

**Thermo Scientific**  
**Heracell VIOS 250i**  
**Heracell VIOS 250i LK**  
**Heracell Vios 250i CR**  
**Forma Steri-Cycle i250**  
**Forma Steri-Cycle i250 LK**  
**Forma Steri-Cycle i250 CR**

*CO<sub>2</sub> Incubators*

**Supplement to Operating Instructions 50144132, 50162744,  
50163129, 50144829, 50162762 and 50163206**

©2021 Thermo Fisher Scientific Inc. All rights reserved.

#### Trademarks

Heracell VIOS™, Heracell Vios CR™, Steri-Run™, Forma™, Steri-Cycle™, Steri-Cycle CR™, iCan™, THRIVE™ and Cell Locker™ are registered trademarks of Thermo Scientific.

G-Rex® is the Trademark of Wilson Wolf.

Thermo Scientific is a brand of 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

Germany

Thermo Electron LED GmbH is a subsidiary company of:

Thermo Fisher Scientific Inc.

168 3rd Avenue

Waltham, MA 02451

USA

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. These operating instructions are protected by copyright. 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 operating instructions manual may change at any time and without any prior notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

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 of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

# Table of Contents

<b>Preface.....</b>	<b>1</b>
General Notes.....	1
Instruction of the Operating Personnel.....	1
Applicability of the Instructions.....	1
Original Document for Translation .....	1
Instructions to be used .....	1
Explanation of Safety Information and Symbols.....	2
Safety Information and Symbols used in the Operating Instructions .....	2
Design of the equipment .....	2
.....	2
<b>Chapter 1 Scope of Delivery .....</b>	<b>1-1</b>
<b>Chapter 2 Installation .....</b>	<b>2-1</b>
<b>Chapter 3 Operation .....</b>	<b>3-1</b>
<b>Chapter 4 Performance Hints .....</b>	<b>4-1</b>
<b>Chapter 5 Parts overview.....</b>	<b>5-1</b>
<b>Chapter 6 Technical Data.....</b>	<b>6-1</b>

# Preface

## General Notes

## Instruction of the Operating Personnel

These operating instructions describe the shelf kit for the CO<sub>2</sub> incubators *Heracell Vios 250i*, *Heracell Vios 250i CR*, *Forma Steri-Cycle i250* and *Forma Steri-Cycle i250 CR*. The shelf kit has been manufactured in keeping with the latest technological developments and is operationally safe. However, the device may present potential hazards, particularly if the device is operated by inadequately trained personnel or if it is not used in accordance with the intended purpose. Therefore, the following must be observed to prevent accidents:

- The CO<sub>2</sub> incubator must be operated only by trained and authorized personnel.
- For any operation of this device, the operator must prepare clear and concise written instructions in the language of the operating and cleaning personnel based on these operating instructions, applicable safety data sheets, plant hygiene guidelines, and technical regulations, in particular:
  - which disinfection measures are to be applied for the device and accessories,
  - which protective measures apply when specific agents are used,
  - the measures to be taken in case of accidents,
  - which precautions and rules of conduct are necessary when entering and working in a clean room.
- Repairs to the device must be carried out only by trained and authorized expert personnel.

## Applicability of the Instructions

### Original Document for Translation

- The contents of the supplement instructions are subject to change without further notice.
- In case of conflicting translations into foreign languages the German-language version of these supplement instructions shall be binding.

### Instructions to be used

- Keep these supplement instructions and the related operating instructions close to the device so that safety instructions and important information are always accessible.

Should you encounter problems that are not detailed adequately in these supplement instructions and the related operating instructions, please contact Thermo Scientific immediately for your own safety.

# Explanation of Safety Information and Symbols

## Safety Information and Symbols used in the Operating Instructions



**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.



**CAUTION** Indicates a situation which, if not avoided, could result in property damage.

**Note** Is used for applicational hints and useful information.

## Design of the equipment

The shelves kit is designed for G-Rex 500 container.

