



# Thermal Rocker®

Model No.

4637Q  
4637JPNQ  
4637-1CEQ  
4637-1CECN

---

# Table of Contents

Description .....	3
Safety Information .....	4
Alert Signals.....	4
Specifications .....	5
Power Requirements .....	5
Rocking Speed.....	5
Temperature Range .....	5
Overall Dimensions .....	5
Platform .....	5
Shipping Weight.....	5
Declaration of Conformity .....	6
Units Environmental Operating Conditions .....	6
Unpacking and Installation .....	7
Shipping Carton.....	7
Location .....	7
Electrical Requirements.....	7
Operation .....	9
Temperature Controlled Operations.....	9
Auto Tune .....	10
For Best Results .....	10
Auto Tuning Procedure .....	11
Temperature Calibration.....	11
Rocking Motion Only Applications .....	13
General Notes .....	13
Maintenance.....	14
Cleaning Spills .....	14
Parts Replacement .....	14
3-Position Rocker Switch.....	14
Rocker Motor.....	14
Motor Brushes.....	15
Temperature Controller .....	15
Changing Angle of Motion.....	16
Replacement Parts.....	17
Ordering Procedures .....	18
Warranty .....	19

---

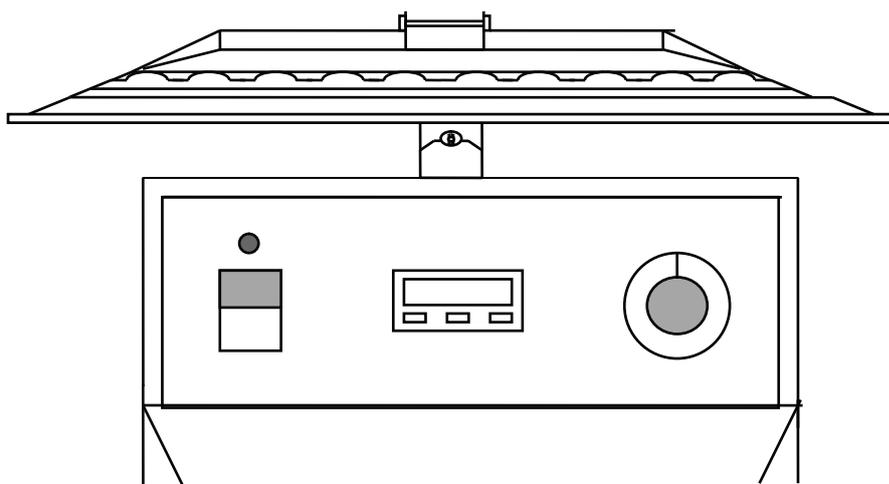
# Description

The Thermo Scientific Lab-Line Thermal Rocker provides the versatility of a variable speed rocking platform in a temperature-controlled environment. This makes it ideal for hybridizations, cell culture studies and blood procedures.

Rocking speed can be set from 0 to 100 cycles per minute and the angle of rotation is adjustable from 10° to 15° from horizontal. Microprocessor-based temperature controller maintains temperature under the polycarbonate cover from ambient +5°C to 70°C. Automatic resetting thermal cutoff terminates power to the heaters in the event temperature set point is exceeded.

The 14-3/16" square (36 cm) platform has a rubber mesh pad that prevents vessels from slipping and is removable for easy cleaning. A 1/2" (1.3 cm) ledge on all sides prevents vessels from sliding off. Rubber feet provide stability and keep the unit from creeping during operation.

The enclosed motor and rugged, all-steel, powder coated body are designed to provide years of virtually maintenance-free service. Convenient, panel mounted, switch allows operation with or without heat.



---

# Safety Information

---

## Alert Signals

**Warning**

Warnings alert you to a possibility of personal injury.

**Caution**

Cautions alert you to a possibility of damage to the equipment.

**Note**

Notes alert you to pertinent facts and conditions.

**Hot Surface**

Hot surfaces alert you to a possibility of personal injury if you come in contact with a surface during use or for a period of time after use.

**Risk of Electric Shock**

Disconnect unit from power source before removing cover.

Your Thermo Scientific Lab-Line Thermal Rocker® has been designed with function, reliability, and safety in mind. It is your responsibility to install it in conformance with local electrical codes. It is most important that the user follow installation instructions exactly as written. Failure to do so is likely to lead to improper operation, erroneous calibrations and possible damage to the equipment. Do not attempt operation without this information.

