

iQue[®] 8-Peak Validation Beads

Product Information

This product contains a mixture of 3.2 µm Rainbow Particles in 8 different fluorescent intensities. Every Rainbow particle contains a mixture of fluorophores that enable their excitation with the blue laser (488 nm) to validate B/Green (FL1), B/Yellow (FL2), and B/Red (FL3) channels of the instrument.

Product

8-Peak Validation Beads (B/Green, B/Yellow, B/Red)
Cat. Number: 90296

Packaging

1-5 mL vial

Concentration

1×10^7 particles/mL

Storage Buffer

Deionized water with 0.02% Sodium Azide
and 0.01% NP40

Storage

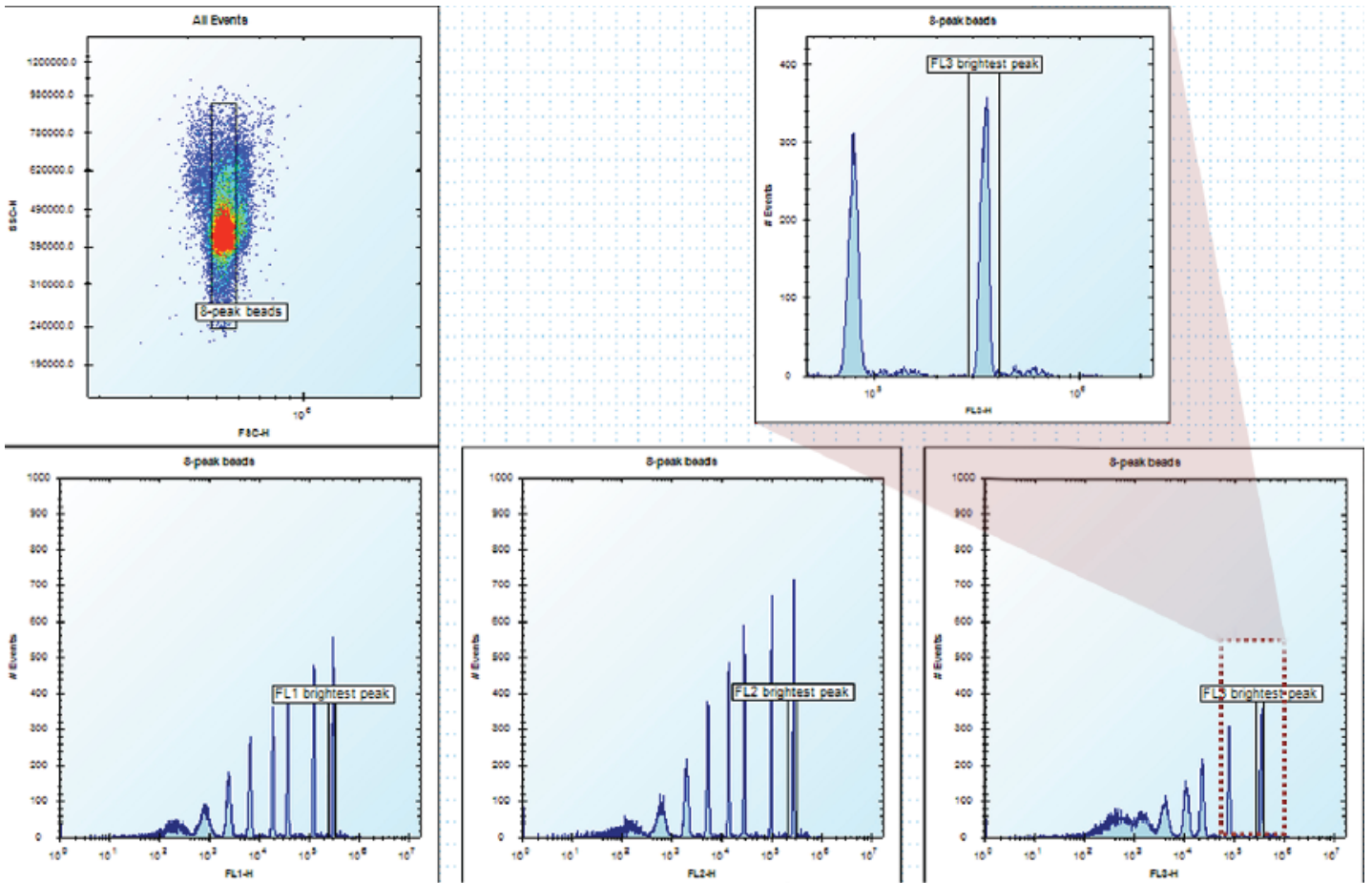
If unopened, stable at room temperature for 3 years.
Store at 2–8°C after opening; expires 1 year after
opening. Do not freeze. Protect from light.

Use

Run daily for instrument performance validation. Dilute
4 drops into 500 µL deionized, 0.22 µm filtered water.
Perform validation as described in the instrument's
User Guide.

Notes

Shake vigorously or vortex before each use. Diluted
bead solution can be stored at 4°C for repeated use.
Preparation of a new stock each week is recommended.



Procedure

1. Set a gate on beads, and collect ~50,000 events.
2. Create 3 histograms and set 1 each for FL1-H, FL2-H, and FL3-H on the x-axis. Draw a gate on the brightest peak in each histogram
3. Create a statistics panel and display the FL-H CV for the brightest peak in each channel. The CV value should be < 5%. If any CV is greater than 5%, the system is out of specification. Refer to instrument's User Guide for detailed troubleshooting 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. #10582 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