

Digital Microplate Shaker 88882005 & 88882006

In the United States:

To fax an order, use 1-800-926-1166 To order online: thermofisher.com

In Canada:

For customer service, call 1-800-766-7000 For customer service, call 1-800-234-7437 To fax an order, use 1-800-463-2996 To order online: thermofisher.ca

Find out more at thermofisher.com



Operating Manual Revision A . 09 03 2019



Contents

Section 1 Important Information	1
Disposing of Product	2
Section 2 Inspection and Installation	3
Packing List	3
Connections	4
Structure Diagram	4
Section 3 Overview	5
Specifications	5
Environmental Conditions	5
Safety Instructions	6
Capacity and Speed	7
Section 4 Operation	8
Control Panel	8
Installation	9
Settings	9
Accessory Installation	10
Section 5 Safety Tips and Maintenance	11
Section 6 Troubleshooting	12
Section 7 Optional Accessories	13
Section 8 Warranty Information	15

Section 1 Important Information

Ignoring the following warnings could cause serious injuries or even fatal accidents

Check the voltage, phase and capacity of power supply on the ID plate before installation. Connect properly.

Power supply must be properly grounded. Abnormal grounded connection causes serious damage. Grounded connection must not be on the water pipe and gas pipe.

Use provided power cord. Power cord: Wall outlet with grounded terminal power cord 250V 10A.

Do not install the product in a place that gas could leak. Do not use in a place that has industrial oil smoke or metallic dust. It causes fire or electric shock. Do not use the machine near to places where explosion could happen due to organic evaporating gases.

Explosive materials: acid, esther, nitro compound.

Inflammable materials: salt peroxides, inorganic peroxide, salt acids.

Check equipment for permissible environmental conditions when using inside of Temperature and Humidity Chamber or Incubator. It can be the cause of fire or trouble by stirrer electricity, electronic, and damage of motor.

Shaker's permissible environmental condition. Temperature 5°C to 40°C, Maximum relative humidity 80%.

Unplug if there is a strange sound, smell and/or smoke from the product. Stop

operating and request the service.

Keep out of the direct sunlight. It may influence product life and proper operation.

Do not use the machine at places where moisture is high and flooding can happen.

Do not assemble, repair, modify on your own. The product may not work well and electric shock is possible with changes in the efficiency of the product. Also this will void the warranty.

Indicates a hazardous situation which, if not avoided, may result in minor or moderate injury.

Do not put heavy things on the power cord. Do not put the machine on the cord. It may take off the wire coating and cause electric shock or fire.

Do not touch it with wet hands and place the main plug correctly. It could cause the electric shock or injuries.

Installing power outlet near instrument may be convenient

Do not install the stirrer near machinery generating high frequency noise. Avoid installation close to high frequency-welding machine, sewing machine, or mass SCR controller.

Do not inject any liquid and inflammable things inside of product.

Do not pour water or put liquid on the top of the product when cleaning. Disconnect the main power immediately and request the service if water may be in the product.

Do not let the product take any strong shock or vibration. It could cause abnormal operation or trouble. It may deteriorate the ability of the product operation and not obtain correct results.

Do not sprinkle insecticide or flammable spray on the product. Use smooth cloths. Cleaning with solvent can cause fire and deformity.

Power off while product cleaning. It may cause electric shock or fire.

Do not drop or allow the machine to fall. It will cause wrong operation and malfunction.

Disposing of Product

Dispose the unit with separating plastic mold, and motor.

Section 2 Inspection and Installation

Before unpacking the unit, first check for damages in the package of the unit.

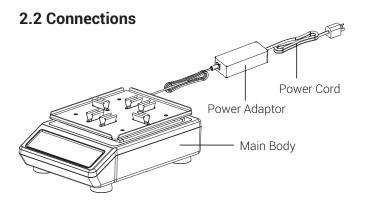
Then unpack the unit. Check carefully to see if there were damages incurred during transit.

After unpacking, check that all unit parts and accessories are as listed below. Contact us or the agent from which you purchased the unit if any components were omitted.