---

# Specifications

---

## Power Requirements

4637Q:	120 V, 50/60 Hz, 3.3 Amps, 400 Watts
4637JPNQ:	100-120 V, 50/60 Hz, 3.0-3.3 Amps, 400 Watts
4637-1CEQ & 4637-1CECN:	220-240 V, 50/60 Hz, 1.7 Amps, 400 Watts

---

## Rocking Speed

0 to 100 cycles per minute

---

## Temperature Range

From ambient +5°C to 70°C (with polycarbonate cover in place) as measured by a thermocouple positioned at approximate geometric center. The temperature in vessels with the polycarbonate cover in place will vary from these values depending on such factors as vessel size and contents, thickness of the glassware used and ambient conditions.

*Measurements conducted under controlled conditions with 30 cm dna tube filled with 10 ml water, ambient at 23°C and temperature equalizes at 42°C.*

---

## Overall Dimensions

16" W x 13-1/4" D x 7-3/4" H\* (41 x 34 x 20 cm)

---

## Platform

14-3/16" W x 14-3/16" D (36 x 36 cm)

---

## Shipping Weight

25 lb. (11 kg)

*\*Polycarbonate cover adds 5" (12.7 cm) to rocker's height*

---

## SPECIFICATIONS

---

### Declaration of Conformity

(For 220 to 240 volt, CE models only)

We hereby declare under our sole responsibility that this product conforms with the technical requirements of the following standards:

EMC:	EN 61000-3-2	Limits for harmonic current emissions
	EN 61000-3-3	Limits for voltage fluctuations and flicker
	EN 61326-1	Electrical equipment for measurement, control, and laboratory use; Part I: General Requirements
Safety:	EN 61010-1	Safety requirements for electrical equipment for measurement, control, and laboratory use; Part I: General Requirements
	EN 61010-2-010	Part II: Particular requirements for laboratory equipment for the heating of materials
	EN 61010-2-051	Part II: Particular requirements for laboratory equipment for mixing and stirring

*per the provisions of the Electromagnetic Compatibility Directive 2004/108/EC, and per the provisions of the Low Voltage Directive 2006/95/EC.*

*The authorized representative located within the European Community is:*

Thermo Fisher Scientific  
Electrothermal House  
Unit 12A  
Purdeys Industrial Estate  
Purdeys Way  
Rochford, Essex, SS4 1ND  
United Kingdom  
Tel: +44(0)1702 303350

Copies of the Declaration of Conformity are available upon request.

---

### Unit's Environmental Operating Conditions

Pollution Degree:	2
Installation Category:	II
Altitude:	2000 Meters MSL (Mean Sea Level)
Humidity:	80% maximum, non-condensing
Electrical Supply:	120VAC or 240VAC
Voltage Tolerance:	±10% of normal rated line
Temperature:	15°C to 40°C
Product Usage:	This product is intended for use indoors only

\*Refer to IEC 664-1

---

# Unpacking and Installation

---

## Shipping Carton

This should be inspected upon delivery. When received, carefully examine for any shipping damage before unpacking. If damage is discovered, the delivering carrier should both 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 is found promptly report it to the carrier and request a damage inspection promptly.

**IMPORTANT:** Failure to request an inspection of damage within a few days after receipt of shipment absolves the carrier from any liability for damage: you must call for a damage inspection promptly.

---

## Location

Place the unit on a level surface in a convenient location and near a grounded electrical outlet that matches the requirements listed on the nameplate. Locate the unit away from heat sources that might affect the temperature being maintained by a Thermal Rocker.

---

## Electrical Requirements

120 VAC models require a 120 VAC, 50/60 Hz power source. They are supplied with a 3-wire line cord. It should be plugged into an outlet designed for 3-prong plugs. If an extension cord is used, it also should be the 3-wire grounded type. For an outlet designed to accept 2-prong plugs (ungrounded), it is required that a qualified electrician replace the outlet with a new grounded type.

240 VAC models require a 240 VAC, 50/60 Hz power source and are supplied with a European cord set.

100-120 VAC models require a 100-120 VAC, 50/60 Hz power source. Because of the variety of plug configurations in use worldwide for 100 to 120 VAC power, the unit is furnished with the plug removed. The user must install a plug to conform with local code and configuration requirements.



### Caution

The power cord serves as the main power disconnect. Access to the power outlet should be unobstructed to allow the unit to be unplugged (disconnected) from the power source.



