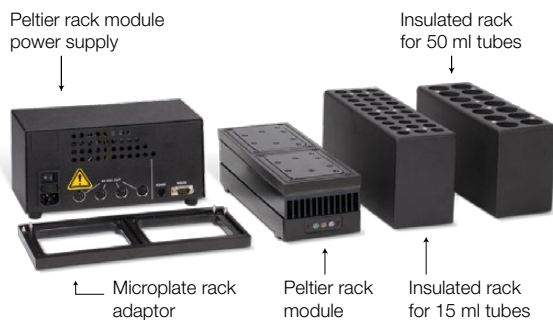


Installing the Peltier Cooling Apparatus

Introduction

The Peltier cooling apparatus enables fractions to be kept cool on the benchtop without the need to place the NGC Fraction Collector in a refrigerator. This quick guide shows you how to install the apparatus on your NGC Fraction Collector.



Installation Instructions

Ensure the NGC Fraction Collector gantry is in the home position and the system is powered down. To install a Peltier rack module (PRM) and insulated rack for 15 or 50 ml tubes, follow the instructions in steps 1–4. To install a microplate rack adaptor on a PRM, follow the instructions in steps 5–9.

- 1 Raise the dispense head to its highest position to avoid contact while placing the PRM and insulated rack on the fraction collector tray.



- 2 Place up to four PRMs on the fraction collector tray with the power cord connectors facing left.



- 3 Connect a power cable to each PRM in place on the NGC Fraction Collector. Use the cable guides on the metal tray insert to direct the cables to the rear of the NGC Fraction Collector, then connect them to the power supply. Position the cables so that movement of the gantry is not obstructed.



- 4 Place the insulated racks on the PRMs and turn on the power supply. The green LED will illuminate, indicating the power supply is on. The blue LED will flash as the PRM is cooling and will become constant once the target temperature has been reached.



- 5 To cool deep-well microtiter plates, a microplate rack adaptor must be installed on the PRM. The screws on the top of the PRM must be removed before installing the adaptor.

- 6 Ensure the PRM is disconnected from the power supply before attempting to remove the screws. Carefully remove the eight retaining screws from the top of the PRM using a flat-head screwdriver.



- 7 Place the screws to one side. Firmly grasp the frame on the PRM and lift it off.
- 8 Place the microplate rack adaptor on the PRM and replace the eight retaining screws.

- 9 Once the microplate rack adaptor is secured, raise the dispense head on the NGC Fraction Collector to its highest position, place the PRM on the tray, and connect it as described in steps 2 and 3.

- 10 Before using the NGC Fraction Collector, ensure the dispense head height is set correctly to avoid any contact between the head and the collection vessels. The optimal height is within 10 mm of the tube/plate to ensure accurate drop collection. When not in use, turn off the power to the Peltier rack module power supply to avoid condensation forming on the PRMs.

4-Port Standard Power Supply (PRPS)

Input: AC Voltage, Frequency 100–240 V, 47–63 Hz
3.6 A at 100 V AC, 1.8 A at 200 V AC

Current Installation Category CAT II (line voltage in appliance and to wall outlet)

Input: DC Voltage, Current 24 V DC, 2 A at each connector
Maximum output current (sum of all connectors): 13 A
Connect only to the supplied Peltier Rack Modules.

Peltier Rack Module (PRM)

Input: DC Voltage, Current 24 V, 2 A 

Installation Category Cat I (Mains isolated)
Use only with the provided power supply.
Connect only with the supplied cable.

Replacing the PRPS Fuse

Replace only with a 5 A 250 V slow-blow 5 x 20 mm fuse (T5AL250V).
Using an incorrect fuse may cause fire or personal injury.



Ordering Information

Catalog #	Description
12003748	Peltier Rack Power Supply , 4 channels
12003749	Peltier Rack Module
12003751	Insulated Rack , for 15 ml tubes
12003759	Insulated Rack , for 50 ml tubes
12003747	Microplate Rack Adaptor

Visit bio-rad.com/web/NGCFractionCollector for more information.

Manufactured by Teledyne CETAC Technologies.

사용자안내문

Warning Statement

EMC Registration is done on this equipment for business use only.
It may cause interference when the product would be used in home.

This warning statement applies to a product for business use.

사용자안내문

이 기기는 업무용 환경에서 사용할 목적으로 적합성평가를 받은 기기로서 가정용 환경에서 사용하는 경우 전파관섭의 우려가 있습니다.

사용자 안내문은 "업무용 방출통신기자재"에만 적용한다.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00
France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 972 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 (0)318 540 666
New Zealand 64 9 415 2280 Norway 47 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 (0) 861 246 723 Spain 34 91 590 5200 Sweden 46 08 555 12700 Switzerland 41 026 674 55 05
Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