## Scope of Delivery



1 x 50166349 Bottom shelf

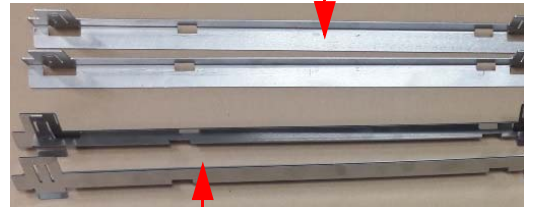


4 x 50166348 Split shelf



4 x 50166347 Guiding rail carrier

2 x 50166346 Guiding rail left



2 x 50166345 Guiding rail right



1 x 50166350 Reinforcement bottom shelf

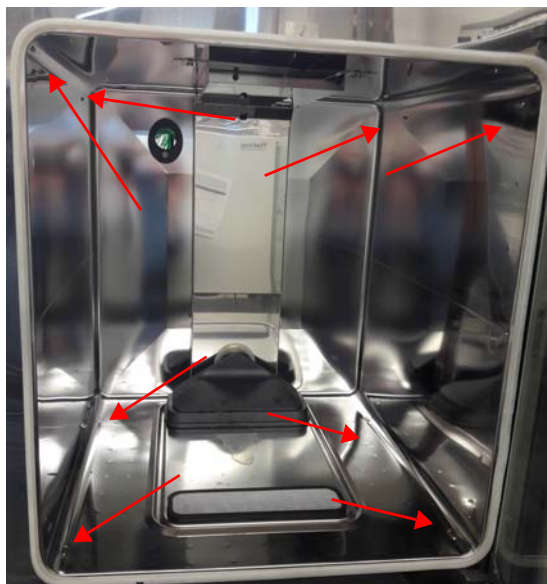
## Installation

**Note**

The parts are delivered with a protective foil. To remove the protective foil more easily warm-up the parts slightly.

**Insert the guiding rail carriers**

- In the inner chamber there are intakes for 4 guiding rail carriers.





- Place the guiding rail carrier in the lower intake.



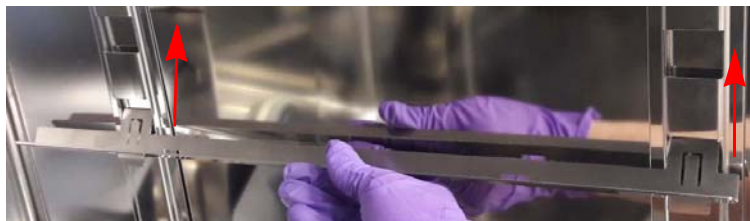
- Snap in or push in the guiding rail carrier in the upper intake.



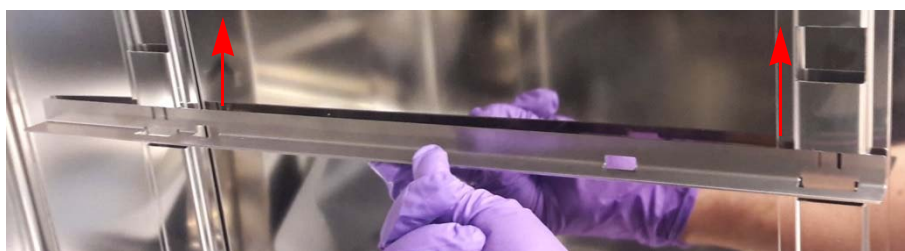


### Insert the guiding rails

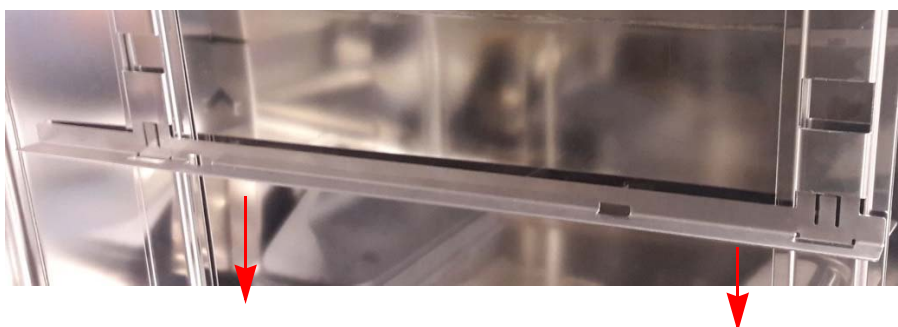
- The positioning of the 4 guiding rails are important. There are 2 left and 2 right side guiding rails. The longer end of the guiding rail should show to the rear side of the chamber.
- Fit in the guiding rail in an angle, in the picture below a right side guiding rail is shown as an example.