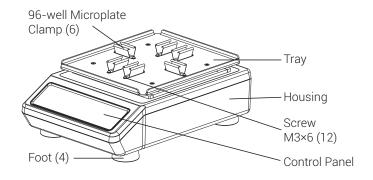
2.1 Packing List

Table -1. Packing List

Description	Catalog Number		Figure
Digital Microplate Shaker	88882005	88882006	
Tray+96-well Microplate Clamp×6 (Installed)	1	1	
General Power Adaptor	1	1	
US Plug	1	N/A	
CN Plug	N/A	1	
EU Plug	N/A	1	
UK Plug	N/A	1	
Screw Driver	1	1	
M3×6 Screw (Installed)	12	12	



2.3 Structure Diagram



Section 3 Overview

3.1 Specifications

Rotation Speed	Speed Range
Load	Maximum Load1kg @1200rpm/ 4kg @≤400rpm(accessory included) Maximum Capacity(Centered on tray)96-well Microplate 4 or 6 depends on different platforms
Time	Timing Range0, 1min.~99h59min.
Size	Overall Dimensions383×260×145mmTray Dimensions275×235×14mmPackaging Dimensions487×387×264mm
Weight	Net Weight
Power Supply	RequirementAC100-240V, 50/60Hz, 72VA
Others	CertificationRoHS, WEEE, cCSAus, CE Mark

3.2 Environmental Conditions

Application Environmental Co	onditions: indoor use
Temperature	5 to 40℃
Altitude	≤2,000 m
Humidity	20% to 85%
Storage Environmental Cond	itions
Temperature	0 to 60°C
Altitude	≤2,000 m
Humidity	20% to 90%, non-condensing

3.3 Safety Instructions

Please read the entire instruction manual before operating the Digital Microplate Shaker.

 Λ

MARNING DO NOT use the Digital Mircopalte Shaker in a hazardous atmosphere or with hazardous materials for which the unit was not designed. Also, the user should be aware that the protection provided by the equipment may be impaired if accessories used are no provided or recommended by the manufacturer, or are used in a manner not specified by the manufacturer.

caution! To avoid electrical shock, completely cut off power to the unit by disconnecting the power cord from the unit or unplug from the wall outlet. Disconnect unit from the power supply prior to maintenance and service. Any spills should be removed promptly. Bio hazard spills should be cleaned using approved liquid promptly. Solvent spills are a fire hazard. Stop the unit immediately, and DO NOT operate until clean up is complete and vapors have dissipated.

DO NOT immerse the unit for cleaning.

DO NOT operate the unit if it shows

signs of electrical or mechanical damage.

Position of Loads

Please place loads symmetrically during operation.

Symmetrical placement













Section 3 | Overview Digital Microplate Shaker

3.4 Capacity and Speed

Load Type	Liquid Capacity	Cat. No.	Max Speed
96-well Microplate	1/2-well deep		1200rpm
96-well Microplate	2/3-well deep		800rpm
Fixed Tube Rack 50ר11mm 1.5ml tube		88882117	500rpm
Fixed Tube Rack 40ר18mm 15ml tube		88882118	450rpm
Fixed Tube Rack 21ר30mm 50ml tube		88882119	400rpm

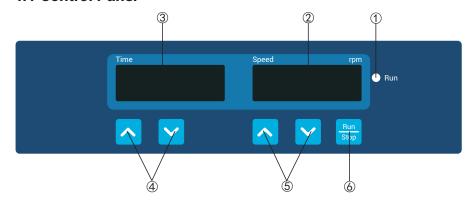
Note: These max. speeds are just for reference. We did the tests on the marble surface. Its friction coefficient is 0.32

Warning: The rotation speed is inversely proportional to the load. When the instrument is running, it is recommended to adjust the rotation speed from low to high step by step and run the instrument at an appropriate speed to avoid spillage of liquids.

Section 4 Operation

This chapter covers the control panel and its operation.

