

Thermo Scientific™ Handheld 8-Channel Decapper

Version 7

June 2022

This page intentionally left blank.

Table of Contents

Chapter 1	Thermo Scientific™ Decapper Overview and Compatibility	1-1
	Instructions for Use.....	1-3
	Hazards and Precautions.....	1-3
	Environmental Operating Conditions.....	1-3
Chapter 2	Parts.....	2-1
Chapter 3	Basic Operation	3-1
	Power Requirements.....	3-1
	Decapping Technique	3-1
	Charging the Decapper	3-2
	Decapping.....	3-2
	Capping/Recapping	3-3
	Ejecting Caps.....	3-3
Chapter 4	Maintenance and Service.....	4-1
	Decontamination	4-1
	Service/Maintenance.....	4-1
Chapter 5	Disposal.....	5-1
Chapter 6	Warranty	6-1
Chapter 7	Regulatory Information.....	7-1
Chapter 8	Contact Information	8-1

Chapter 1

Decapper Overview and Compatibility

4105MAT

Thermo Scientific™ Matrix™ 2D ScrewTop Tubes (4105MAT)

- 4906: Matrix™ ScrewCap Tray, empty, 5 trays/case
- 4477: Matrix™ ScrewCap Tray, colorless caps, 5 trays of 96 caps/case
- 3740: 1.0mL Matrix™ 2D ScrewTop Tubes, bulk, 480 tubes/case
- 3741: 1.0mL Matrix™ 2D ScrewTop Tubes, 5 racks of 96 tubes/case
- 3742: 1.0mL Matrix™ 2D ScrewTop Tubes, uncapped, 5 racks of 96 tubes/case
- 3743: 500µL Matrix™ 2D ScrewTop Tubes, bulk, 480 tubes/case
- 3744: 500µL Matrix™ 2D ScrewTop Tubes, 5 racks of 96 tubes/case
- 3745: 500µL Matrix™ 2D ScrewTop Tubes, uncapped, 5 racks of 96 tubes/case
- Please contact your Thermo Fisher sales representative for additional product options including barcoded racks and colored cap options.

4105NUN

Thermo Scientific™ Nunc™ Universal, Nunc™ Cryobank™, Nunc™ Bank-IT™, and Matrix™ 200µL Tubes

- 3747: 200µL Matrix™ 2D ScrewTop Tubes, bulk, 480 tubes/case
- 3748: 200µL Matrix™ 2D ScrewTop Tubes, 5 racks of 96 tubes/case
- 374074: Nunc™ 0.5mL Internally-threaded Bank-It Tubes, 2D, 10 racks of 96 tubes/case
- 374078: Nunc™ 1.0mL Internally-threaded Bank-It Tubes, 2D, 10 racks of 96 tubes/case
- 374086: Nunc™ 0.5mL Internally-threaded Cryobank Tubes, 2D, 10 racks of 96 tubes/case
- 374088: Nunc™ 1.0mL Internally-threaded Cryobank Tubes, 2D, 10 racks of 96 tubes/case