- Lift it upwards.



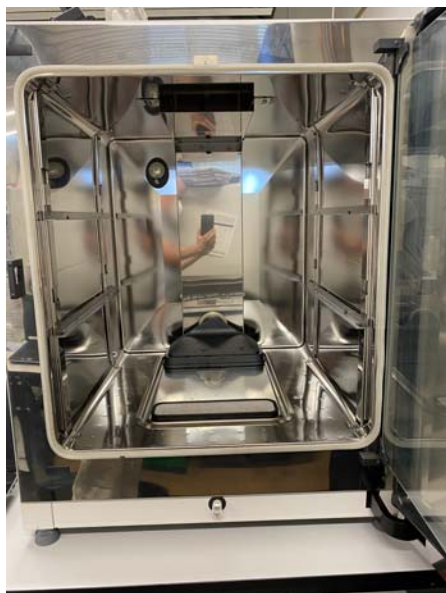
- Press it downwards.



- The tongue/latch must snap in.



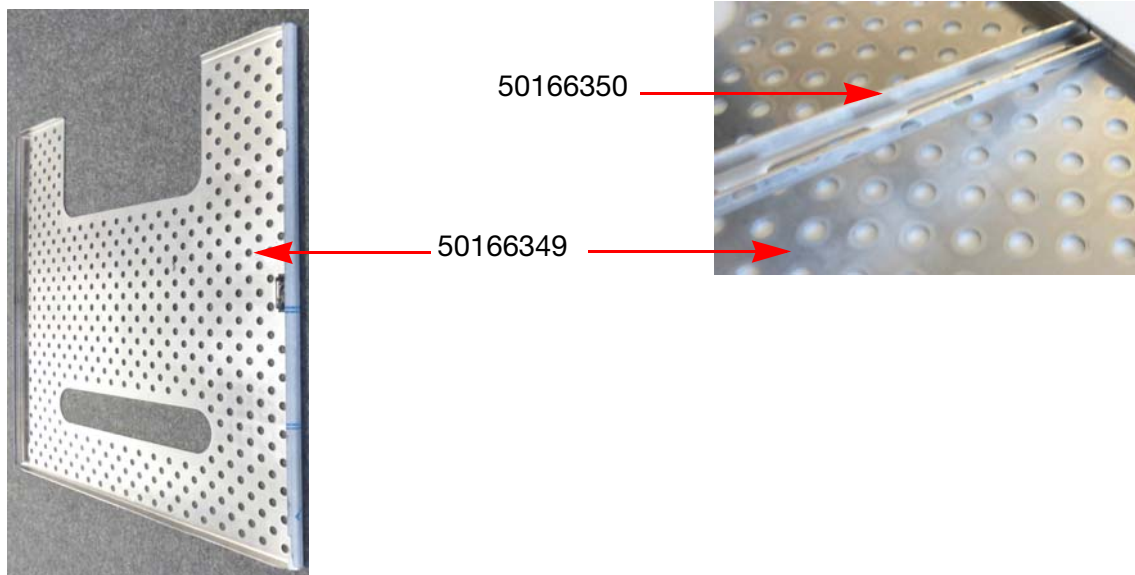
- Repeat the above step for the other 3 guiding rails.