### Note

Leave rotator disconnected when not in use.

---

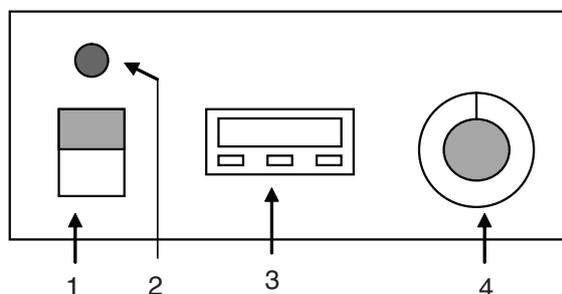
## UNPACKING AND INSTALLATION

If a plug must be installed, use only the 3-prong grounded type, rated for the unit load requirements and matching the power outlet. Make sure the green ground wire is secured to the plug ground terminal.

The user shall be made aware that the protection provided by the equipment may be impaired if the equipment is used with accessories not provided or recommended by the manufacturer, or used in a manner not specified by the manufacturer.

# Operation

Control Panel



1. **ROCKER SWITCH:** 3-position switch enables platform to: rock with heat, rock without heat, or heat only without motion.
2. **STATUS LAMP:** Illuminates when rocker's heaters are activated.
3. **TEMPERATURE CONTROLLER:** Microprocessor-based controller maintains temperature and provides display of either set point or actual temperature.
4. **SPEED CONTROL:** Adjusts rocking speed. Markings on dial are for reference purposes only.



## Warning

Do not use in the presence of flammable or combustible materials or explosive gases. Do not use in the presence of pressurized or sealed containers—fire or explosion may result, causing death or severe injury.



## Warning

The outer edges and bottom of the platform can become very hot at elevated temperatures. To avoid burns do not touch or get too close to the platform when heating at higher temperatures. Allow unit to cool down before removing vessels or touching unit.

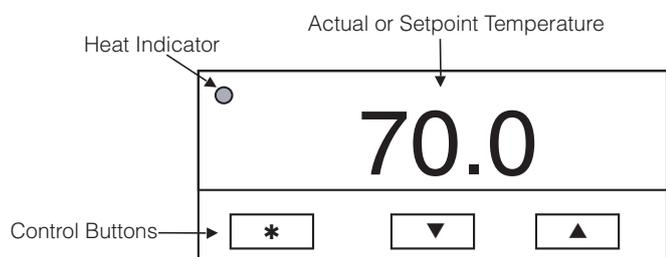
---

## Temperature Controlled Applications

- The polycarbonate cover must be in place for those applications where temperature control and rocking action are combined.
- Make certain that the 3-position rocker switch is in **OFF** position.
- Connect power cord to electrical outlet meeting requirements stated on unit's nameplate.
- Load platform with vessels to be shaken. Allow ample space between vessels so that accidental contact will be avoided.

The user is responsible for determining safe operation in the use of vessels at various rocking motion speeds and for checking that the vessels are secure while in motion.

- Turn the rocker switch to the **ROCK N HEAT** position to rock and heat simultaneously. Status light over switch will illuminate indicating power is supplied to the heaters. Adjust the temperature and speed control dials to desired settings.



1. **Controller Self-Test:** When the rocker is powered up, the controller will display **8888** along with the three decimal points and the heat **ON** indicator lamp. The display will then blank out for 2 seconds before showing the temperature.
2. **Heat **On** Indicator:** The heat **ON** indicator lamp is lit when the rocker's heaters are receiving power. The lamp will normally flash when the temperature is at set point.
3. **Set Point Adjustments:** The controller normally displays the temperature. To view or change the temperature set point proceed as follows:

Press	Controller
*	View setpoint
*▼	Decrease setpoint
*▲	Increase setpoint

- A. Press and hold the star key and use either the up or down arrow key to adjust the set point to the desired temperature. Release the star key.
- B. Allow at least 30 minutes for the chamber temperature to stabilize.

## Auto Tune

The auto tune program automatically adjusts the controller's parameters to achieve optimal temperature control. It is not necessary to run the auto tune program when setting up the rocker. However, if the temperature appears to be unstable, the auto tune program can be run using the procedure shown below:

### For Best Results

- Set the usual set point temperature and use normal load conditions.
- Allow the rocker to stabilize at set point for at least 30 minutes.