- 374220: Nunc™ 5.0mL Internally-threaded Cryobank Tubes, 2D, 10 racks of 48 tubes/case
- 374220LOCK: Nunc™ 5.0mL Internally-threaded Cryobank Tubes, 2D, Lock Rack, 10 racks of 48 tubes/case
- 374221: Nunc™ 2.0mL Internally-threaded Cryobank Tubes, 2D, 10 racks of 48 tubes/case
- 374221LOCK: Nunc™ 2.0mL Internally-threaded Cryobank Tubes, 2D, Lock Rack, 10 racks of 48 tubes/case
- 374222: Nunc™ 5.0mL Internally-threaded Cryobank Tubes, 10 racks of 48 tubes/case
- 374223: Nunc™ 2.0mL Internally-threaded Cryobank Tubes, 10 racks of 48 tubes/case
- 374223LOCK: Nunc™ 2.0mL Internally-threaded Cryobank Tubes, Lock Rack, 10 racks of 48 tubes/case
- 374258: Nunc™ 2.0mL Internally-threaded Cryobank Tubes, 2D, Bulk, 960 tubes/case
- 374259: Nunc™ 2.0mL Internally-threaded Cryobank Tubes, Bulk, 960 tubes/case
- 374260: Nunc™ 5.0mL Internally-threaded Cryobank Tubes, Bulk, 528 tubes/case
- 374261: Nunc™ 5.0mL Internally-threaded Cryobank Tubes, 2D, Bulk, 528 tubes/case
- 374500: Nunc™ 1.8mL Externally-threaded Universal Tubes, 2D, White Patch and Graduations, 10 racks of 48 tubes/case
- 374501: Nunc™ 1.8mL Externally-threaded Universal Tubes, White Patch and Graduations, 10 racks of 48 tubes/case
- 374502: Nunc™ 1.8mL Externally-threaded Universal Tubes, 2D, White Patch and Graduations, Bulk, 480 tubes/case
- 374503: Nunc™ 1.8mL Externally-threaded Universal Tubes, White Patch and Graduations, Bulk, 480 tubes/case
- 374510: Nunc™ 2.0mL Internally-threaded Universal Tubes, 2D, 10 Racks of 48 tubes/case
- 374511: Nunc™ 2.0mL Internally-threaded Universal Tubes, 10 Racks of 48 tubes/case
- 374512: Nunc™ 2.0mL Internally-threaded Universal Tubes, 2D, Bulk, 480 tubes/case
- 374513: Nunc™ 2.0mL Internally-threaded Universal Tubes, Bulk, 480 tubes/case
- Please contact your Thermo Fisher sales representative for additional product options including barcoded racks and colored cap options.

Instructions for Use

The Thermo Scientific™ Handheld 8-Channel Decapper is intended to be used for capping and decapping Matrix™ and Nunc™ storage tubes during experimentation and storage.

These products are intended for general laboratory use only. They have been tested and found to comply with the limits for a Class B residential device.

Hazards and Precautions

There are no known hazards associated with the Thermo Scientific™ Handheld 8-Channel Decapper when it is operated in accordance with the instructions in this manual.

If these products are used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.

Environmental Operating Conditions

Operating Temperature: 10-35°C (50-95°F)

Relative Humidity: 10-95% Non-Condensing

Be sure to follow the rules listed below while charging the battery. Failure to do so may cause the battery to become hot, explode, or ignite and cause serious injury.

1. When Using the Battery**DANGER**

(1) Misusing the battery may cause the battery to get hot, explode, or ignite and cause serious injury. Be sure to follow the safety rules listed below:

- Do not place the battery in fire or heat the battery.
- Do not install the battery backwards so that the polarity is reversed.
- Do not connect the positive terminal and the negative terminal of the battery to each other with any metal object (such as wire).
- Do not carry or store the batteries together with necklaces, hairpins, or other metal objects.
- Do not penetrate the battery with nails, strike the battery with a hammer, step on the battery, or otherwise subject it to strong impacts or shocks.
- Do not solder directly onto the battery.
- Do not expose the battery to water or salt water, or allow the battery to get wet.

(2) Do not disassemble or modify the battery. The battery contains safety and protection devices which, if damaged, may cause the battery to generate heat, explode or ignite.

(3) Do not place the battery on or near fires, stoves, or other high temperature locations. Do not place battery in direct sunshine, or use or store the battery inside cars in hot weather. Doing so may cause the battery to generate heat, explode, or ignite. Using the battery in this manner may also result in a loss of performance and a shortened life expectancy.

(4) Do not insert the battery into equipment designed to be hermetically sealed. In some cases hydrogen or oxygen may be discharged from the cell which may result in rupture, fire or explosion.

