

Operating Instructions

Cell Locker

50155007 Revision C May 2021



Preface

© 2021 Thermo Fisher Scientific Inc. All rights reserved.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions.

In case of conflicting translations into foreign languages the German-language version of these operating instructions shall be binding.

Trademarks

CELL LOCKER® is a registered trademark of Thermo Scientific.

Thermo Scientific is a brand owned by Thermo Fisher Scientific, Inc.

All other trademarks mentioned in the operating instructions are the exclusive property of the respective manufacturers.

Thermo Electron LED GmbH
Robert-Bosch-Straße 1
D - 63505 Langenselbold
Deutschland

Die Thermo Electron LED GmbH is an affiliate to:

Thermo Fisher Scientific Inc.
168 Third Avenue
Waltham, MA 02451
USA

Table of Contents

Chapter 1 General notes	2
1.1. Explanation of safety information and symbols	2
1.2. Safety notes	3
1.3. Operational Safety Rules	3
1.4. Warranty	3
1.5. Intended Purpose	4
1.5.1. Correct Use	4
1.5.2. Incorrect Use	4
Chapter 2 Delivery of the device	5
2.1. Scope of delivery	5
2.2. Acceptance inspection	5
Chapter 3 Description of the device	6
3.1. Description	6
3.2. First use	7
3.3. Operation	7
3.4. Application versions	7
Chapter 4 Cleaning and disinfection	9
4.1. Cleaning	9
4.2. Wipe/spray disinfection	10
4.3. Autoclaving	11
Chapter 5 Service	12
5.1. Membrane filter replacement	12
5.2. Cover gasket replacement	14
Chapter 6 Spare and wear parts	15
Chapter 7 Disposal	16
Chapter 8 Technical Data	17

1. General notes

1.1 Explanation of safety information and symbols



Warning

Indicates a hazardous situation which, if not avoided, could result in death or serious injuries.



Caution

Indicates a hazardous situation which, if not avoided, could result in minor or moderate injuries.

Notice

Indicates a situation which, if not avoided, could result in damage to equipment or property.

	In chapters of the instruction manual which deal with safety, this icon appears under the title of the chapter. Displayed on the equipment, this icon denotes that special attention must be paid to the information given in the instruction manual or accompanying documents.
	Is used for application hints and useful information
	Harmful liquids!
	Fire hazard!
	Toxic substances!
	Wear safety gloves!
	Wear safety goggles!
	Wear breathing protection!

1.2 Safety notes

The *CELL LOCKER* may only be used by instructed personnel.

The present operating instructions, applicable safety data sheets, hygiene guidelines and the corresponding technical rules issued by the operator shall be used to create written procedures targeted at personnel working with the subject matter device, detailing:

- the decontamination measures to be employed for the *CELL LOCKER* and the accessories used with it,
- the safety precautions to be taken when processing specific agents,
- wearing protective equipment when handling e.g. microbiological and biological samples,
- the measures to be taken in case of accidents,
- drying time after decommissioning, as Transport 2 hours,
- working according to GLP - Good Laboratory Practice.

1.3 Operational Safety Rules

Observe the sample weight limits as a whole and its shelving in particular.

Check the door seal every six months for proper sealing performance and possible damage.

Do not process any samples containing hazardous chemical substances that may be released into the ambient air through defective seals or may cause corrosion or other defects.

1.4 Warranty

Thermo Fisher Scientific warrants the operational safety and functions of the *CELL LOCKER* for one year only under the condition that:

- the *CELL LOCKER* is operated and serviced exclusively in accordance with its intended purpose and as described in these operating instructions,
- the device is not modified,
- only original spare parts and accessories that have been approved by Thermo Fisher Scientific are used (third-party spares without Thermo Fisher Scientific approval void the limited warranty),
- inspections and maintenance are performed at the specified intervals,
- an operation verification test is performed after each repair activity.

The warranty is valid from the date of delivery of the *CELL LOCKER* to the customer.

