

iQue[®] 6-Peak Validation Beads

Product Information

This product contains 2 vials of a mixture of 3.0–3.4 μm beads that are excited by the instrument's red laser and are detected in the R/Red (FL4) channel. Unlabeled beads are in the vial labeled Peak 1 and a mixed population of beads exhibiting 5 different fluorescent intensities are in the vial labeled Peaks 2–6.

Product

6-Peak Validation Beads (R/Red)
Cat. Number: 90295

Packaging

2 x 5 mL (1 vial for peak 1; 1 vial for peaks 2–6) and a USB stick with the bead template on it.

Concentration

1×10^7 particles/mL

Storage Buffer

Deionized water with 0.02% Sodium Azide and 0.01% NP40

Storage

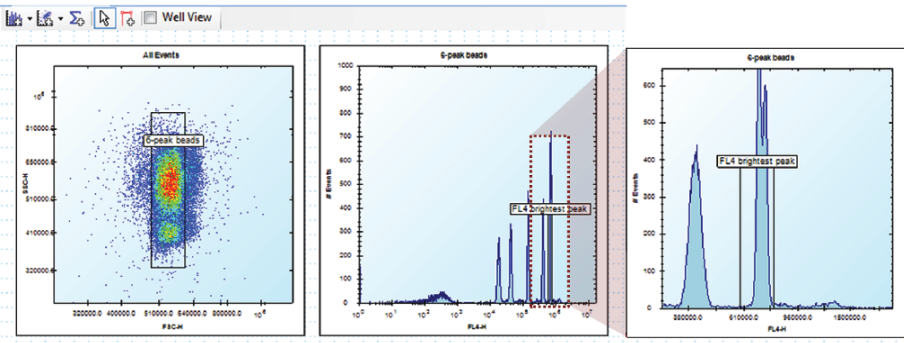
If unopened, stable at room temperature for 3 years. Store at 2–8°C after opening; expires 1 year after opening. Do not freeze. Protect from light.

Use

For daily validation of the instrument performance. Dilute 4 drops from the Peak 1 vial and 4 drops from the Peaks 2–6 vial into a single tube containing 500 μL deionized, 0.22 μm filtered water. Perform validation as described in the instrument's Hardware Manual.

Notes

Shake vigorously or vortex before each use. The diluted bead solution can be stored at 2–8°C for repeated use up to 1 week. Discard solution after 1 week.



Procedure

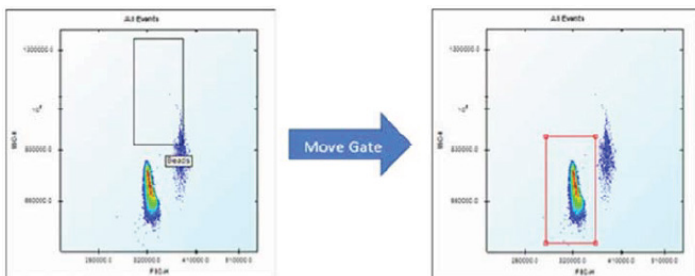
1. Install the Bead Lot Template (iQue Forecyt® 5.1 or newer)

- 1.1 Launch iQue Forecyt®.
- 1.2 Insert the enclosed USB stick.
- 1.3 Click File Import Templates.
- 1.4 Click the Add...button on the right side of the screen.
- 1.5 Browse the USB directory and select the .IET file that applies to your instrument.
- 1.6 Click OK. The default template is now installed.

2.0 Customize the Template for your Instrument (iQue Forecyt® 5.1 or newer)

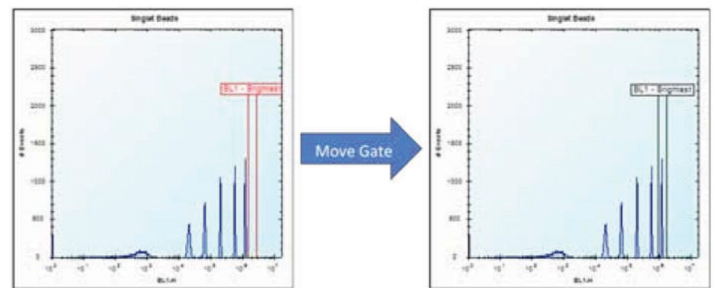
Prior to first use, optimize the template for the performance characteristics of your specific instrument.

- 2.1 Click File New Experiment.
- 2.2 Choose, Use Template and Browse...select the template that came with these beads.
- 2.3 Give the experiement a name and click Save.
- 2.4 Place the preparation of the Validation Beads on the detector inlet.
- 2.5 Click the Run button in the Controller Console and wait for the beads to be acquired.
- 2.6 Move the Bead gate to encompass the primary bead population (see insert below).

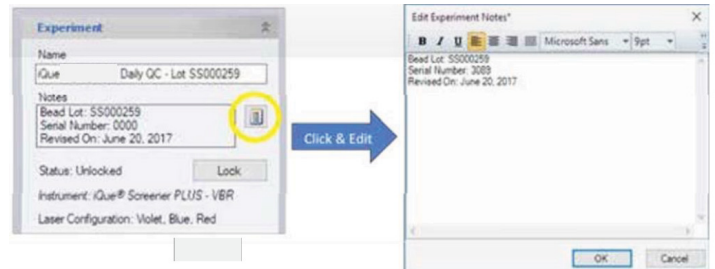


- 2.7 Move the "FL4 - Brightest" gate to encompass the top (right-most) peak and the FL4 histogram.

Note: Do not resize the gate. Just click and drag the label to move it sideways.



- 2.8 Switch to the Design tab. Click on the Notepad icon next to the Notes field. Edit the notes to include your instrument's serial number and the date you revised the template and click OK.



- 2.9 Click save.
- 2.10 Click File Save as Template. Give your template a new name and click save. Your cusotmized template is now ready to use for future Daily QC runs.

Metrics

1. Go to metrics tab and review 6 peak tube experiment results.
2. Make sure all boxes read 1 (pass). If a 0 appears (fail) refer to the instrument hardware manual for detailed trouble shooting procedures.

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. #10581 Rev E.

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