cabinet by Velcro strips. Grasp the frame by the corners and pull to remove. There are also small indents located along the edges of the panel to accommodate a flat screwdriver blade. Figure 4-2 illustrates the underside of the frame.



Figure 4-2. Spare Fuse Location

Spare Fuses (continued)

Three fuse holders are located on the left side of the relay tray located in the lower part of the console cabinet. Figure 4-3 shows the location of the fuses. Refer to Table 4-2 for a list of their electrical ratings, part numbers, and applications.





Access to the relay tray is made by removing the grille from the front of the cabinet. It is held in place by six press-in type retainers and is easily removed by grasping the edges of the grille and pulling it off.

To remove the panel beneath the grille, remove six Phillips screws; three on the bottom of the panel and three on the top. The two Phillips screws on the left side of the foot pedal will need to be loosened to allow the panel to slide outward.

Fuses, SHKE435HP		
Rating	Application	Part Number
0.25 amp	Main Power Relay Board	230144
1.6 amp	Drive Motor	230145
0.1 amp	Recorder (opt.)	30107

Fuses, SHKE436HP		
Rating	Application	Part Number
0.15 amp	Main Power Relay Board	230142
0.8 amp	Drive Motor	230141
0.1 amp	Optional Recorder	230107

Caution Do not substitute! Replace these fuses with fuses of identical electrical ratings only. ▲

Circuit Boards Warning Only qualified service personnel should perform this procedure.

Four circuit boards control the Orbital Shaker. Three boards are located in the relay tray compartment, the fourth is behind the LCD display. Figure 4-4 identifies the circuit boards and other major components in the relay tray. Refer also to the relay tray wiring diagrams in Section 7.

To access the panel, unplug the shaker and move it to a sturdy location that will allow the back of the cabinet to swing down and lie flat. Remove the screws indicated by the arrows in Figure 4-2 and lower the back panel.

Components in the electronics panel are identified in Figure 4-4. Refer to the parts list and the electrical schematics in the back of this manual for part numbers.



Figure 4-4. Electrical Component Locations, Relay Tray Compartment

Temperature Sensors

Two temperature sensors are located behind a perforated cover plate on the right side of the chamber. Refer to Figures 4-1 and 4-5. To access these sensors, pull outward on the two black press-in fasteners on the top edge of the cover and lift the cover upward. The cover is held in place with four metal clips.



Figure 4-5. Chamber Right Side with Perforated Cover Plate Removed

To replace the cover, make sure all four clips engage the metal edges of the chamber and the two fasteners are firmly seated in their holes. Press the top of the fastener in until a "click" is heard. See Figure 4-6.



Figure 4-6. Fastener

Blower Fan Motors and Heating Elements

Locations of the Blower Motor and Ambient Fan Motor are shown in Figure 4-7. Removal of the relay tray is necessary to service the Blower Motor or the Heating Elements. Access to the heaters is through an access port on the front of the air plenum.



Figure 4-7. Air Plenum with Heating and Air Moving Components

Heater Element Circuit Breaker

Warning Only qualified service personnel should perform this procedure.

Warning Remove and lock-out electrical power when working on or near the relay control tray and heating element connectors. Allow sufficient time for the heating elements to cool before Heating Element reaching into that area. \blacktriangle

A manual reset circuit breaker is located between the heating element electrical connectors on the side of the air plenum. (Figures 4-7 and 4-8) The breaker can be reset by removing the front grille and reaching over the relay tray.



Figure 4-8. Connectors and Reset

Tune the Cabinet After the console cabinet is in place and leveled, with the platform installed, turn the unit on and set the speed to 300RPM. Kneeling in front of the console, lightly touch the lower left and right corners of the cabinet. If one side seems to vibrate more than the other, raise or lower the corner support leg using the 3/4" open end wrench supplied in the parts bag. Continue this "fine tuning" until the vibrations are reduced as much as possible.

Service the Drive Belt



Figure 4-9. Front View of Cabinet with Grille Removed

The motor drive pulley, large mechanism pulley, belt and motor mounting bolts are visible after removing the grille and belt/pulley tray (Figures 4-9 and 4-10).

Loosening the three 7/16" bolts at the base Figure 4-10. Drive Belt of the drive motor allows the belt to be changed or tension applied to the belt.

Tighten the three bolts and torque to 10 ft. lbs.



Service the Drive Belt (continued)

To remove the pulleys from their shafts, use a 1/8" Allen wrench to remove two set screws from the belt groove of the larger pulley; use a 5/16" Allen wrench to remove the single set screw from the belt groove of the smaller pulley.