---

## Auto Tuning Procedure

- A. Enter the program mode by pressing and holding BOTH the up and down arrow keys for 3 seconds.
- B. Release BOTH arrow keys when **tunE** is displayed.
- C. The controller display should now be alternating between **tunE** and **oFF**.
- D. Press and hold the "**STAR**" (\*) key. Press and release the up arrow key until **At.SP** is displayed. Release the "**STAR**" (\*) key.
- E. After one minute has elapsed, the controller display will begin to alternate between showing the **chamber temperature, tunE** and **At.SP**.
- F. Allow the program to run until the display again shows only the temperature.

---

## Temperature Calibration

- A. Place a calibrated thermometer near the approximate geometric center of the platform in a position that would allow it to be read through the glass door.
- B. Press and hold the "**STAR**" (\*) key and using the up or down arrow key, adjust the set point to the desired temperature.
- C. Allow the unit to run for at least 30 minutes.
- D. Controller display should now be indicating the set point temperature. Make note of the thermometer reading.
- E. Press and hold both arrow keys until the controller display indicates **tunE**. Release the arrow keys. Press and release the down arrow key, the

display should now indicate **LEUL**. Press and hold the "**STAR**" (\*) key and using the up arrow key adjust the display to read **3**. Release the "**STAR**" (\*) key. Press and release the up arrow key until the display indicates **Zero**. The display should now alternate between **Zero** and a numerical value.

- F. Using the examples shown below and the thermocouple value obtained in step above, enter the correct **Zero** value into the controller by pressing the "**STAR**" (\*) key and using the up or down arrow keys. If there is already a **Zero** value present then add the new value to the one already present.

Thermometer	= 70 °C	Thermometer	= 70 °C
Controller Reading	= 75 °C	Controller Reading	= 65 °C
Subtract	= -5 °C	Subtract	= +5 °C
Enter <b>Zero</b> value of	-5 °C	Enter <b>Zero</b> value of	+5 °C

- G. When the correct **Zero** value has been entered, press and hold the two arrow keys together until the display again indicates the chamber temperature. If the procedure was done correctly, the controller display should now agree with the thermometer reading to within  $\pm 0.5^{\circ}\text{C}$ .
- H. Allow the unit to run for at least 30 minutes.
- I. Re-check the thermometer reading, the controller display and the thermometer should agree to within  $\pm 0.5^{\circ}\text{C}$ . If not repeat steps D, E and F above.

---

## Rocking Motion Only Applications

- Turn the rocker switch to the **ROCK** position for rocking action only.
- Adjust the speed control to the desired rocking speed.
- Load platform with vessels to be shaken. Allow ample space between vessels, so that accidental contact will be avoided.

The user is responsible for determining safe operation in the use of vessels at various rocking motion speeds and for checking that the vessels are secure while in motion.

---

## General Notes

During rocking action, observe vessels in motion to determine that contents are not escaping or spilling onto the platform. If this should occur, turn temperature and speed control dials to **0** and the 3-position rocker switch to the **OFF** position. Wait for platform to cool down before attempting clean up.

In the event of electrical overload, the unit is protected by resettable, pushbutton circuit breaker on the back. If overload occurs, the breaker will trip. To reset, disconnect rocker from power source and push button in.

However, if breaker trips repeatedly, unplug the unit and have the shaker serviced.

It is best to leave unit disconnected from power source when not in use.

---

# Maintenance



## Note

Make no attempt to service or repair a Thermo Scientific product under warranty before consulting your distributor. After the warranty period, such consultation is still advised, especially when the repair may be technically sophisticated or difficult.

If assistance is needed beyond what the distributor can provide, please call the customer service department at 1-800-553-0039. No merchandise, however should be returned without prior approval.



## Caution

Disconnect plug from electrical outlet before attempting any maintenance or repair of this unit.

---

## Cleaning Spills

In the event of any spills on the surfaces of the unit, carry out appropriate neutralization procedures and/or wash off with a cleaning solution of warm water and a mild detergent. Rinse off with clean water and dry thoroughly. Do not use abrasive cleaners on surfaces.

---

## Parts Replacement

---

### 3-Position Rocker Switch

To access the rocker switch, remove the cover. Turn the unit over and remove 6 Phillips screws holding the bottom. Remove the 5 wires attached to the switch. Note attachment sequence and numbers imprinted on switch terminal. Compress spring release on sides of switch and withdraw through the front of the panel. Insert the replacement switch. Reconnect wires to the new switch in the correct locations.

