

iQue® 3 & iQue® Screener PLUS Validation Beads

Note: To set up your Validation Beads, please follow these instructions to install and customize the template. Also, the lot number shown in the template examples may be different than your lot number of Validation Beads. Please insert the lot number from the beads you received into the templates.

Product Information

Kit Contents

- One vial of Validation Bead stock
- USB Stick with template (.IET or .BLT) files

Applies To

- iQue® 3 & iQue® Screener PLUS family of instruments

Software Requirements

- iQue Forecyt® Software 5.4 R1 or Newer

Procedure

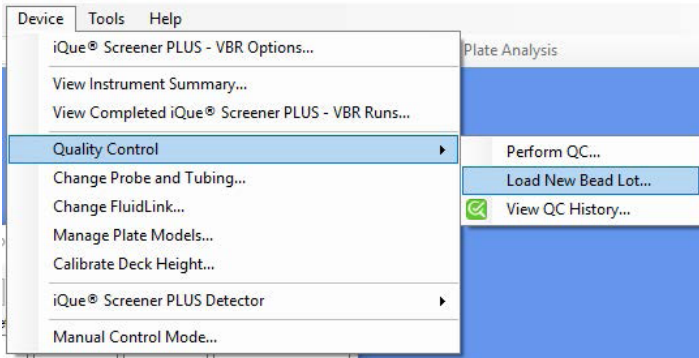
1.0 Prepare the Beads

- 1.1 Add 300 μ L of iQue® Screener Buffer or iQue® QSol Buffer to a 1.5 mL microcentrifuge tube.
- 1.2 Vortex the vial of Validation Beads.
- 1.3 Add two drops of Validations Beads stock to the buffer.
- 1.4 Vortex for 5 seconds.

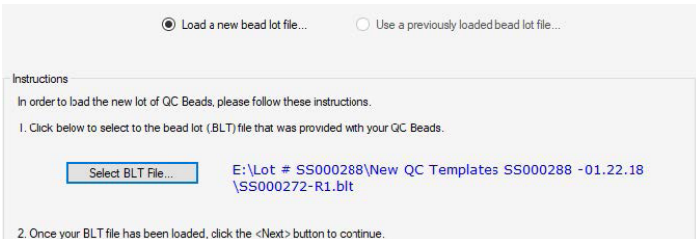
2.0 Install the Bead Lot Template (iQue Forecyt® 6.2 and newer)

The following screenshots demonstrate the workow to update your validation bead lot template.

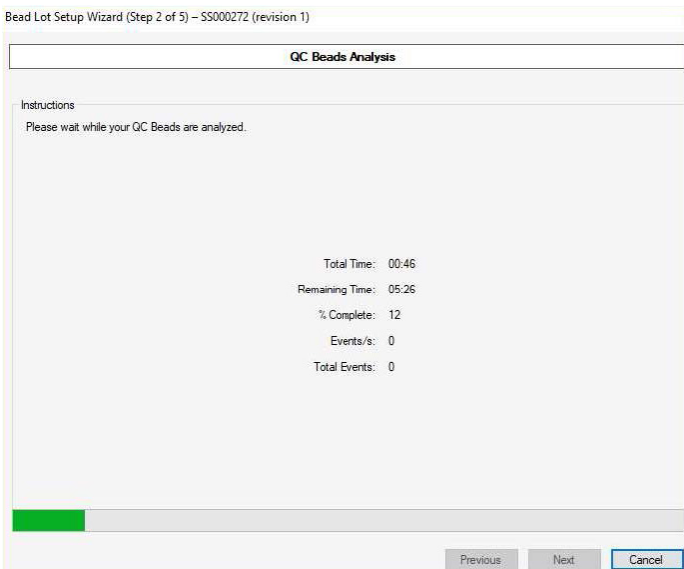
- 2.1 Launch iQue Forecyt® software.
- 2.2 Insert the enclosed USB stick.
- 2.3 Click Device | Quality Control | Load New Bead Lot...



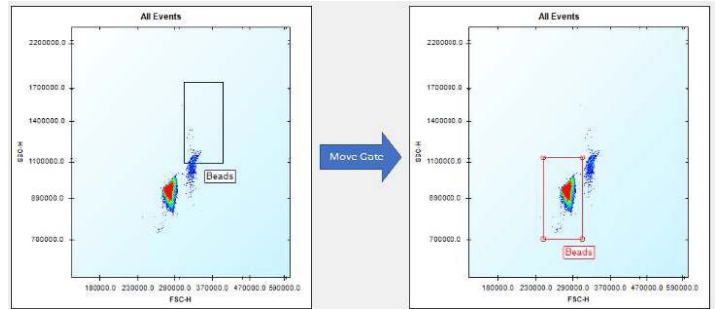
- 2.4 Click the Select .BLT File... button in the Bead Lot Setup Wizard.
- 2.5 Browse the USB directory and select the .BLT file.



