

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

- 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

2

Chapter 2

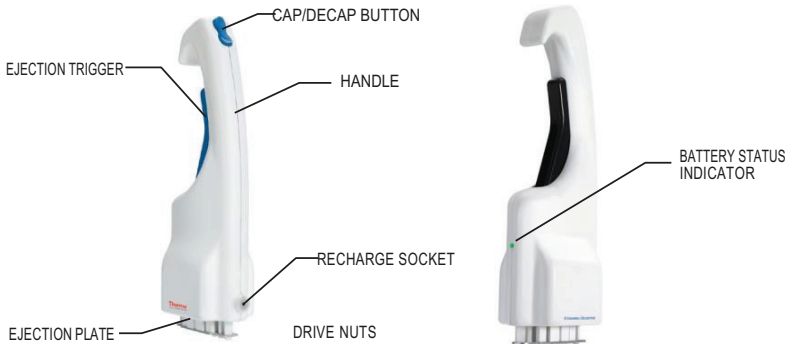
Parts

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



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.

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