When replacing the pulleys, seat the larger pulley completely against its baseplate. The smaller pulley, however, is installed with 0.300" space between it and the baseplate.

Calibrate Speed (RPM)

An external calibrated speed measuring device can be used to adjust the actual platform speed so that the unit setpoint speed matches the external device measurement. Calibration of the platform speed is performed at the factory at 250 RPM. To change the actual platform speed, open the Configuration menu by pressing the down arrow, up arrow, and the

Silence button, and then the down arrow twice, in the sequence shown at right.

The screen at right appears on the display.



Then press the Speed button beneath RPM.

The value shown on this screen is the present Speed setpoint. Using the up

and down arrows, increase or decrease the platform speed until the reading on an independent, accurate speed measuring device matches the Speed setpoint.



When finished, press any of the three buttons (Time, Speed, or Temp) to save the new setpoint value and return to the previous screen, or press nothing for about 15 seconds to save the new setpoint value and return to the Operating Screen.

Calibrate the Temperature

- 1. Place a 250ml Erlenmeyer flash (filled with approximately 100ml of liquid) in approximately the geometric center of the platform.
- 2. Suspend an independent temperature measuring device of known accuracy into the flask. The sensor should be submerged in the liquid but not in contact with the bottom or sides of the flask.
- 3. Adjust the shaker temp setpoint at desired calibration temperature.
- 4. Set the shaking speed setpoint to 75 RPM.
- 5. Start the unit and allow a minimum of 2 hours stabilization of cabinet and flask liquid.
- 6. Enter Calibration mode by pressing the down arrow, the up arrow, the Silence button. Then press the down arrow twice. The screen at right will appear on the display:
- 7. Press the Temperature button Setpoints beneath Temp.
- 8. Using the up and down arrows, increase or decrease the temperature value to match the independent, accurate temperature measuring device.



- 9. When complete, press the Time, Speed, or Temp button to save the setting. The display will return to the Calibrate RPM Temp screen. (Or, if nothing is pressed for about fifteen seconds, the display will return to the Operating Screen and the setting will be automatically saved to memory.)
- 10. Allow the cabinet to re-stabilize for 1 hour. Recheck temperature. If necessary, return to Step 6 until no additional adjustments are needed.
- 11. Temperature calibration is now complete. Remove calibration equipment and resume use.

Alarm Message	Alarm Criteria	Alarm Delay*	Alarm Ringback*	System State	Corrective Action
Over Temp Shutdown	Temperature at the over temp sensor is about 1° over shut down set point	None	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters off	Press SILENCE to silence audible alarm Check for air intake blockage Over temperature probe malfunction Sensor connector unplugged Heater circuit not cycling Main circuit board failure Call Forma's Service Department
Main Temp Sensor	Sensor circuit is open or shorted beyond the expected resistance range in either direction	30 sec.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters off	Press SILENCE to silence audible alarm Check board connector Check sensor circuit Replace sensor Call Forma's Service Department
Over Temp Sensor	Sensor circuit is open or shorted beyond the expected resistance range in either direction.	30 sec.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters on	Press SILENCE to silence audible alarm Check board connector Check sensor circuit Replace sensor Call Forma's Service Department
Temperature is High	Temperature is above control set point by temperature tracking limit	None	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters on	Press SILENCE to silence audible alarm Check temperature tracking limit Check sensor circuit Replace main temperature sensor Call Forma's Service Department
Temperature is Low	Temperature is below control set point by temperature tracking limit	30 min.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters on	Press SILENCE to silence audible alarm Check if lid is completely closed Check temperature tracking limit Check sensor circuit Replace main temperature sensor Call Forma's Service Department
	Electrical power has been disrupted	Upon power up	None	Not affected	Warning notice only Press SILENCE to silence audible alarm