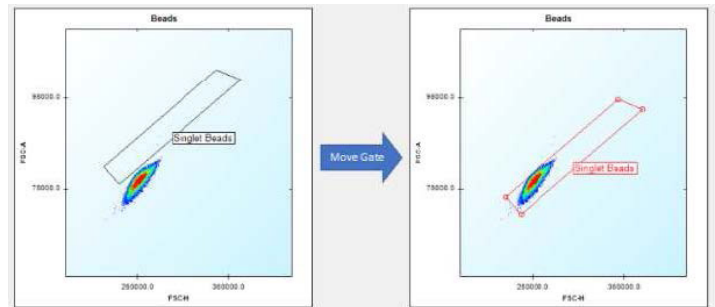
- 2.6 Place the preparation of the Validation Beads in the microcentrifuge tube holder.
- 2.7 Click the Next button to begin QC Beads Analysis.



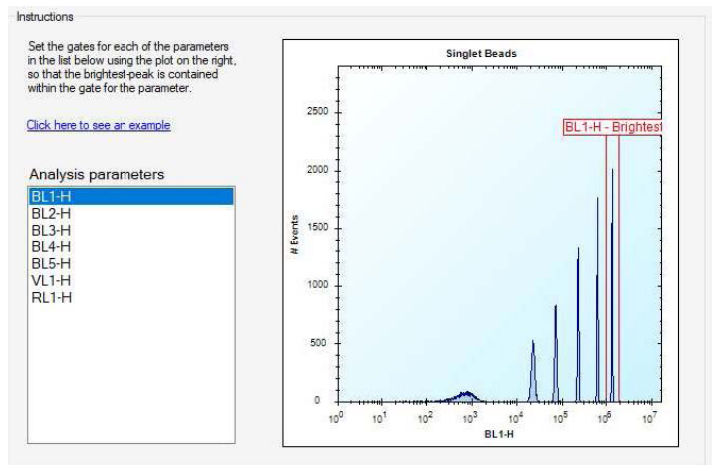
- 2.8 Move the Bead Gate to encompass the primary bead population then press the Next button.



- 2.9 Move the Singlet Bead gate to encompass the singlet bead population then press the Next button.

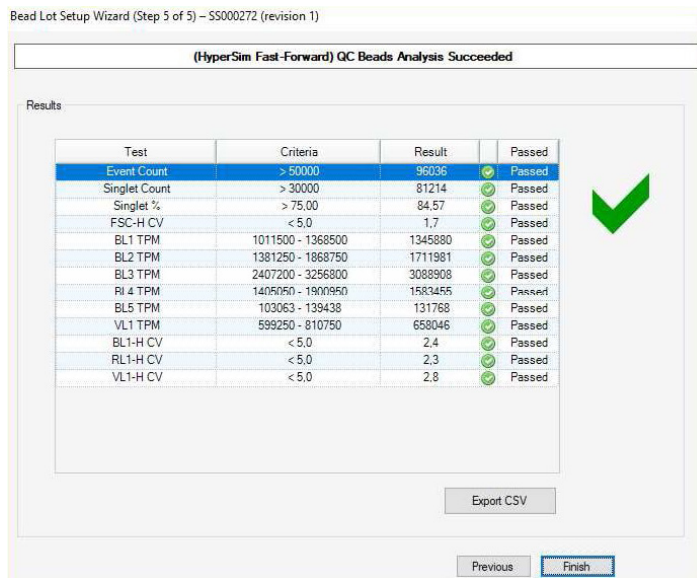


- 2.10 Move the "BL1-H - Brightest" gate to encompass the top (right-most) peak on the histogram and press the Next button.



- 2.11 Repeat Step 2.10 for the other Analysis parameters.

2.12 Click Finish to complete the Bead Lot Setup for the instrument.



3.0 Install the Bead Lot Template (iQue Forecyt® 6.1 and earlier)

1. Launch iQue Forecyt® software.
2. Insert the enclosed USB stick.
3. Click File | Import Templates.
4. Click the Add... button on the right side of the screen.
5. Browse the USB directory and select the .IET file that applies to your instrument.
6. Click OK. The default template is now installed.

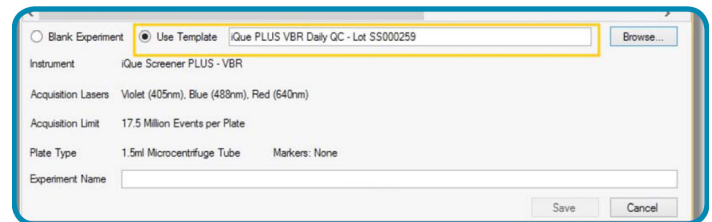


Name	Date modified	Type	Size
<input type="checkbox"/> iQue PLUS BR Daily QC - Lot SS000259.i	6/23/2017 2:26 PM	IET File	9 KB
<input type="checkbox"/> iQue PLUS VBR Daily QC - Lot SS000259.i	6/23/2017 2:25 PM	IET File	10 KB
<input type="checkbox"/> iQue PLUS VYB Daily QC - Lot SS000259.i	6/23/2017 2:25 PM	IET File	10 KB