Warning

- 1) Immediately discontinue use of the battery if, while using, charging, or storing the battery, the battery emits an unusual smell, feels hot, changes color, changes shape, or appears abnormal in any other way. Contact your sales location or Thermo Scientific if any of these problems are observed.
- (2) Do not place the batteries in microwave ovens, high-pressure containers, or on induction cookware.
- (3) In the event that the battery leaks and the fluid gets into one's eye, do not rub the eye. Rinse well with water and immediately seek medical care. If left untreated the battery fluid could cause damage to the eye.

Caution

(1) If the device is to be used by small children, the caregiver should explain the contents of the user's manual to the children. The caregiver should provide adequate supervision to insure that the device is being used as explained in the user's manual.

(2) When the battery is worn out, insulate the terminals with adhesive tape or similar materials before disposal.

Danger

Be sure to follow the rules listed below while charging the battery. Failure to do so may cause the battery to become hot, explode, or ignite and cause serious injury.

When charging the battery, either use a specified battery charger or otherwise insure that the battery charging conditions specified by ThermoFisher are met.

Do not attach the batteries to a power supply plug or directly to a car's cigarette lighter.

Do not place the batteries in or near fire, or into direct sunlight. When the battery becomes hot, the built-in safety equipment is activated, preventing the battery from charging further, and heating the battery can destroy the safety equipment and can cause additional heating, breaking, or ignition of the battery.

Warning

Do not continue charging the battery if it does not recharge within the specified charging time. Doing so may cause the battery to become hot, explode, or ignite.

2. While Charging

Caution

The temperature range over which the battery can be charged is 0°C to 45°C. Charging the battery at temperatures outside of this range may cause the battery to become hot or to break. Charging the battery outside of this temperature range may also harm the performance of the battery or reduce the battery's expectancy.

3. When Discharging the Battery

Danger

Do not discharge the battery using any device except for the specified device. When the battery is used in devices aside from the specified device it may damage the performance of the battery or reduce its life expectancy, and if the device causes an abnormal current to flow, it may cause the battery to become hot, explode, or ignite and cause serious injury.

Caution

The temperature range over which the battery can be discharged is -10°C to 60°C. Use of the battery outside of this temperature range may damage the performance of the battery or may reduce its life expectancy.

2

Chapter 2

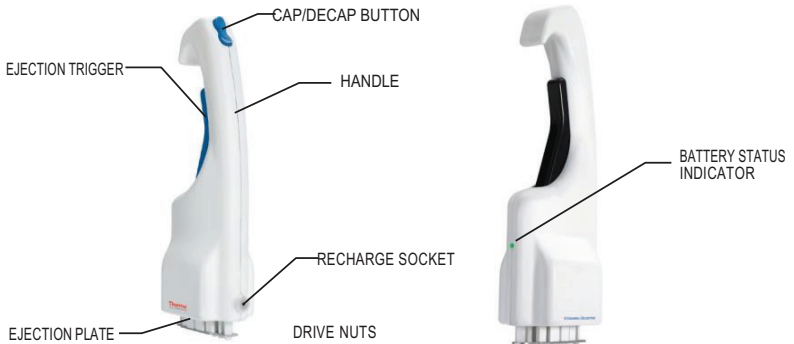
Parts

Included in this packaging should be the following:

- 1) Thermo Scientific™ 8-Channel Decapper (Part Numbers 4105MAT, 4105NUN)
- 2) Power Supply with Adapters for International Use (Part Number 4109)
- 3) Benchtop Stand (Part Number 4106)
- 4) Rack of Matrix™ or Nunc™ Caps (depending upon model ordered)
- 5) Spacer for use with Part Numbers 3747, 3748, 374074, 374078, 374086, 374088, 374500, 374501, 374502, 374503 (Only included with 4105NUN Model)
- 6) Operation Manual
- 7) Certificate of Warranty and Warranty Registration Card

If any of the above parts are missing, please contact your Thermo Fisher Scientific representative or their authorized dealer

