

# iQue® Enhanced Cartridge Kit

For use in iQue®3



## Product Information

iQue® Enhanced Cartridge Kit includes one each of the standard cartridges. The 4-Pack contains: (1) iQue® QSol cartridge, (1) Flush cartridge, (1) Clean cartridge, and (1) Rinse cartridge.

### List of Available Products

Product Name	Cat. No.	Format
iQue® Enhanced Cartridge Kit	97000	1 Pack
iQue® Enhanced Cartridge Kit	97001	4 Pack
iQue® Enhanced Cartridge Kit	97002	8 Pack

**Table 1.** List of Available Products

### Volume

Each cartridge contains 37 mL, sufficient to run >100 384-well plates. Run time is dependent on protocol.

### Storage

Store iQue® QSol cartridge at room temperature. Once opened, store at 2–8°C. Store Rinse, Clean and Flush cartridges at room temperature.

### Notes

Enhanced Cartridge Kits are only compatible on iQue®3 platforms running iQue Forecyt® versions 7.1 or higher. Cartridges expire 30 days after opening.

## Procedure

### 1. Installing Cartridges on iQue®3 Platform

- 1.1 DO NOT shake the cartridges. Open the blue seal before installing the cartridge in the station.
- 1.2 Open the seal by pulling up on the blue tab. Do not use a tool.
- 1.3 Install the cartridge in the appropriate color-coded slot.

### 2. Uninstalling and Storing Cartridges

- 2.1 Remove the cartridge from the enhanced rinse station.
- 2.2 Close the cartridge after removing it from the station by pressing the seal closed. Store the iQue® QSol at 2–8°C, the rest of the cartridges (Rinse, Clean and Flush) are stored at room temperature. To reuse the iQue® QSol cartridge, remove it from 2–8°C and allow it to equilibrate to room temperature for at least 20 minutes before starting the next experiment.

## Sales and Service Contacts

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

**Sartorius BioAnalytical Instruments, Inc.**

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

© 2021. All Rights Reserved.  
Sartorius BioAnalytical Instruments, Inc.  
iQue® is a Sartorius brand. Intellicyt®, iQue®3, iQue® Screener PLUS, iQue Forecyt®, iQue QBead® and all names of iQue® products are registered trademarks and the property of Sartorius unless otherwise specified.  
Publication No. #17092 Rev B.

#### North America

Sartorius Corporation  
300 West Morgan Road  
Ann Arbor, Michigan, 48108  
Telephone +1 734 769 1600  
E-Mail: [AskAScientist@sartorius.com](mailto:AskAScientist@sartorius.com)  
Online Store: [shop.intellicyt.com](http://shop.intellicyt.com)

#### Europe

Sartorius UK Ltd.  
Units 2 & 3 The Quadrant  
Newark Close  
Royston Hertfordshire  
SG8 5HL  
United Kingdom  
Telephone +44 (0) 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)