1.5 Intended Purpose

1.5.1. Correct Use

The *CELL LOCKER* is intended for use in laboratory equipment, which among others is used for the preparation and cultivation of cell and tissue cultures for research applications.

The maximum Autoclave temperature is 15 minutes at 121°C (250 °F) with fast exhaust.

1.5.2. Incorrect Use

Do not use tissues, substances or liquids that:

- are easily flammable or explosive;
- release vapor or dust that forms combustible or explosive mixtures when exposed to air;
- release poisons;
- release dust;
- exhibit exothermic reactions;
- are pyrotechnical substances.

Do not pour any liquids on the bottom of the interior surface or into a collecting basin inside the unit.

The *CELL LOCKER* is not for use in Medical Devices.

2. Delivery of the device

2.1 Scope of delivery

Description	Pcs.
<i>CELL LOCKER</i>	1
Membrane filter	2
Work tray with cover	1
Consumables kit	1
Operating instructions	1
Summarized safety instructions	1

2.2 Acceptance inspection

After the *CELL LOCKER* has been delivered, check the delivery immediately for:

- completeness,
- possible damage.

If components are missing or damage is found on the device or the packaging, please contact the carrier, as well as the customer service, immediately.

3. Description of the device

3.1 Description

The *CELL LOCKER* consists of a polycarbonate chamber with two membrane filters, which are permeable to water vapor and gas, but are impermeable to microorganisms.

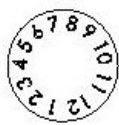


Fig. 1. *CELL LOCKER*

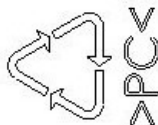
Position	Name
1	work plate
2	membrane frame with filter
3	gasket
4	latch
5	cover
6	handle knob



Production
year



Production
month

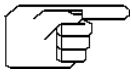


Polycarbonate
Makrolon 2528

Fig. 2. Bottom marks

3.2 First use

When using the first time, clean and disinfect the *CELL LOCKER*.



Note

Clean the *CELL LOCKER* before inserting the membrane frame with filter.



Warning

The *CELL LOCKER* must be removed before the incubator is sterilized.

3.3 Operation

The *CELL LOCKER* can be placed into the incubator with 6-door gas tight screen with or without cover.

When inserting into a laboratory device make sure that the *CELL LOCKER* is not twisted.



3.4 Application versions

Depending on user preference, the *CELL LOCKER* can be modified to attach or remove the transport cover, using the consumables kit:

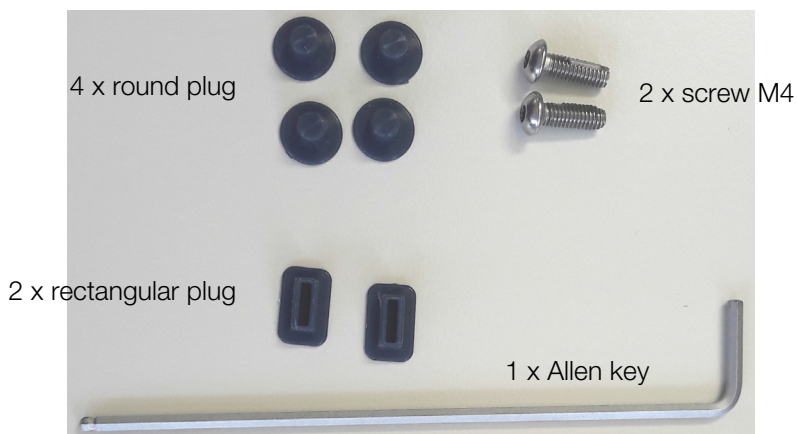


Fig. 3. Consumables kit

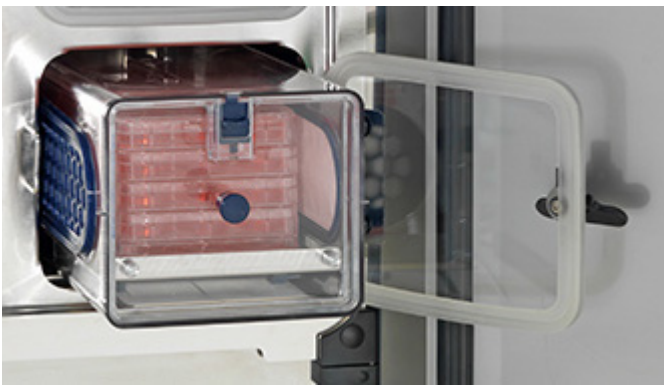
Cover can be attached to work tray with the Allen key, allowing protected transport.