Critical parts of the Thermo Scientific™ 8-Channel Decapper are identified below



3



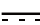
Chapter 3

Basic Operation

Charging the Decapper

- Before first-time use, charge the decapper by plugging the power supply into the recharge socket at the base of the handle.
- Simply plug the other end into a standard 100-240 volt, 50-60 Hz wall socket and allow the unit to charge up to 2 hours at which time the unit will be fully charged
- For added ease and convenience, the decapper features:
 - A bi-color LED on the front of the handle that indicates the level of battery charge:
 - LED alternatively blinking **Red** & **Green** = Charging
 - Solid **Green** = In operation or charge complete
 - Blinking **Red** = Low battery
 - Special circuitry makes it impossible to overcharge the battery and allows the decapper to be charged or partially charged at any convenient time
 - For long-term storage, leave the decapper attached to power supply

Power Requirements

- Power supply input: INPUT: 100-240V ~ 50/60Hz 0.3A
- Power supply output: OUTPUT: 12.0V @ 1.0A, constant voltage
- Explanation of symbols on power supply:
 - Indoor use only = 
 - Class 2 power supply = 
 - DC Symbol = 

Decapping Technique

- Holding the Decapper loosely with a 75° angle between the forearm and upperarm will minimize potential muscle strain





Decapping

- Place the decapper over row of tubes to be decapped, making sure that the ejector plate is parallel with tops of caps




TIP: Line up the cap driver closest to the recharge socket with the tubes that are closest to you. The other drivers and tubes will then line up

- While lightly pushing down on caps, toggle the “” button on the handle to make sure all of the drive nuts have been properly seated into the caps
- Once all of the drive nuts are seated, push the “” button on handle until the tubes are decapped

Ejecting Caps

- Once the tubes have been decapped, line up the decapper with an empty row on the cap tray
- Insert caps into cavities in cap tray
- Squeeze ejection trigger

Capping/Recapping

- Apply the caps from the cap tray onto the drive nuts. Line up the decapper with the row of tubes to be recapped, making sure the ejection plate is parallel with the tubes
- Push the “” button to cap. DO NOT push caps on. Let the caps screw onto the tubes on their own
- Once the caps are securely on the tubes, squeeze the ejection trigger to remove the caps from the drive nuts

4

Chapter 4

Maintenance and Service

Decontamination

- The entire 8-Channel Decapper can be cleaned with most common laboratory disinfectants but must not be submerged in liquid.
- This instrument contains electrical components that can not be autoclaved
- Below are the steps to disinfect:
 - Wet a cloth with the desired disinfectant
 - Wipe down decapper making sure not to drip disinfectant into the inner cavity
- Suggested Disinfectants:
 - Alcohol wipes
 - 5% Bleach solution
 - Thermo Scientific™ RNase AWAY™ disinfectant
 - Thermo Scientific™ DNA Away™ disinfectant
 - Phenolic cleansers

Service/Maintenance

- Service the decapper after each year of use.

5

Chapter 5

Disposal

Regulations and laws pertaining to the recycling and disposal of lithium ion batteries vary from country to country as well as by state and local governments. The European governments have more strict regulations on the disposal of rechargeable batteries than the USA and Canada. You will need to check the laws and regulations where you live.

6

Chapter 6

Warranty

The Thermo Scientific™ Handheld 8-Channel Decapper comes with a 12-month warranty. Please visit www.thermofisher.com/global-registration to register your product and activate your warranty. Preventive maintenance and service plans are available from Thermo Fisher Scientific. Email technicalsupport@thermofisher.com for details.

