

# iQue® QSol Buffer 100x and 1000x Concentrate

## Product Information

iQue® QSol Buffer is designed to reduce sample build-up and improve sampling on iQue® Platform. 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 a concentrate for custom usage in assays.

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

The 100x Concentrate formulation provides users with the flexibility to generate different quantities of buffer for use in assays. Additionally, this 100x and 1000x concentrate can be directly diluted in sample buffer or assay media and used in an assay, such as by supplementing microwells with the 1000x iQue® QSol.

## Product

Product Name	Cat. No.	Volume	Format
iQue® QSol 100x Concentrate	91304	5mL	1 vial
iQue® QSol 100x Concentrate	91161	25mL	1 vial
iQue® QSol 1000x Concentrate	91363	5mL	1 vial

## Packaging

Cat. No.	Volume	Sufficient to prepare
91304	5mL	500mL
91161	25mL	2.5L
91363	5mL	not apply

## Storage

Store at room temperature. Protect from prolonged exposure to light. Do not freeze. Once iQue® QSol is diluted, store for 2 months at room temperature.

## Procedure

### 1. Prepare 1x Working Solution.

1.1 Before use, gently hand mix or vortex the reagents to ensure a homogenous solution.

### 2.0 Determine how much volume of solution you are trying to prepare, and dilute concentrate by 100x or 1000x to achieve a 1x working solution.

2.2 If preparing 1x working solution, dilute iQue® QSol concentrate into a solution containing PBS + 0.1% BSA.

2.3 If preparing 1x working solution of iQue® QSol in sample buffer or assay media – prepare buffer as normal and add appropriate volume of iQue® QSol concentrate.

## 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)

© 2023. 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: 10|2023

#### North America

Sartorius Corporation  
300 West Morgan Road  
Ann Arbor, Michigan, 48108  
Telephone +1 734 769 1600  
E-Mail: [LPS.OPM.NA@Sartorius.com](mailto:LPS.OPM.NA@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)