with attached transport cover



without transport cover



with cover to remove *CELL LOCKER* with cultures



without cover to remove cultures with single door access



samples access inside the chamber with transport cover

Note:
Attach cover for protected transport to biosafety cabinet or to another incubator or lab. Use of cover in the chamber may require two hand opening.

Fig. 4. *CELL LOCKER* options

4. Cleaning and disinfection

Before starting any work, the following safety hints should be followed:



Warning
Health hazard!



The surfaces of the work space may be contaminated. Contact with contaminated cleaning liquids may cause infections. Disinfectants may contain harmful substances. When cleaning and disinfecting, always observe the safety instructions and hygiene guidelines.



Wear safety goggles!



Wear safety gloves!



Wear breathing protection!

Observe the safety instructions of the disinfectant's manufacturer and the hygiene supervisor!

4.1 Cleaning



Caution
Incompatible cleaners!



Some of the **CELL LOCKER** components are made of plastic.

Solvents may attack plastics.

Strong acids or alkaline solutions may cause corrosion of plastics.



Caution
Moisture-sensitive components!

Do not spray cleaning agent onto the membrane filter of the **CELL LOCKER**. When wiping the membrane filter frame make sure that no moisture contacts the membrane filter.

Remove dirt residues and deposit thoroughly using a solution of lukewarm water and mild detergent.

Wipe the surfaces and work plate with a clean cloth and clear water. Then, wipe the surfaces dry using a clean cloth.

4.2 Wipe/spray disinfection



Caution

Alcoholic disinfectants!

Disinfectants having a alcohol content of more than 10% may form, in combination with air, easily combustible and explosive gas mixtures.



When using such disinfectants, avoid open flames or exposure to excessive heat during the entire disinfection process.

Use such disinfectants only in adequately ventilated rooms.

After the disinfectant has been allowed to react, wipe the cleaned **CELL LOCKER** components thoroughly dry.

Observe safety regulations to avoid fire and/or explosion hazard caused by alcohol-containing disinfectants.



Caution

Chloride-containing disinfectants!



Do not use chloride-containing disinfectants.

Use only disinfectants that do not affect stainless steel.



Warning

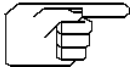
Decontamination or cleaning agents!

Thermo Electron LED GmbH or his agent is to be consulted, if there is any doubt about the compatibility of decontamination or cleaning agents.

1. Remove all samples from the work space and store them in a safe place.
2. Spray disinfectant onto the surfaces of the work space and of the accessories or wipe the surfaces clean using disinfectant. Do not spray disinfectant onto the membrane filters.
3. Allow time for the disinfectant to act as specified by the manufacturer.
4. Wipe off the chamber surfaces and the work tray using lukewarm water mixed with standard rinsing agents. Rinse completely using warm water.
5. Re-rinse the cleansed surfaces 3 to 5 times with autoclaved water to completely remove any cleaning agent residues.
6. After this, wipe the cleansed surfaces, air baffle and internals dry with a soft, sterile cloth.

4.3 Autoclaving

The *CELL LOCKER* can be autoclaved up to 15 minutes at 121°C with fast exhaust. Autoclaving at higher temperature is not recommended since the material cannot withstand higher temperature and its composition would be negatively affected.



Note

Remove the membrane frame with filter before autoclaving the *CELL LOCKER*.

5. Service

Any service work should be recorded.

5.1 Membrane filter replacement

Periodically check the membrane frame filters and at least once a year replace them.



Warning
Health hazard!



Wear safety gloves!