7

Chapter 7

Regulatory Information

Conformity Statements

Europe:

We herewith declare under our sole responsibility that the above mentioned products meet the provisions of the following EC Directives and harmonized standards:

- 2014/30/EU (EMC Directive)**
- 2014/35/EU (LVD Directive)**
- 2006/42/EC (Machinery Directive)**
- 2012/19/EU (WEEE Directive)**
- 2011/65/EU (RoHS Directive)**

United Kingdom:

We herewith declare under our sole responsibility that the above-mentioned products meet the provisions of the following UK Regulation(s):

- Electromagnetic Compatibility Regulations 2016**
- Electrical Equipment (Safety) Regulations 2016**
- Supply of Machinery (Safety) Regulations 2008**
- The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012**

United States of America:

FCC Compliance Statement products subject to **Part 15B**

This device complies with Part 15 of the FCC Rules. Operation is subject to the following two conditions: (1) This device may not cause harmful interference, and (2) this device must accept any interference received, including interference that may cause undesired operation

Canada:

This device complies with ICES-003 (Issue7) of the ISED Rules

CAN ICES-003(B) / NMB-003(B)

WEEE Compliance:

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on Thermo Fisher Scientific's compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive via email: technicalsupport@thermofisher.com

Conformité WEEE:



Ce produit doit être conforme à la directive européenne (2012/19/EU) des Déchets d'Équipements Électriques et Electroniques (DEEE). Il est marqué par le symbole suivant: Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles par email: techsupport.labproducts.eu@thermofisher.com

WEEE Konformität:



Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2012/19/EU erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet: Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs- / Entsorgungs- anlagen in allen EU-Mitgliedstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden.

Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Fisher Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Scientific Produkte zu identifizieren, die unter diese RoHS Anweisung fallen, bekommen Sie per email: techsupport.labproducts.eu@thermofisher.com

China RoHS Declaration

Below is the China EEP Hazardous Substances Information Table

产品中有害物质的名称及含量 **China EEP Hazardous Substances Information**

部件名称 Component Name	Thermo Scientific Handheld 8-Channel Decapper					
	有害物质 Hazardous Substances					
	铅 (Pb)	汞 (Hg)	镉 (Cd)	六价铬 (Cr(VI))	多溴联苯 (PBB) Polybrominated biphenyls	多溴二苯醚 (PBDE) Polybrominated diphenyl ethers
金工 Metal Parts	0	0	0	0	0	0
塑料 Plastic Parts	0	0	0	0	0	0
显示器 Display	0	0	0	0	0	0
电缆 Cabling	0	0	0	0	0	0
电路板 PCBA's	0	0	0	0	0	0
标签 Labels	0	0	0	0	0	0
本表格依据 SJ/T 11364 的规定编制 This table is compiled according to SJ/T 11364 standard.						
0: 表示该有害物质在该部件所有均质材料中的含量均在 GB/T 26572 规定的限量要求以下。 Indicates that the concentration of the hazardous substance in all homogeneous materials for the part is below the relevant threshold of the GB/T 26572 standard.						
X: 表示该有害物质至少在该部件的某一均质材料中的含量超出 GB/T 26572 规定的限量要。 Indicates that the concentration of the hazardous substance in at least one homogenous material of the part is above the relevant threshold of the GB/T 26572 standard.						

The China EEP Hazardous Substances Information table is available at <https://www.thermofisher.com/in/en/home/technical-resources/rohs-certificates.html>



Chapter 8

Contact Information

For technical support in North America and Europe, please contact us at the following email address for your region:

North America:

info.LH@thermofisher.com

Europe:

techsupport.labproducts.eu@thermofisher.com

For technical support outside of these regions, and for all other product inquiries, please visit www.thermofisher.com, and go to the Contact Us page under the Services menu. After choosing a support option, search for your product, and select your country/region to find the most up to date information on:

Product Information

Contact Phone Numbers

Contact Email Addresses

Regulatory Information

Instrument Service

This page intentionally left blank.

This page intentionally left blank.

Thermo Fisher Scientific Oy
Ratastie 2
01620 Vantaa
Finland

thermofisher.com/samplestorage

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries unless otherwise specified. 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
scientific