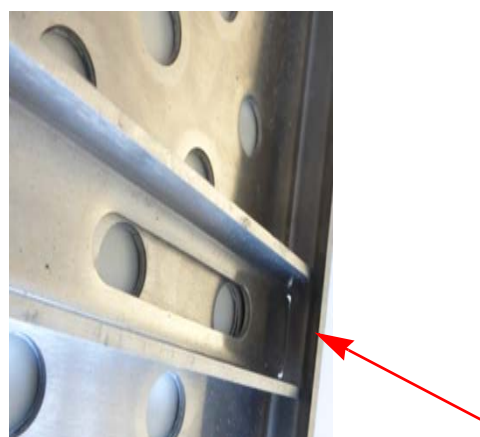
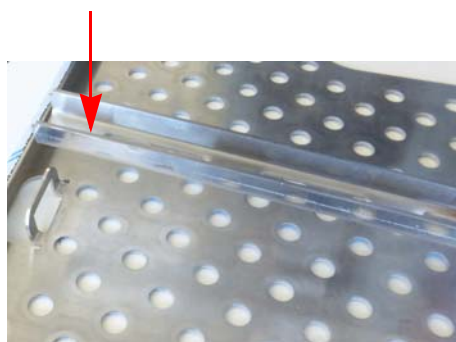
The inner chamber with installed guiding rails.

**Preparing the bottom shelf**

- The bottom shelf has a reinforcement.
- Install the reinforcement into one slot.

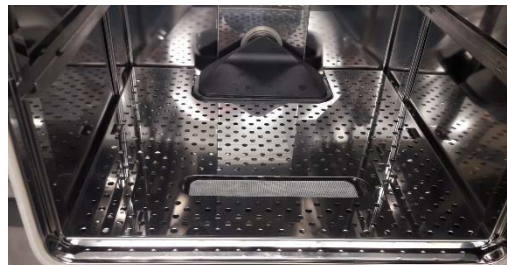
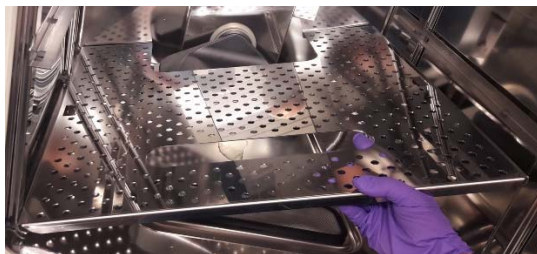


- Press the reinforcement into the other slot, so it snaps in.



### Inserting the shelves

- Use the bottom shelf for the lowest shelf position and 2 split shelves for upper shelf positions.
- Insert the bottom shelf in an angle, so the shelf fits in.

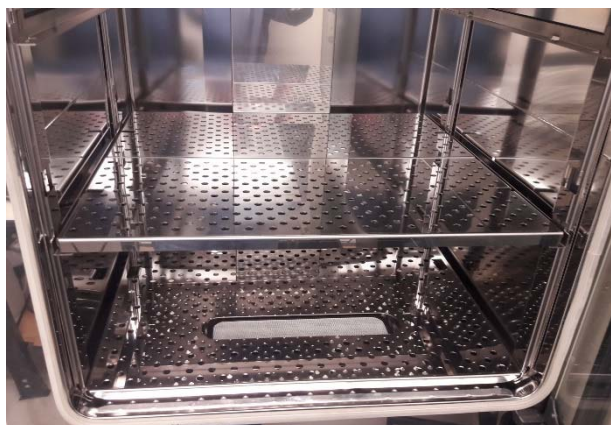
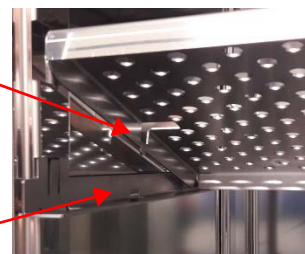


- Insert the split shelves in the middle guide rails, so the pin fits in the slot.



pin

slot



- Insert the split shelves in the upper guiding rails.





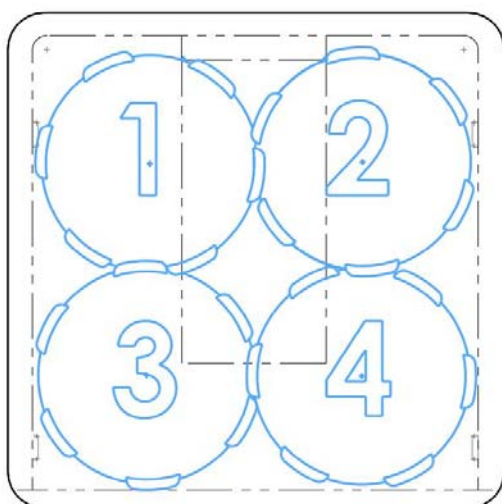
The fully installed shelves.