* Alarm Delay and Ringback times are approximate

CHART.cdr

Alarm Message	Alarm Criteria	Alarm Delay*	Alarm Ringback*	System State	Corrective Action
Cycle Complete	Count-down time has reached zero	None	None	Alarm light on Audible alarm on Blower fans on Shaker motor off Heaters on	Advisory notice only Press SILENCE to silence alarm
RPM High	RPM is above control set point by tracking limit	2 min.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters on	Press SILENCE to silence audible alarm Check platform loading Check RPM tracking limit setting Shut the unit off and call Forma's Service Department
RPM Low	RPM is below control set point by tracking limit	2 min.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters on	Press SILENCE to silence audible alarm Check for overloaded platform Check for obstruction to edges of platform Check for low input AC mains voltage. Shut the unit off and call Forma's Service Department
Check Belt	Rotation sensor circuit sees no mechanical rotation or excessive belt slippage	None	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on Heaters on	Press SILENCE to silence audible alarm Shut the unit off and check the belt If alarm persists, call Forma's Service Department
Audible is disabled!	Operator has turned off the audible alarm	None	None	Normal operation	Lower half of the LCD display will show this warning as long as the audible alarm remains turned off
Platform Stalled	Motor tries to start but platform is obstructed	15 sec.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor on/off/on Heaters on	Press SILENCE to silence audible alarm Check for overloaded platform Check for platform edge obstruction Shut the unit off and check the belt If alarm persists, call Forma's Service Department
Check Fuse Power Failure	Primary drive motor fuse blown	15 sec.	15 min.	Alarm light on Audible alarm on Blower fans on Shaker motor off Heaters on	Press SILENCE to silence audible alarm Check/replace drive motor fuse If alarm persists, call Forma's Service Department

* Alarm Delay and Ringback times are approximate

Section 5 Specifications

Shaking

Range 25-525 RPM	
Accuracy±1 RPM	
Motion One inch/orbital	
IndicatorLCD in 1 RPM increments	
Temperature	
Range . 5° C (41°F) above ambient to 60° C (140°F)	
Control±0.1°C	
Uniformity±0.2°C (in flask)	
IndicatorLCD in 0.1°C increments	

Timer

Range . . Programmable from 5 minutes to 200 hours, or continuous operation

Indicator LCD in 5 minute increments

Run Time . . LCD counts down for a timed run or counts up in a "Hold" function in 1 minute increments/decrements

Alarms

Temperature . . Adjustable tracking high/low temp Speed Adjustable tracking high/low RPM Time Cycle complete Power Failure Loss of input power

Safety

Platform Speed Software independent speed control circuit

Platform Stall Software independent motor overcurrent protection circuit

LCD (Liquid Crystal Display)

Top line displays actual elapsed run time, speed, and temperature. Bottom line displays user time, speed and temperture setpoints alternating with any active alarm messages.

Drive

Triple counterbalanced. Compensates for unbalanced platform loads

Drive Motor

1/3 HP brushless DC, permanently-lubricated ball bearing

Lid

Counterbalanced, hand or foot operated, with tempered thermal pane window and key lock.

Automatic Restart

Microprocessor retains all programming in non-volatile memory. In the event a power outage, the shaker restarts automatically.

Construction

Interior Stainless steel with coved corners

Exterior Cold rolled steel

Finish Powder coated for a durable, easily maintained surface

Platform Anodized brushed aluminum

Dimensions

Exterior 45.0" W x 39.0" H x 30.0" F-B

Exterior (lid open) . 45.0" W x 39.0" H x 30.5" F-B

Interior 34.3" W x 18.8" H x 21.1" F-B

Electrical

SHKE435HP

Nominal: 120VAC, 60Hz, 1 PH, 8.6 FLA

SHKE436HP

Nominal: 230VAC, 50/60Hz, 1 PH, 3.5 FLA

Remote Alarm Contacts . . Cycle Completion, Speed, Temperature and Power Failure Alarms as selected by user

Certifications

Declaration of Conformity available on request

Ambient Operating Conditions

Temperature $18^{\circ}C$ (64°F) to 40°C (104°F)

Humidity 80% RH at or below 31°C, decreasing linearly to 50% RH at 40°C

Capacity

Flasks From (91) 25ml up to (4) 6L

Weights

Optional Platforms

Size 29.5" x 18" (74.9cm x 45.7cm)

Clips 25ml, 50ml, 125ml, 250/300ml, 500ml

1L, 2L, 4L, 6L and 2800ml Fernbach sizes available

Racks . . Adjustable angle test tube holder with rack, 13-30mm

Filter

HEPA: Rated efficient at 0.3 microns

Size: 18" x 16" x 2" (45.7cm x 40.6cm x 5.1cm)

Intended Use

Orbital shakers are designed to provide increased aeration in a stable temperature environment

Unintended Use

1) Not intended for use in Class I or II applications as defined in 21 CFR

2) Not intended for mixtures of flammable materials

Sound Level

Not to exceed 85db

Safety Specifications

Indoor Use Only

Humidity . . 80% RH at or below 31°C, decreasing linearly to 50% RH at 40°C

Fluorinated Greenhouse Gases

Compliant with REGULATION (EU) No 517/2014 OF THE EURO-PEAN PARLIAMENT AND OF THE COUNCIL on fluorinated greenhouse gases.