4.1 Control Panel



- ①. Run indicator: The light is on when the instrument is running and off when the instrument is in standby.
- ②. Speed display window: The window displays set speed (when the instrument is in standby) or current speed (when the instrument is running).
- ③. Time display window: The window displays cumulative time (in continuous mode) or remaining time (in timer mode).
- ④. Time setting buttons: UP/DOWN arrow buttons are used to increase/ decrease the set time of the instrument.
- (5). Speed setting buttons: UP/DOWN arrow buttons are used to increase/ decrease the set speed of the instrument.
- **(6**). Run/Stop button: Start or stop the instrument.

4.2 Installation

- 1. Connect all the components according to the figures shown on page 4 of this manual. Use grounded power outlet.
- 2. Press the power switch "I" side and switch on the instrument.

4.3 Settings

Time Settings

1. Continuous mode

Press the "\[\times \]" or "\[\times \]" arrow button below the Time display window. When the number shown on the Time display window starts flashing, press "\[\times \]" arrow button to decrease the time to 00:00 and then release the button. The time setting is finished after the number shown on the Time display window flashed twice.

2 Timer mode

Press the "\(\)" or "\(\)" arrow button below the Time display window. When the number shown on the Time display window starts flashing, press "\(\)" or "\(\)" arrow button to increase or decrease the time value. Release the button when the time shown on the Time display window reaches the set value. The time setting is finished after the number shown on the Time display window flashed twice.

Speed Settings

Press the "\sqrt " or "\sqrt " arrow button

below the Speed display window. When the number shown on the Speed display window starts flashing, press "\sigma" or

"
" arrow button to increase or
decrease the speed value. Release the
button when the speed shown on the
Speed display window reaches the set
value. The speed setting is finished after
the number shown on the Speed display
window flashed twice.

Note: Press the "\(^{"}\)" or "\(^{"}\)" arrow button for a longer time to accelerate the setting.

Run and Stop

Press "" button and the instrument will start running with the specified settings and the Run indicator light will be on.

The Time display window will show the cumulative time (Continuous Mode), or remaining time (Timer Mode) and the Speed display window will show the current speed.

Press " button again and the instrument will slow down until it stops. The instrument will then be in standby and the two display windows will show the set values.

Note: To ensure shaking operation smooth and steady, it may take 1 minute for the microprocessor control system to accelerate the tray to the set speed.

Finish Operation

After the operation is finished, press the power switch at the back right side of the instrument and put it into the "O" state. Unplug the instrument and store the instrument according to the storage guide.

Alarm System

Exceed the speed limit: After entering the operation mode for 5s, if the actual speed of the instrument is 0 or more than 1200rpm, the instrument will alarm immediately.

Err1: After entering the operation mode for 10s, and the instrument runs steadily (running stability means that the actual speed of the instrument is within the set value of ±10rpm). If the speed exceeds the set speed of ±10rpm, the instrument stops running and buzzes to alarm, Run indicator light goes off, and "Err1" is displayed in the Speed display window.

End of timer: The instrument buzzing alarm, the "End" is displayed in the Speed display window.

When the instrument alarms, press any key, the instrument is back to the standby mode.

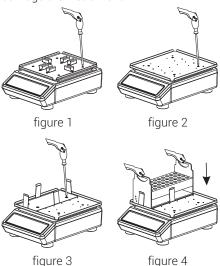
Power Recovery

If the power supply is cut off suddenly while the instrument is in operation, the unit will automatically run at the previously set parameter upon power restoration. The display windows will flash. Press any button to stop flashing.

4.4 Accessory Installation

- 1. Loosen the screws of the tray with a cross screwdriver (figure 1), and replace the standard tray with platform large (figure 2).
- 2. Split the Fixed Tube Rack and the holder (88882117, 88882118, 88882119), put the holder on tray and fix it with the 4 sunk screws (figure 3).
- 3. Vertically clip the Fixed Tube Rack into the holder and make sure there is no gap between the rack and the holder (figure 4).
- 4. Insert the test tube.

Note: When using single test tube, it is recommended not to set the speed too high because the imbalanced load may damage the instrument.



