

Bottle/Tube Roller

Catalog #88881-003 and 88881-004

Operation Manual 9240-11-022



MANUAL NUMBER 9240-11-022

0		11/12/13	Original	CCS
REV	ECR/ECN	DATE	DESCRIPTION	Ву

Thermo Scientific Bottle/Tube Roller

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 Direct

1-800-438-4851 Toll Free, U.S. and Canada

1-877-213-8051 FAX

service.led.marietta@thermofisher.com Tech Support Email Address www.unitylabservices.com 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 401 Millcreek Road, Box 649 Marietta, OH 45750

International customers, please contact your local Thermo Scientific distributor.

Thermo Scientific Bottle/Tube Roller

Table of Contents

Section 1	Unpacking and Installation	1 -1
	Packing List	1-1
Section 2	Overview	.2 -1
Section 3	Operation	.3 -1
Section 4	Troubleshooting	.4 -1
Section 5	Packing List	.5 -1
Section 6	Warranty Information	6-1

iii

Section 1 Unpacking and Installation

When unpacking, follow the steps below.

- 1. Unpack and inspect for any damage that may have occurred during shipping. Call the seller if damage is found.
- 2. Verify contants against packing list and call the seller if anything is missing.
- 3. Place the unit on level and sturdy surface to avoid vibration and noise.
- 4. Check if the specified instrument voltage is identical with local power supply before powering on.

Packing List

Description		Catalog Number	Qty.	Remark
Bottle\Tube Roller		88881003 US 88881004 GBL	1	
Roller bar		88881010	6	
Power cables			1	
Operation Manual		9240-11-022	1	
O-Ring (23.7X4.1mm)		88881011	20	
O-Ring (13.8X3.1mm)		88881012	20	
Control box		88881014	1	
Stacking	Stacking Rods	88881013	4	
Accessories (Optional)	Knob		4	
	Allen Wrench		1	

Thermo Scientific Bottle/Tube Roller 1-1

Installation

1. Control box installation

This unit is equipped with a separate control box. Fasten the nut on the connection plug of control box to the unit.

Note Keys on the panel of the unit will become invalid when control box is connected. One control box controls one unit only. ▲

2. Roller bars removal and installation

Roller bars can be removed and installed for different sized tubes.

2.1 Roller bars removal

Counterclockwise twist the knob to be loose on the unit end(see Figure 1-1) and then uplift the pressure sheet-metal. Uplift the roller bar (see Figure 1-2) and remove it. Press the sheet metal to the original position and clockwise twist the knob for fastening.





Figure 1-1. Loosen the Knob

Figure 1-2. Lift Roller Bar

2.2 Roller bars installation

Counterclockwise twist the knob to be loose on the unit end (see Figure 1-1) and then uplift the pressure sheet-metal. Aim hexagon hole of roller at hexagon shaft of the unit and press the other head into semicircle orifice. Press the sheet metal to the knob position and clockwise twist the knob for fastening.

3. O-Rings

When rolling test tubes, tubes, tubes may roll off the bars due to the larger outer diameter of the tube cap. Place o-rings on the other end of the tube (see Figure 1-3) to prevent the tube from falling off. Two o-rings for each tube are recommended.



Figure 1-3. O-rings

1-2 Bottle/Tube Roller Thermo Scientific

Installation (continued)

4. Installation for stacked units

For stacking two units, optional accessories are necessary. The two units can be stacked together.

Loosen the inner hexagon screw on the bottom of the unit, pull the round head of the inner hexagon screw out as the arrow direction (see Figure 1-4) until a support base is pulled out (see Figure 1-5), then tighten the inner hexagon screw. Repeat Step a for other three support bases to be extended.





Figure 1-4. Loosen Inner Hexagon Screw

Figure 1-5. Support Base Out

- b. Locate the four stainless pillars in the stacking accessories. Screw the head with the M5 screw thread into extended support base in Step a (see Figure 1-6).
- Screw the other head with M4 screw thread into support base hole of the stacked unit and fasten the knob (see Figure 1-6. Pillar Figure 1-7). Repeat Step c to install remaining three support bases (see Figure 1-8).





