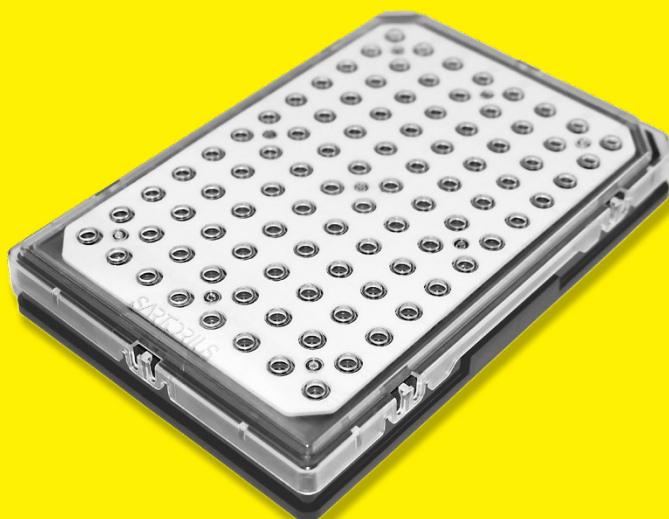


Instructions for Use

# Octet® AE Microplate Evaporation Cover

Cover to Minimize Sample Evaporation in Microplates  
For Octet® BLI systems with flat microplate sample stages



1000147423



SARTORIUS



# Contents

<b>1</b>	<b>About These Instructions.....</b>	<b>4</b>
1.1	Validity.....	4
1.2	Related Documents.....	4
1.3	Symbols Used.....	4
<b>2</b>	<b>Safety Instructions .....</b>	<b>4</b>
2.1	Intended Use .....	4
<b>3</b>	<b>Product Description.....</b>	<b>4</b>
3.1	Product Overview.....	4
3.2	Filling trough.....	4
3.3	Hydration .....	4
<b>4</b>	<b>Process Preparation .....</b>	<b>5</b>
4.1	Scope of Delivery .....	5
4.2	Unpacking .....	5
4.3	Hydrating the Product .....	5
<b>5</b>	<b>Usage .....</b>	<b>5</b>
5.1	Placing the Product on a Microplate.....	5
<b>6</b>	<b>Disposal .....</b>	<b>5</b>
<b>7</b>	<b>Technical Data.....</b>	<b>6</b>
7.1	Dimensions and Weights.....	6
7.2	Storage Conditions .....	6
7.3	Devices and Microplates.....	6
7.4	Liquids .....	6
7.5	Temperatures .....	6
7.6	Average Sample Volume Remaining .....	6
<b>8</b>	<b>Trademark Information.....</b>	<b>6</b>

# 1 About These Instructions

## 1.1 Validity

These instructions are part of the product; they must be read in full. These instructions apply to the following versions of the product:

Product	Item number
Octet® AE microplate evaporation cover	PN 18-5152

## 1.2 Related Documents

- In addition to these instructions, observe the following documents: Octet® BLI Discovery User Guide for the appropriate Octet® BLI system, e.g. Octet® R8e

## 1.3 Symbols Used

- Required action: Describes activities that must be carried out. The activities in the sequence must be carried out in succession.
- ▷ Result: Describes the result of the activities carried out.

# 2 Safety Instructions

## 2.1 Intended Use

The product is intended to reduce sample evaporation in biolayer interferometry experiments (BLI experiments). The product must be hydrated before use.

The product is intended for pharmaceutical and biotechnological research. The product is **not** suitable for diagnostic procedures, therapeutic procedures, or patient-related applications.

The product must only be used with approved analysis devices (for suitability, see chapter “7.3 Devices and Microplates”, page 6).

The product is designed for single-use and must be disposed of after use.

The product is intended exclusively for use in accordance with these instructions. Any other use is considered **improper**.

### Operating conditions for the product

Only use the product indoors. The product may only be used with the equipment and under the operating conditions described in the Technical Data section of these instructions.

