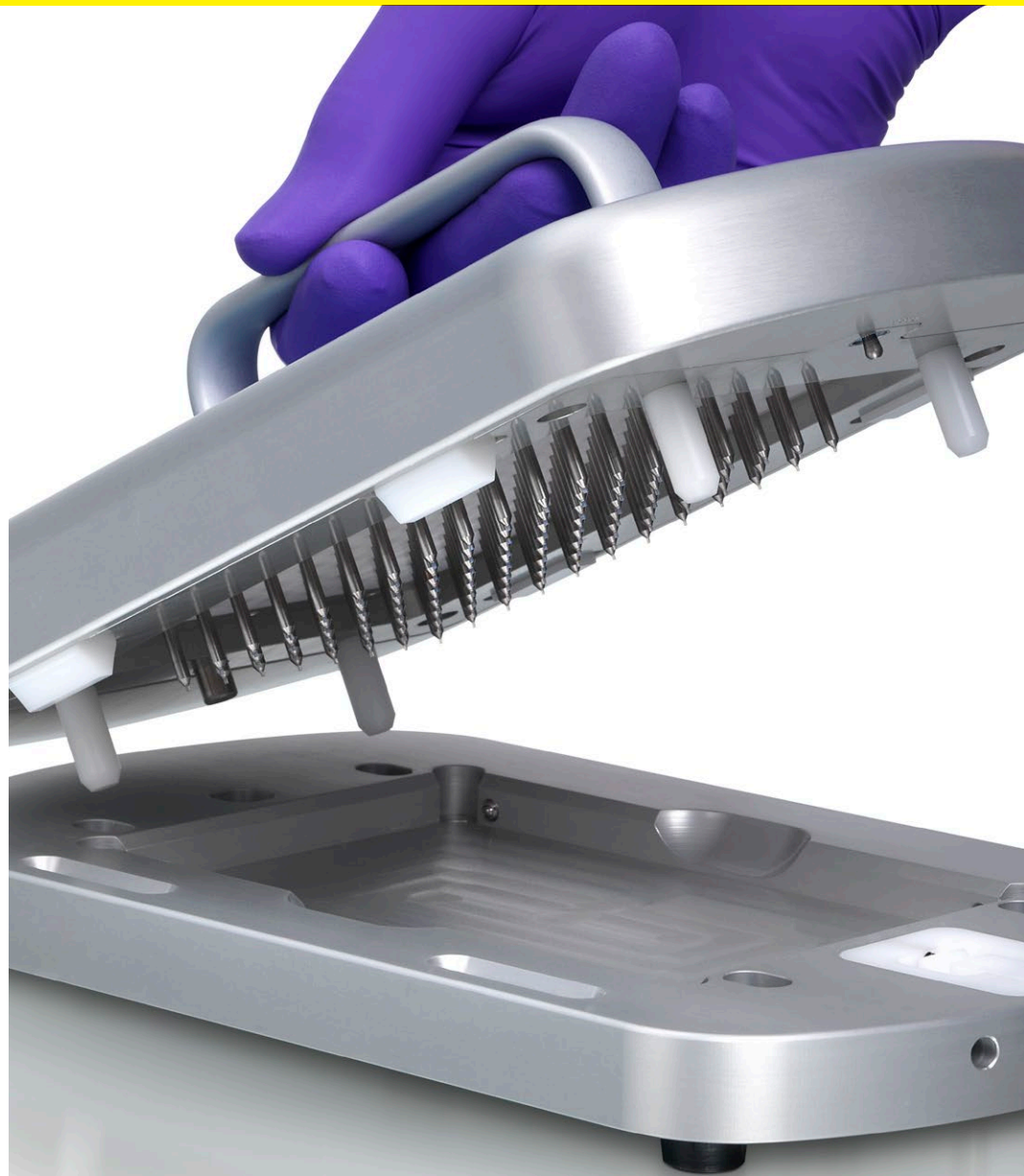


## Incucyte® 96-well Woundmaker Tool Care Instructions

### General Guidelines

- Never place the pin block directly onto the bench surface.
- Never turn the pin block over—this can damage the pins.
- Never use the Woundmaker with a dry or partially dry plate. Make sure each well contains media or PBS before wounding.
- Never use the Woundmaker with more than 100  $\mu$ L of media per well.
- Do not push down on the pin block when making wounds.
- Store the Woundmaker in a clean, dust-free and safe location, such as a biological safety cabinet, when not in use.
- Clean the Woundmaker according to the wash protocol after every use.
- Never put more than 45 mL of any solution into the wash boat. This will prevent damage to the pin mechanism.



# Incucyte® 96-well Woundmaker Tool Care Instructions

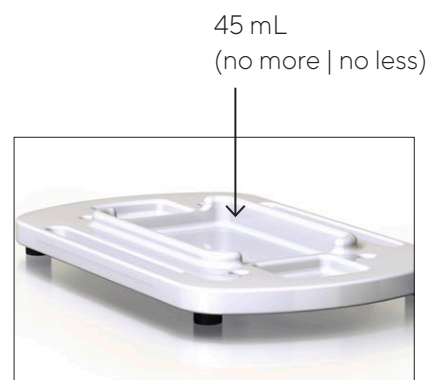


## Wash Instructions

1. Prior to use after storage (5 minute washes)
  - Sterile distilled water (45 mL)
  - 70% ethanol (45 mL)
2. Between plates of the same cell line (5 minute washes)
  - Sterile distilled water (45 mL)
3. After last experiment of the day or between plates containing different cells lines (5 minute washes)

Note: Do not use bleach when cleaning the Woundmaker as it will damage the tool, causing issues creating accurate wounds.

  - 0.5% Alconox (45 mL)
  - Sterile distilled water (45 mL)
  - 70% ethanol (45 mL)



## Sales and Service Contacts

For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)

### Essen BioScience, A Sartorius Company

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

E-Mail: [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)

**North America**  
Essen BioScience Inc.  
300 West Morgan Road  
Ann Arbor, Michigan, 48108  
USA  
Telephone +1 734 769 1600  
E-Mail:  
[orders.US07@sartorius.com](mailto:orders.US07@sartorius.com)

**Europe**  
Essen BioScience Ltd.  
Units 2 & 3 The Quadrant  
Newark Close  
Royston Hertfordshire  
SG8 5HL  
United Kingdom  
Telephone +44 1763 227400  
E-Mail:  
[euorders.UK03@sartorius.com](mailto:euorders.UK03@sartorius.com)

**APAC**  
Sartorius Japan K.K.  
4th Floor Daiwa Shinagawa North  
Bldg.  
1-8-11 Kita-Shinagawa  
Shinagawa-ku, Tokyo  
140-0001  
Japan  
Telephone: +81 3 6478 5202  
E-Mail: [orders.US07@sartorius.com](mailto:orders.US07@sartorius.com)

Specifications subject to change without notice.

© 2020. All rights reserved. Incucyte, Essen BioScience, and all names of Essen BioScience products are registered trademarks and the property of Essen BioScience unless otherwise specified. Essen BioScience is a Sartorius Company. Publication No.: 8000-0126-D00

Status: 09|2020