

ProLine™ Calibration Beads

Catalog #145-1081

Intended Use

ProLine calibration beads are optimized to verify the alignment performance and determine the sort drop delay on the S3™ cell sorter. Each bottle contains a single population of ready-to-use fluorescent beads.

Contents

3 x 5 ml bottles

Concentration

1 x 10⁶ beads/ml

Bead Size

6 µm

Storage Buffer Composition

Deionized water with 0.02% sodium azide and 0.01% NP-40

Storage

Store at 2–8°C.

Important: Do not freeze. Protect from light.

Bead Lot Specifications

Specifications for each bead lot can be found in the certificate of analysis. Download the certificate of analysis from the Support section in Bio-Rad's website, www.bio-rad.com.

Directions for Use

Note: The supplied concentration will provide the adequate number of beads/µl for use.

Note: If the S3™ ProSort™ software is not open, double click the **ProSort icon**, log in, and select **Start-Up** in the Setup and Maintenance toolbar.

Important: Dilution of these beads is not required.

Preparation of Beads

1. Vortex the 5-ml bottle containing the ProLine beads vigorously.
2. Using the bottle dropper, release ten drops of solution (approximately 0.5 ml) into a 5-ml tube.

Quality Control Procedure

1. Place the 5-ml tube containing the ProLine calibration beads on the loading stage.
2. Move the loading stage to the run position.

3. Click **Run QC** in the Setup and Maintenance toolbar (Figure 1).

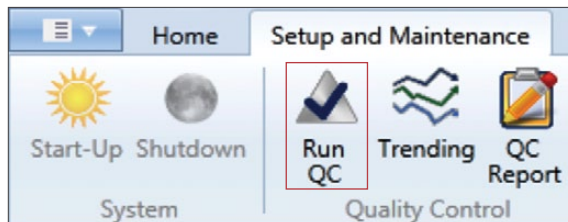


Fig. 1. Setup and Maintenance toolbar with Run QC selected.

4. Select your Bead Lot in the Event Based Alignment Settings window and click the **checkmark** (Figure 2). The software will perform both event-based alignment and drop delay setting.

Note: If your bead lot is not available in the dropdown menu, please see the S3 cell sorter manual at www.bio-rad.com/cellsorter for instructions on uploading a bead lot file.

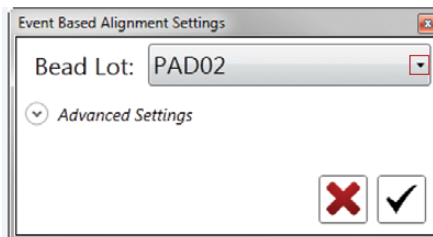


Fig. 2. Event Based Alignment Settings window with bead lot dropdown arrow selected.

5. After the drop delay setting is completed, click the **checkmark** in the Drop Delay Alignment Status window (Figure 3).

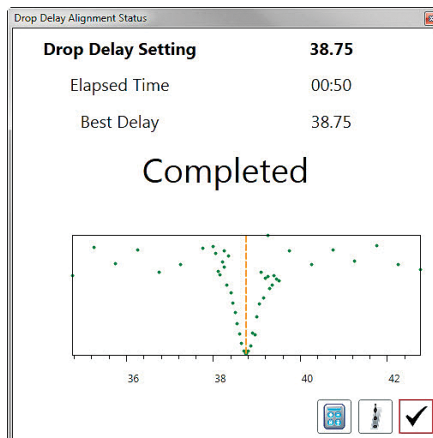


Fig. 3. Drop Delay Alignment Status window with checkmark selected.

6. The QCReportWindow will appear with a passed/failed QC Status (Figure 4).

Passed QC Status: The S3 cell sorter is ready to use. Close the QCReportWindow and start your run.

Failed QC Status: Rerun the quality control procedure up to two more times. If a failure continues, contact your local technical support team at [www. bio-rad.com](http://www.bio-rad.com).

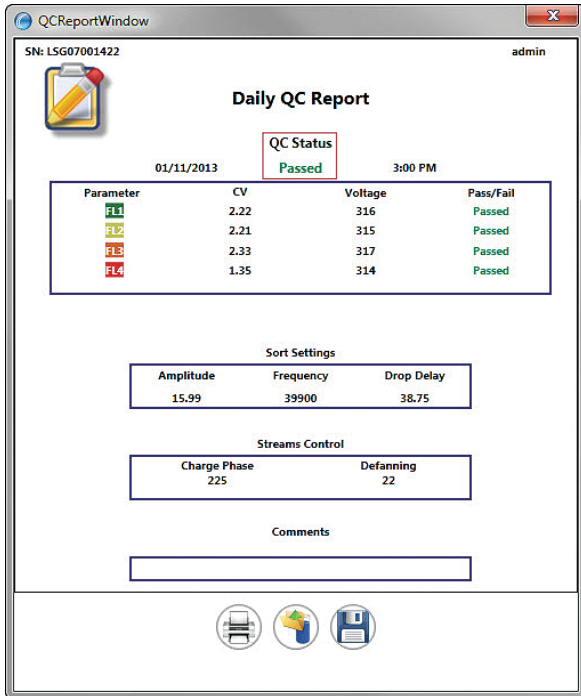


Fig. 4. QCReportWindow with passing QC Status.

Caution: This reagent is for research use only.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000