This product contains foam blown with fluorinated greenhouse gas, R-245fa.

Mains Supply Fluctuations . . ±10%

Installation Category II¹

Pollution Degree 2²

Class of Equipment I

1 Installation category (overvoltage category) defines the level of transient overvoltage which the instrument is designed to withstand safely. It depends on the nature of the electricity supply and its overvoltage protection means. For example, in CAT II which is the category used for instruments in installations supplied from a supply comparable to public mains such as hospital and research laboratories and most industrial laboratories, the expected transient overvoltage is 2500V for a 230V supply and 1500V for a 120V supply.

2 Pollution Degree describes the amount of conductive pollution present in the operating evironment. Pollution Degree 2 assumes that normally only non-conductive pollution such as dust occurs with the exception of accasional conductivity caused by condensation.

Section 6 Parts List

*Refer to Section 4 for electrical fuse information.

SHKE435HP

Part No
129024 Pneumatic Spring, 80 lbs.
138009 Heater, Wirewound 450W 115V/230V
156089Motor, Brushless 24V
191535Motor Drive Board
191734Display/Keypad Replacement Kit
190525 Relay Board
192589 Control Board Replacement
290160 Probe, 2252 Ohm/25°C, 1/8 x 2 (2)
300275 Relay, DPDT 20A 120V
400113 Thermostat - Opens at 200°F
420064 Transformer, 130VA
420085 Transformer, 25VA
800040V-Belt, A x 43 1/2" x 45"
435051 Screwdriver, Phillips 8-3/4"
443020Wrench, 5/32" Hex with T-handle
194046 Spare Part Screw Bag, (for platform and clips)
900113 Tubeaxial Fan, 665 CFM 115V
900092 230 CFM Blower 115V 60Hz
760164 HEPA Filter
760167Air Filter 9.5" x 23.625"
230107 100mA Fuse (for optional recorder) T.D. 5mm x 20mm
230144
230145 1.6A Fuse T.D. 5mm x 20mm
443021
107003 Lid Glass Window
990024

SHKE436HP

Part No
129024Pneumatic Spring, 80 lbs.
138009 Heater, Wirewound 450W 115V/230V
156089Motor, Brushless 24V
191535
191545 Temp Control Board
191734 Display/Keypad Replacement Kit
192589 Control Board Replacement
290137 Probe, 2252 Ohm/25°C, 1/8 x 2 (2)
300276 Relay, DPDT 20A 240V
400113 Thermostat
420064 Transformer, 130VA
420085 Transformer, 25VA
800040V-Belt, A x 43 1/2" x 45"
435051 Screwdriver, Phillips 8-3/4"
443020Wrench, 5/32" Hex with T-handle
194046 Spare Part Screw Bag, (for platform and clips)
900149 Tubeaxial Fan 547 CFM 230V
900093 Blower 230 CFM 220V 50Hz
760164 HEPA Filter
760167Air Filter 9.5" x 23.625"
230107 100mA Fuse (for optional recorder) T.D. 5mm x 20mm
230144
230145 1.6A Fuse T.D. 5mm x 20mm
443021 ¾" Open End Wrench
107003 Lid Glass Window
990024