### Loading the G-Rex Container

- At each of the splitted shelves 2 container can be loaded. At the bottom shelf 2 container.
- The lugs of the container's must fit in ones each other.

**Note** The order to place the G-Rex container is very important!

- Load the G-Rex container beginning at position 1 at the rear side of the inner chamber.
- Then position G-Rex container number 2, so the flap is beside the air duct.
- Load G-Rex container 3 and 4, so the flaps match as shown in the picture below.





The fully loaded incubator.

**Note**

Due to the limited space, the loading of the unit with 10 vessels need to be done properly according the arrangement shown in the pictures. Even if you load according to the picture, it might be needed to fine-tune the arrangement to close the glass door. The containers should not touch the glass door. Please close the glass door in the first step very carefully to check if the arrangement fits and the glass door can be closed and secured by the glass door lock.

## Operation

**Note**

Shelves are not designed to be pulled out.

Containers must be taken out individually.



## Performance Hints

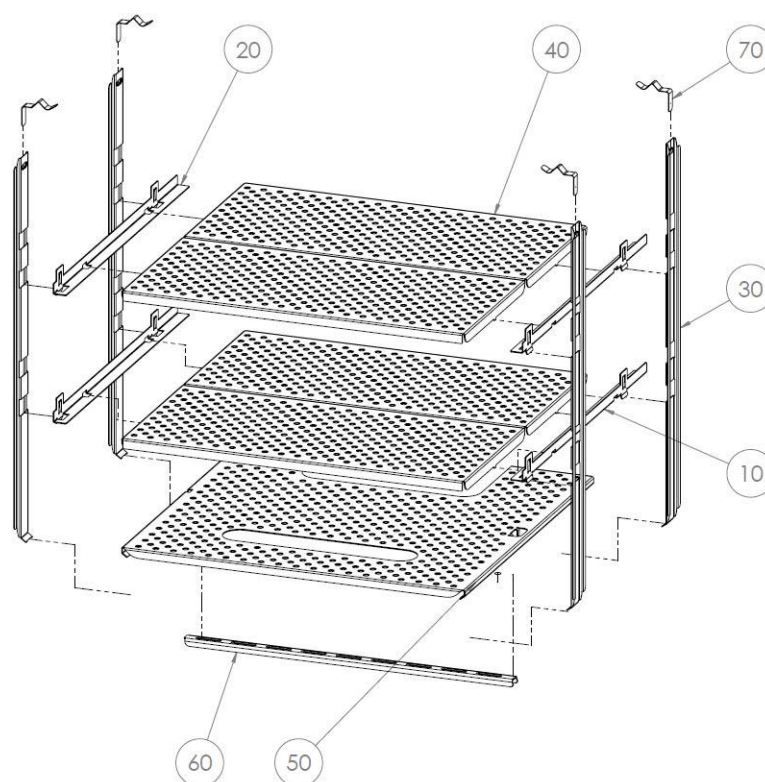
**Note**

Use only pre-heated containers with media to avoid long process stabilization time.

The temperature of the preheated container with medium must not be above the set incubator temperature, otherwise the usable space temperature may overshoot and a temperature alarm may occur.

The customer must verify his application process himself.

## Parts overview



Pos.	Part-No.	Description
10	50166345	Support bar right VIOS
20	50166346	Support bar left VIOS
30	50166347	Supporting profile G-Rex VIOS
40	50166348	Small shelf G-Rex VIOS
50	50166349	Base shelf bottom G-Rex VIOS
60	50166350	Reinforcing strut VIOS
70	50166351	Spring for supporting profile VIOS

## Technical Data

<b>Shelf kit G-Rex VIOS</b>	<b>Unit</b>	<b>Value</b>
Maximum load capacity small shelf G-Rex VIOS, equal distribution on the shelf	kg / lbs	13 / 28.7
Maximum load capacity layer, equal distribution on the shelf	kg / lbs	26 / 57.4
Maximum load capacity base shelf bottom G-Rex VIOS	kg / lbs	13 / 28.7
Maximum overall device loading capacity, equal distribution on the shelves	kg / lbs	65 / 143



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

**thermo**scientific