# 3 Product Description

## 3.1 Product Overview

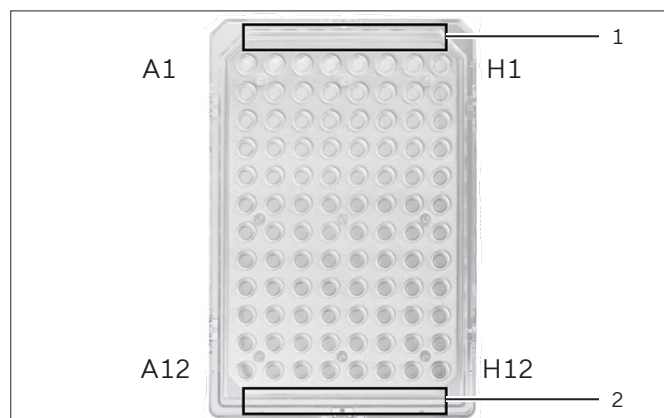


Fig. 1: Evaporation cover (back)

Pos.	Name
1	Top filling trough
2	Bottom filling trough

## 3.2 Filling trough

The filling trough allows liquid to flow into the absorbent layer.

## 3.3 Hydration

The hydration of the absorbent layer creates a saturated vapor layer over the well, which ensures that the sample concentration remains constant.

## 4 Process Preparation

### 4.1 Scope of Delivery

Item	Quantity
Octet® AE microplate evaporation cover	10

### 4.2 Unpacking

#### Procedure

- ▶ If the product is stored before use: Observe the storage conditions (see chapter “7.2 Storage Conditions”, page 6).
- ▶ Unpack the product.

### 4.3 Hydrating the Product

The deionized water must be pipetted into the top and bottom filling troughs. This will hydrate the absorbent layer.

The pipette tips must be held at a slight angle in the filling trough. This will ensure that the pipette tip openings are **not** blocked by the filling trough walls.

Material:     Recommended for even dispensing:  
                     8-channel pipette, set to 150 µL  
                     Liquid: 1.2 mL of deionized water (8 x  
                                  150 µL) per trough filling; 7.2 mL in total

#### Procedure

- ▶ Place the product on the laboratory bench with the bottom facing upwards. This makes it possible to access the top and bottom filling troughs.
- ▶ When pipetting, proceed as follows:
  - Dispense the liquid slowly so that it does **not** overflow into the adjacent wells.
  - Do **not** tilt the product. The liquid flow must **not** be increased.

- ▶ Repeat the following 2 hydration steps 3 times in total; so that 3.6 mL is dispensed per trough:



- ▶ Pipette 1.2 mL of liquid into the top filling trough.
- ▶ Pipette 1.2 mL of liquid into the bottom filling trough.

- ▶ When the entire 7.2 mL of deionized water has been dispensed: Ensure that the absorbent layer is fully saturated.
- ▶ To remove excess liquid in the top or bottom filling trough: Tilt the product slightly and absorb the excess liquid with a lint-free cloth, e.g., Kimwipes™.

## 5 Usage

### 5.1 Placing the Product on a Microplate

#### Procedure

- ▶ Turn the product over and place it on a prepared microplate.
- ▶ Push down on the product until it engages.
- ▶ The product is ready for assays.

## 6 Disposal

#### Procedure

- ▶ Dispose of the product in accordance with local government regulations.
- ▶ Dispose of the packaging in accordance with local government regulations. Utilize local recycling options.

## 7 Technical Data

### 7.1 Dimensions and Weights

	Unit	Value
Dimensions (L x W x H)	mm	127.8 x 85.9 x 9.8
Weight	g	27

### 7.2 Storage Conditions

Storage at room temperature

Dry, in original packaging

### 7.3 Devices and Microplates

	Unit	Value
Approved devices		
Octet® R8e   RH16   RH96		
Biosensor offset, set to at least	mm	3
Approved microplate		
Greiner 96-well microplate (catalog no. 655209)		

### 7.4 Liquids

Deionized water

**No** organic solvents

### 7.5 Temperatures

	Unit	Value
Temperature in assays	°C	+15 – +40

### 7.6 Average Sample Volume Remaining

	Unit	Value
When using a minimum volume of 200 µL in 96-well microplate, after 16 hours at 25 °C	%	90

## 8 Trademark Information

Kimwipes is a registered trademark of Kimberly-Clark Worldwide, Inc.

Sartorius BioAnalytical Instruments Inc.  
47661 Fremont Blvd.  
Fremont, CA, 94538  
USA

[www.sartorius.com](http://www.sartorius.com)

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote all genders.

Copyright notice:

These instructions, including all components, are protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

Last updated:

01 | 2025

© 2025

Sartorius BioAnalytical Instruments Inc.  
47661 Fremont Blvd.  
Fremont, CA, 94538, USA

NT | Publication No.: WIQ6008-e250102