Section 7 Electrical Schematics





Section 7 Electrical Schematics

	WIDE #							
77	WIRE #	GALIGE		WIRE #	10			
/8	2	14	BLUE	25	18	BREWN		
79	3	14	GRN/YEL	33	24	RED		
80	3A-3C	18	GRN/YEL	34	24	BLACK		
81	3E-3K		SHIELDS	35	24	RED		
01	4	14	BLACK	36	24	BLACK		
82	4A	14	BLACK	37	24	BLACK		
83	5	14	WHITE	38	24	REU		
84	JA 5B	14	WHITE	29	24	WHITE		
05	6	14	BROWN	41	24	BROWN		
00	6A	14	BROWN	42	24	BLACK		
86	8	24	BLACK	43	24	RED		
87	9	24	PURPLE	44	24	GREEN		
88	10	18	BR⊡₩N	45	24	WHITE		
00	11	18	RED	46	24	GREEN		
89	12	18	ORANGE	47	24	BLACK		
90	13	18	YELLUW	48	24	RED		
91	14	18		49 50	24	PED		
	16	18	BLUE	51	24	GREEN		
92	17	18	RED	52	24	WHITE		
93	18	18	DRANGE	53	24	BROWN		
94	19	18	RED	54	18	RED		
95	22	18	YELLOW	55	18	WHITE		
95	23	18	BROWN	56	18	BLACK		
90	23A	18	BRUWN	57	24	BLACK		
97	24	24 24		58	24 19	BLACK		
98	25	24	RL ACK	594	18	BROWN		
00	27	24	BLACK	82A	18	PURPLE		
33	28	24	RED	86	14	BROWN		
100	29	24	BLACK	87	14	BLUE		
101	30	24	ORANGE					
102								
100								
104								
105								
106								
107								
107								
				2 DS-751	04-03-14 GLS GL	S CCS CHG MIC	RD BD FRDM 191542	LEIGCILICOL 7C
				1 DS-423	09-01-11 GSW SAU	DRP 900149		
				0 07-5/8		CUN RELEASE		
		THIS DOCU	IENT CONTAINS PROPR	REV ECN ND.	DATE BY CAU PART NAME: 436 C	D APPO DESCRIP DNSDLE INCLUBATOR SHA	TION OF REVISION	– 436 T Console Inc
AT	TENTION	THIS DOCU INFORMATION BE DISCLOSED USED FOR MA	MENT CONTAINS PROPR AND SUCH INFORMATION I O TO DTHERS FOR ANY PUR INUFACTURING PURPOSES		DATE BY CAU PART NAME: 436 C TLE: ELECTRICAL	CLUN RELEASE D APPD DESCRIP DNSDLE INCUBATOR SHA SCHEMATIC		– 436 Console Inc – Shaker





THERMO FISHER SCIENTIFIC DIGITAL SHAKER WARRANTY USA
The Warranty Period starts two weeks from the date your equipment is shipped from our facility. This allows shipping time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the warranty period.
During the first 24 months, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, <u>labor included</u> . For an additional 3 years, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, <u>labor excluded</u> . In addition, the Orbital Shaker mechanism is warranted for 10 years, parts only, F.O.B. factory. The mechanism is defined as the bearing assemblies. The warranty will be void if the equipment is altered without written authorization from Thermo. Installation and calibration is not covered by this warranty agreement. The Technical Services Department must be contacted for warranty determination and direction prior to performance of any repairs. Expendable items, i.e., glass, filters, light bulbs and lid gaskets are excluded from this warranty. Extended warranties are dependent on the units being maintained regularly as stated in the operation and service manuals.
Replacement or repair of components parts or equipment under this warranty shall not exceed the warranty to either the equipment or to the component part beyond the original warranty period. The Technical Services Department must give prior approval for return of any components or equipment.
THIS WARRANTY IS EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER WRITTEN, ORAL, OR IMPLIED. NO WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY. Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.
Your local Thermo Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction manuals carefully detail equipment installation, operation, and preventive maintenance.
If equipment service is required, please call your Technical Services Department at 1-800-438-4851 (USA and Canada) or 1-740-373-4763. We're ready to answer any questions on equipment warranty, operation, maintenance, service and special applications. Outside the USA, contact your local distributor for warranty information.
Rev. 2 6/2015

THERMO FISHER SCIENTIFIC INTERNATIONAL DIGITAL SHAKER WARRANTY
The Warranty Period starts two months from the date your equipment is shipped from our facility. This allows shipping time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the warranty period.
During the first 24 months, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, <u>including labor</u> . For an additional 3 years, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, <u>excluding labor</u> . In addition, the Orbital Shaker drive mechanism is warranted for 10 years, parts only, F.O.B. factory. The mechanism is defined as the bearing assemblies. The warranty will be void if the equipment is altered without the written authorization from Thermo. Installation and calibration is not covered by this warranty agreement. The local Thermo Fisher Scientific office must be contacted for warranty determination and direction prior to performance of any repairs. Expendable items, i.e., glass, filters, light bulbs and lid gaskets are excluded from this warranty. Extended warrants are dependent on the units being maintained regularly as stated in the operation and service manuals.
Replacement or repair of component parts or equipment under this warranty shall not exceed the warranty to either the equipment or to the component part beyond the original warranty period. The local Thermo Fisher Scientific office must give prior approval for return of any components or equipment.
THIS WARRANTY IS EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER WRITTEN, ORAL, OR IMPLIED. NO WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY. Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.
Thermo International Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction manuals carefully detail equipment installation, operation, and preventative maintenance.
If equipment service is required, please call your local Thermo Fisher Scientific office. We're ready to answer your questions on equipment warranty, operation, maintenance, service and special applications.
Rev. 2 6/2015 Rev. 2 6/2015

thermoscientific.com

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Thermo Fisher Scientific 401 Millcreek Road Marietta, Ohio 45750 United States