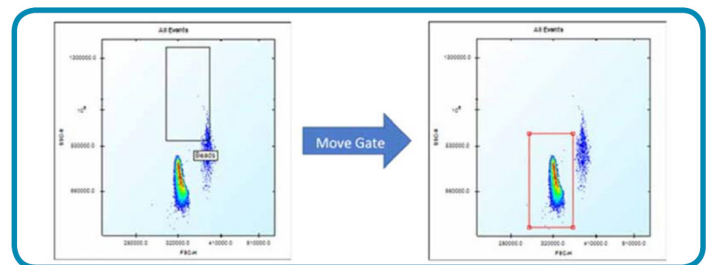
4.0 Customize the Template for your Instrument (iQue Forecyt® 6.1 and earlier)

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

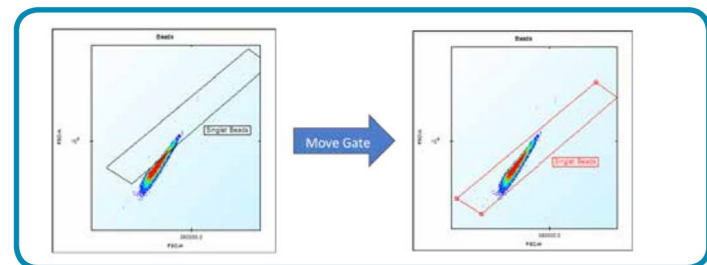
1. Click File | New Experiment.
2. Choose Use Template and Browse...select the template that came with these beads.



3. Give the experiment a name and click Save.
4. Place the preparation of the Validation Beads in the microcentrifuge tube holder.
5. Click the Run button in the Controller Console and wait for the beads to be acquired (approx. 5 min)
6. Move the Bead gate to encompass the primary bead population.

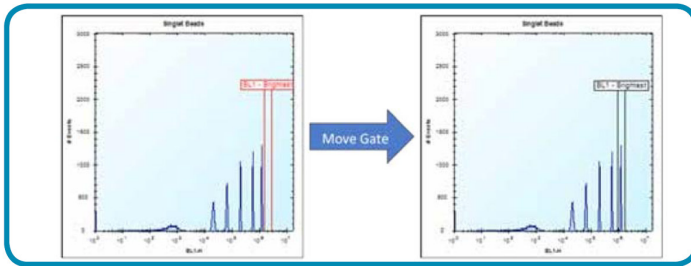


7. Move the Singlet Bead gate to encompass the singlet bead population.

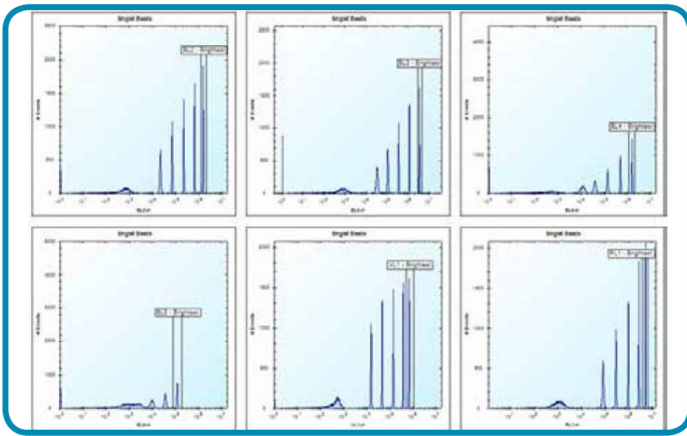


4.8 Move the "BL1-H-Brightest" gate to encompass the top (right-most) peak on the BL1-H histogram.

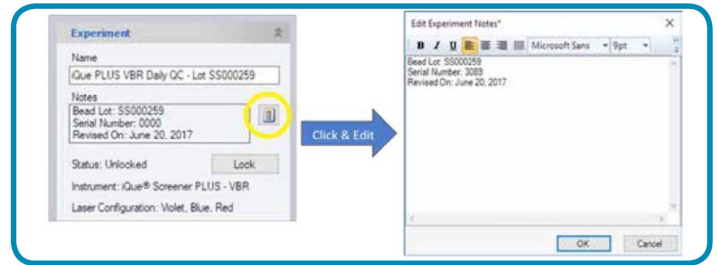
Note: Do not resize the gate. Just click and drag the label to move it sideways.



4.9 Repeat **Step 4.8** for the other six fluorescence histograms: BL2, BL3 BL4, BL5, VL1 and RL1-H.



4.10 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.



4.11 Click Save.

4.12 Click File | Save as Template. Give your template a new name and click Save. Your customized template is now ready to use for future Daily QC runs.

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. #11368 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