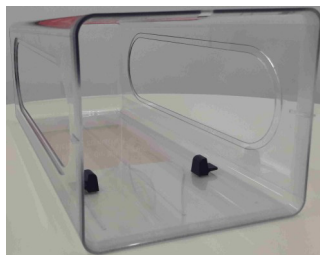
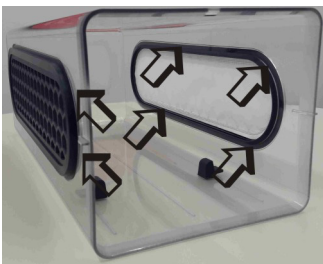


Wear breathing protection!

1. Pull out the work tray and take it out.



2. Press and remove the membrane frame from the inside.



3. Put the membranes into a plastic bag, autoclave them and dispose as hazardous waste.



4. Unpack the new membranes.



5. Insert the new membrane frame and press firmly in the radial direction.

6. Visually inspect from the inside, whether the gasket lip is correctly seated, if necessary press from the outside.



7. Insert the second membrane according to 5. and 6.



Warning Contamination hazard!

Damaged or improperly mounted membranes can contaminate the user or the environment!

8. Insert the work tray over the tilting stop and push into the interior.



5.2 Cover gasket replacement

Inspect the cover gasket periodically, e.g. monthly, for integrity, no damage and replace as needed.

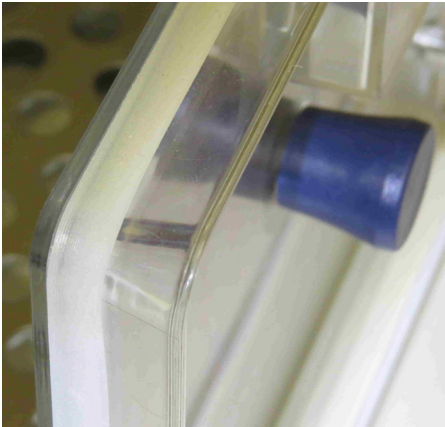
1. Pull out the work tray and take it out.



2. Remove cover gasket from the notch.



3. Insert a new gasket in the notch and take care that the gasket lip is outside.



Warning Contamination hazard!

Damaged or improperly mounted membranes can contaminate the user or the environment!

4. Insert the work tray and push into the interior.



6. Spare and wear parts

Description	Material-No.
Membrane frame's	50153148
Cover seal (6 pieces)	50156433
Work tray, stainless steel	50155808
Work tray, copper	50155809
Consumables kit	50161834
Cover <i>CELL LOCKER</i>	50156434

Only original spare parts and accessories that have been approved by Thermo Fisher Scientific used, third-party spares without Thermo Fisher Scientific approval void the limited warranty.

7. Disposal



Warning

Contamination hazard!

The device may have been used for treating and processing infectious substances. Therefore, the device and device components may have been contaminated. Prior to disposal, all device components must be decontaminated!

Clean the unit components thoroughly, then disinfect ,decontaminate or autoclave them (depending on application).

Attach a declaration of non-objection with details of performed decontamination measures to the items that are to be disposed of.

All device components can be disposed of after they have been decontaminated properly. Membrane filters must be disposed of in accordance with applicable national legislation and directives on the handling of hazardous waste.

Overview of the materials used

Component	Material
<i>CELL LOCKER</i> housing and cover	Polycarbonat, Makrolon 2858
Work tray stainless steel	1.4301+2B / Steel 10088-2
Work tray copper	Copper
Membrane frame	ABS
Membrane filter	PES Supor-200R Membrane 0,2 µm PALL

8. Technical Data

Description	Unit	Value
Outer dimensions (W x H x D)	mm	185 x 156 x 380
Inner space volume	l	approx. 5.4
Weight	kg	1.8
Loading capacity per work tray	kg	5
Operating temperature	°C / °F	4 - 55 / 39 - 131
Autoclave cycles		max. 12
Autoclave temperature	°C / °F	max. 121 / 250

Find out more at [thermofisher.com](https://www.thermofisher.com)

thermoscientific