Section 4 | Operation Digital Microplate Shaker

Section 5 Safety Tips and Maintenance

Safety Tips

- 1. Use independent power supply.
- 2. Check if the local power supply voltage is suitable for use.
- 3. Do not drag the power supply cable when unplugging.
- 4. Do not use non-specified power cable or damage cable.
- 5. Service should only be performed by a qualified professional.
- 6. The power supply must be unplugged under the following situations:
- (1). When the unit is moved
- (2). When the electrical cabinet or the moving component is opened
- (3). When the equipment is malfunctioning
- (4). When the equipment is not in use

Maintenance

- a. This instrument uses brushless DC motor. It is maintenance free and has a long service time, high quality, and low noise level.
- b. Surface can be cleaned with a mild detergent and water.

Clean Spill

If accidental spillage of liquids caused by mishandling or contained breakage occurs on the surface of the instrument, please shut down the instrument and clean up the liquid immediately. If the liquid has already spilled into the unit, cut off the power supply first and immediately clean up the liquid at the surface of the instrument. Place the instrument in a ventilated and dry environment for 24 hours before reuse. If the instrument is not functioning after drying for 24 hours, please contact the manufacturer.

Warning: Disassembling/Assembling without a qualified professional's guidance may cause malfunctioning of the instrument.

Section 6 Troubleshooting

Please refer to the following table to troubleshoot if any malfunction occurs.

If the problem still exists, contact your local sales representative.

Error	Cause	Solution	
Cannot start instrument, display window off	Power disconnected	Connect the power	
	Power switch off	Switch on power	
	Power adaptor failure	Replace power adaptor	
No shaking of the tray	Over-weighted or unbalanced load	Adjust the weight and position of load,decrease rotation speed	
	Electrical malfunction	Contact Thermo Scientific	
	Mechanical malfunction	Contact Thermo Scientific	
Loud noise	Microplate loose	Adjust position of the microplate clamp	
Loud Hoise	Tray loose	Fasten screws	
Other	Keep record for maintenance		

Note:

Err1-Speed alarm

If Err1 occurs, please contact Thermo Scientific Customer Service for solutions.

Section 7 Optional Accessories

Description	Cat. No.	Dimensions	Max. Qty.	Figure
Platform Large (For 6 x Microplates)	88882114	292x275x14mm	1	
96-well Microplate clamp	88882115	54x19x19mm (7pcs/pack)	7	
Platform Large Assemble (For 6 x Microplates)	88882116	292x275x25mm	1	
Fixed Tube Rack 50ר11mm 1.5ml tube	88882117	197x86x102mm	3	N.
Fixed Tube Rack 40ר18mm 15ml tube	88882118	262x112x140mm	2	
Fixed Tube Rack 21ר30mm 50ml tube	88882119	262x112x140mm	2	
General Power Adaptor w/US plug	88870126	125VAC 10A 1.8n	n 1	
General Power Adaptor w/AUS, CN plug	88870127	250VAC 10A 1.8n	n 1	
General Power Adaptor w/EU plug	88870128	250VAC 16A 1.8n	n 1 '	
General Power Adaptor w/UK plug	88870129	250VAC13A 1.8m	1	

Description	Cat. No.	Dimensions	Max. Qty.	Figure
Screw GB/T819.1 M4×8 (Platform installation)	88882130	4ea/pack	/	
Screw GB/T819.1 M5×10 (Tube rack installation)	88882131	4ea/pack	/	
Screw GB/T9074.4 M3×6 (96-well Microplate clamp installation)	88882132	2ea/pack	/	

13 Section 7 | Optional Accessories Digital Microplate Shaker 14

Section 8 Warranty

THERMO FISHER SCIENTIFIC STANDARD PRODUCT WARRANTY

The Warranty Period starts two weeks from the date your equipment is shipped from our facility. This allows for 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 first year warranty period.

During the first two (2) years, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, labor included. Installation and calibration are 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, glass, filters and gaskets are excluded from this warranty.

Replacement or repair of components parts or equipment under this warranty shall not extend 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. At Thermo's option, all non-conforming parts must be returned to Thermo Fisher Scientific postage paid and replacement parts are shipped FOB destination.

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 1-866-984-3766, option number 2. We're ready to answer your questions on equipment warranty, operation, maintenance, service and special application. Outside the USA, please contact local Thermo Technical Services Department or local distributor for warranty information.

15 Section 8 | Warranty Digital Microplate Shaker 1