Reverse procedure and reattach the bottom cover.

---

### Rocker Motor

- To access rocker motor, remove bottom plate.
- Remove Allen screw from the end of the eccentric. Loosen the 2 set screws holding the eccentric to the motor shaft.
- Slide the eccentric off the end of the motor shaft. Remove the 4 screws holding the motor to bracket.
- Disconnect wires to the motor.
- Install new motor and reverse the procedure.

---

## Motor Brushes

- Under normal conditions of use, the DC motor brushes should be checked once a year and replaced if worn.
- Disconnect plug from outlet. Remove bottom plate.
- Unscrew brush cap from side of motor.
- Remove spring and brush. Replace brushes if worn to less than 1/4" (0.64 cm) in length. It is important that BOTH brushes be replaced, even though only one may be worn.

---

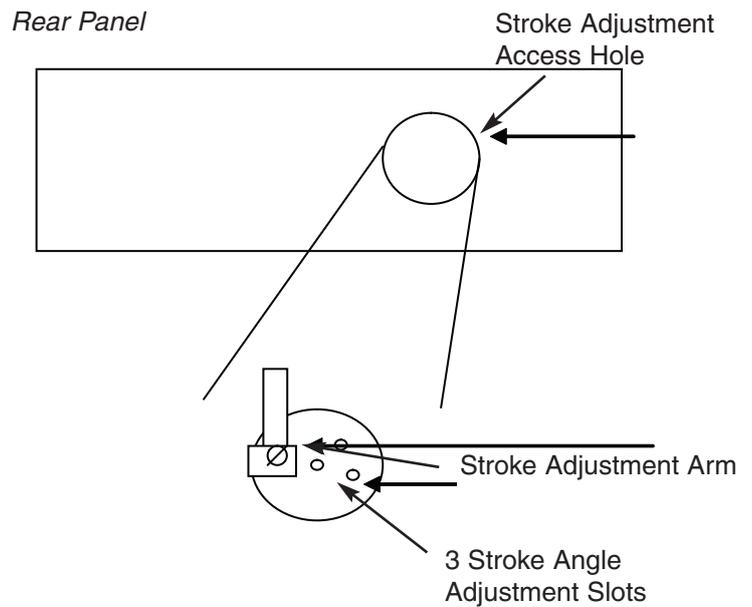
## Temperature Controller

1. Place ON/OFF switch in **OFF** position.
2. Unplug rocker from outlet power supply.
3. To remove controller from control housing:
  - use both hands to firmly grip each side of the controller bezel
  - press on the bezel side grips until the bezel tabs release
  - slowly pull controller from housing
4. To install new, factory configured controller:
  - Carefully slide new controller into controller housing.
  - Press controller bezel into controller housing until bezel tabs securely lock controller into place
5. Plug rocker into outlet power supply.
6. Place ON/OFF switch in **ON** position.

**Note**

PCB contacts at rear of controller fit into contacts at rear of controller housing.

## Changing Angle of Motion



1. Pry off plug from stroke adjustment accesses hole located on the right side of rocker's rear panel.
2. Insert screwdriver and remove the stroke adjustment screw that holds stroke adjustment arm in place.
3. Reposition stroke adjustment arm over one of the 2 remaining stroke adjustment slots.
4. Tighten into position with screw removed in step 1.
5. Snap plug into place over access hole.
6. Repositioning stroke adjustment allows angle of rotation to be adjusted from 10° to 15° from horizontal.

---

# Replacement Parts

<b>Description</b>	<b>Part Number</b>
Circuit Breaker, 1 Amp:	330-158-00
Circuit Breaker, 5 Amp:	330-118-00
Cordset for 4637Q and 4637JPNQ:	470-187-00
Cordset for 4637-1CEQ:	CRX77
Cordset for 4637-1CECN:	CRMX10
Cover:	720-505-00
Rubber Feet (4):	790-214-00
Heater, 100/120V, 200 Watts:	340-002-00
Speed Control Knob:	560-225-00
Mat, Rubber:	790-380-00
Motor:	370-341-00
RPO Motor Brushes:	370-341-01
Lamp Base, Green:	360-249-01
Lens, Green:	360-238-00
Speed Control, 120V:	227-841-00
Speed Control, 240V:	229-419-00
Switch, Rocker:	440-373-00
Configured Temperature Controller:	485-360-12
RTD Temperature Sensor:	410-632-00
Solid State Relay:	400-233-00
Wiring Schematic, 100/120V Unit:	229-235-00
Wiring Schematic, 240V Unit:	229-299-00

---

# Ordering Procedures

Please refer to the Specification Plate for the complete model number, serial number, and series number when requesting service, replacement parts or in any correspondence concerning this unit.