Figure 1-7. Tighten Knob after Installation



Figure 1-8. 4 Pillars into Support Bases

Thermo Scientific Bottle/Tube Roller 1-3

Section 2 Overview

Features

- Wide speed range: 0.5-80 rpm
- Timing function (control box connected)
- Remote control (control box connected)
- Stackable function

Specifications

Roller Bars	6		
Roller Speed	0.5 ~ 80rpm		
Roller Bars Dimensions	Ø28mm×259mm		
Available Tubes or Bottles	Max. Ø120mm bottles; tubes		
Dimensions(W×D×H)	360mm×402mm×72mm		
Timing Range	0 ~ 999min or continuous run		
Max. Load	6.5Kg		
Power Requirement	10W		
Ambient Temperature	4°C~60°C		
Relative Humidity	≤95%		
Net Weight	6.5Kg		
Stacking Function	Optional		
Power Supply	AC100~240V,50~60Hz		

Notice

- Use the unit in specified environment.
- Verify specified instrument voltage is identical with local power supply.
- DO NOT move the unit while in operation.
- DO NOT use to mix flammable or volatile liquid or operate the unit under hazardous condition.

Thermo Scientific Bottle/Tube Roller 2-1

Section 3 Operation

Interface Instructions

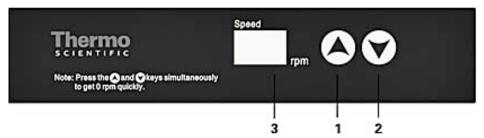


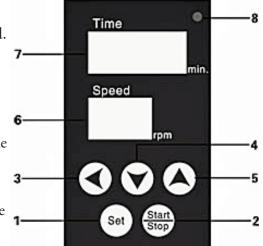
Figure 3-1. Interface Components

- 1. "▲" : To increase value of roller speed.
- 2. "▼" : To decrease value of roller speed.
- 3. Speed display: Show current speed of roller.

The speed of the Mini Roller will drop to 0 RPM when the down arrow key and up arrow key are pressed simultaneously.

Control Box Instructions

- 1. To set value of time and speed.
- 2. To start or stop the unit.
- 3. To move the cursor.
- 4. To decrease value. Hold for 3 seconds to quickly decrease the value.
- 5. To increase value. Hold for 3 seconds to quickly increase the value.



6. Displays current roller speed. Figure 3-2. Control Box Components

- Displays remaining hours.
 When it reads "00", Continuous working Mode is activated. Note
 Timing function is deactivated when the unit is in Continuous
 working Mode.
- 8. Power Indicator.

Thermo Scientific Bottle/Tube Roller 3-1

Section 4 Troubleshooting

Error	Cause	Solution
Unit does not start	Power cable not connected	Check that power cable is firmly plugged into wall outlet
	Power supply	Check that cable is correctly installed into the unit
	Control box has been connected	Remote control box from unit
No display	Power failure	Ensure unit is powered on
	Time cycle completed	Repeat run
LED light on, unit not working	Overloaded	Check and repeat run
Not-responsive keypad	Control box is connected	Remote control box

Thermo Scientific Bottle/Tube Roller 4-1

Section 5 Packing List

Description		Cat. Number	Qty.	Remark
Bottle/Tube Roller		88881003 US 88881004 GLBL	1	
Roller bar		88881010	6	
Power cables			1	
Operation Manual		9240-11-022	1	
O-Ring(23.7X4.1mm)		88881011	20	
O-Ring(13.8X3.1mm)		88881012	20	
Control box		88881014	1	
Stacking	Stacking Rods	88881013	4	
Accessories (Optional)	Knob		4	
(Sprishur)	Allen Wrench		1	

Thermo Scientific Bottle/Tube Roller 5-1

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 1-800-438-4851 (USA and Canada) or 1-740-373-4763. We're ready to answer your questions on equipment warranty, operation, maintenance, service and special application. Outside the USA, contact your local distributor for warranty information.

Rev. 0 9/13

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

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 +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

Thermo Scientific Bottle/Tube Roller 6

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

www.thermofisher.com

