

iQue® Enhanced Buffer Cartridge

For use in iQue®3



Product Information

iQue® Enhanced Buffer Cartridge formula is PBS + 0.1% BSA. It is used to coat the sampling line to promote smooth flow of sample.

List of Available Products

Product Name	Cat. No.	Format
iQue® Enhanced Buffer Cartridge	97003	1 Pack
iQue® Enhanced Buffer Cartridge	97004	4 Pack
iQue® Enhanced Buffer Cartridge	97005	8 Pack

Table 1. List of Available Products

Volume

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

Storage

Store at room temperature.
Once opened, store at 2–8°C.

Notes

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

Procedure

1. Installing Cartridge on iQue®3 Platform

- 1.1 DO NOT shake the cartridge. Open the blue seal before installing the cartridges 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 at 2–8°C. To reuse the 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

Sartorius BioAnalytical Instruments, Inc.

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. #17087 Rev B.

North America

Sartorius Corporation
300 West Morgan Road
Ann Arbor, Michigan, 48108
Telephone +1 734 769 1600
E-Mail: AskAScientist@sartorius.com
Online Store: 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

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