All parts listed herein may be ordered from the **Thermo Scientific** dealer from whom you purchased this unit or can be obtained promptly from the factory. When service or replacement parts are needed we ask that you check first with your dealer. If the dealer cannot handle your request, then contact our Customer Service Department at 563-556-2241 or 800-553-0039.

Prior to returning any materials, please contact our Customer Service Department for a "Return Materials Authorization" number (RMA). Material returned without an RMA number will be refused.

---

# One Year Limited Warranty

This Thermo Scientific product is warranted to be free of defects in materials and workmanship for one (1) year from the first to occur of (i) the date the product is sold by the manufacturer or (ii) the date the product is purchased by the original retail customer (the "Commencement Date"). Except as expressly stated above, the MANUFACTURER MAKES NO OTHER WARRANTY, EXPRESSED OR IMPLIED, WITH RESPECT TO THE PRODUCTS AND EXPRESSLY DISCLAIMS ANY AND ALL WARRANTIES, INCLUDING BUT NOT LIMITED TO, WARRANTIES OF DESIGN, MERCHANT ABILITY AND FITNESS FOR A PARTICULAR PURPOSE.

An authorized representative of the manufacturer must perform all warranty inspections. In the event of a defect covered by the warranty, we shall, as our sole obligation and exclusive remedy, provide free replacement parts to remedy the defective product. In addition, for products sold within the continental United States or Canada, the manufacturer shall provide free labor to repair the products with the replacement parts, but only for a period of ninety (90) days from the Commencement Date.

The warranty provided hereunder shall be null and void and without further force or effect if there is any (i) repair made to the product by a party other than the manufacturer or its duly authorized service representative, (ii) misuse (including use inconsistent with written operating instructions for the product), mishandling, contamination, overheating, modification or alteration of the product by any customer or third party or (iii) use of replacement parts that are obtained from a party who is not an authorized dealer of Thermo Scientific products.

Heating elements, because of their susceptibility to overheating and contamination, must be returned to the factory and if, upon inspection, it is concluded that failure is due to factors other than excessive high temperature or contamination, the manufacturer will provide warranty replacement. As a condition to the return of any product, or any constituent part thereof, to the factory, it shall be sent prepaid and a prior written authorization from the manufacturer assigning a Return Materials Number to the product or part shall be obtained.

IN NO EVENT SHALL THE MANUFACTURER BE LIABLE TO ANY PARTY FOR ANY DIRECT, INDIRECT, SPECIAL, INCIDENTAL, OR CONSEQUENTIAL DAMAGES, OR FOR ANY DAMAGES RESULTING FROM LOSS OF USE OR PROFITS, ANTICIPATED OR OTHERWISE, ARISING OUT OF OR IN CONNECTION WITH THE SALE, USE OR PERFORMANCE OF ANY PRODUCTS, WHETHER SUCH CLAIM IS BASED ON CONTRACT, TORT (INCLUDING NEGLIGENCE), ANY THEORY OF STRICT LIABILITY OR REGULATORY ACTION.

For the name of the authorized Thermo Scientific product dealer nearest you or any additional information, contact us:

North America: USA/Canada +1 866 984 3766 (866-9-THERMO) [www.thermo.com](http://www.thermo.com)

Europe: Austria +43 1 801 40 0, Belgium +32 2 482 30 30, France +33 2 2803 2180, Germany national toll free 08001-536 376,

Germany international +49 6184 90 6940, Italy +39 02 02 95059 434-254-375, Netherlands +31 76 571 4440, Nordic/Baltic countries +358 9 329 100,

Russia/CIS +7 (812) 703 42 15, Spain/Portugal +34 93 223 09 18, Switzerland +41 44 454 12 12, UK/Ireland +44 870 609 9203

Asia: China +86 21 6865 4588 or +86 10 8419 3588, India toll free 1800 22 8374, India +91 22 6716 2200, Japan +81 45 453 9220,

Other Asian countries +852 2885 4613 Countries not listed: +49 6184 90 6940 or +33 2 2803 2180