

iQue® S1 Fluid Station QSol Buffer Cartridge

Product Information

iQue® QSol buffer is designed to reduce sample build-up and improve sampling on iQue® platforms, including the iQue® Screener, iQue® Screener HD, and iQue® Screener PLUS. Reported benefits of the enhanced buffer include:

- Flexibility in sample selection by reducing carryover and protein buildup
- Better data resulting from a more robust/ more consistent sampling
- Faster time to results from a streamlined workflow
- Convenient, ready to use cartridges and cost saving concentrate.

The buffered solution is biologically compatible and contains a nonionic, non-denaturing surfactant that solubilizes extracellular proteins. The buffer has been tested and validated for use with both cell and bead based assays. It is particularly useful for difficult samples such as:

- Primary cells (PBMCs, bone marrow)
- Yeast cells (sugar in media)
- Long incubation assays (3-7 days)
- High density samples

Kit Contents

Product Name	Cat. No.	Format
iQue® S1 Fluid Station QSol Buffer Cartridge	90283	1 Pack
iQue® S1 Fluid Station QSol Buffer Cartridge	90287	10 Pack
P200 Pipet Tip for Puncturing Cartridge	-	1 Tip

Table 1: Kit Contents

Packaging

Each cartridge contains 29 mL, sufficient to run >25 384-well plates using standard protocol. Alternate protocols could achieve more or fewer plates per cartridge.

Storage

Store at room temperature. Buffer cartridges are manufactured under aseptic conditions. Once opened, store at 2-8°C. Protect from prolonged exposure to light. Do not freeze.

Procedure

1. Opening and Installing Cartridges on your iQue® Platform.

- 1.1 Before use, gently rock cartridge back and forth several times to ensure a homogenous buffer solution. Do NOT shake the cartridge as this will cause the solution to foam.
- 1.2 To open a cartridge, puncture the top foil seal at the indicated location. Use of a standard laboratory pipet tip (P200 size or larger) will aid in the puncture and expansion of the opening to the indicated size.
- 1.3 Buffer cartridges must always be installed on the “Buffer” location on the fluid station. Gently place the opened cartridge into the slot and ensure that the cartridge is properly seated.



Figure 1: Buffer should always be installed at the first, left most position.

2.0 Prime the Installed Probe and Tubing in Buffer.

- 2.1 After initial installation of a cartridge, prime the installed probe and tubing with buffer for several minutes to ensure buffer is coated throughout the sample path. This can be done from the controller window, using the “Prime” function. The system should be primed for at least 5 minutes.
- 2.2 Once the probe and tubing are primed, proceed with normal experiment setup and acquisition.

Note: this step should be repeated every time a new probe and tubing assembly is installed on the system.

3.0 Uninstalling and Storing Buffer Cartridges

- 3.1 Remove the cartridge from the rinse station.
- 3.2 Cover the opening with the provided sticker. Store at 2–8°C, taking precautions to prevent evaporation.
- 3.3 To reuse the cartridge, remove it from 2–8°C and allow it to equilibrate at room temperature for at least 20 minutes before starting next experiment. Shake well before each usage, taking precautions to avoid spilling from the opening.

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.
Status: 08|2022

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