AlignFlow[™] Flow Cytometry Alignment Beads

Table 1 Contents and storage

Material	Amount	Concentration	Storage	Stability		
AlignFlow [™] beads	2 × 1.5 mL dropper vials	Suspension in water containing 0.05% Tween [®] 20 and 2 mM sodium azide*	 2-8°C Protect from light Sonicate before use Do not freeze 	When stored as directed, reagents are stable for at least 1 year.†		
* The AlignFlow [™] suspensions contain ~1.2 × 10 ⁸ beads/mL (0.1% solids). † No leaching or degradation is expected.						
Approximate fluorescence excitation and emission maxima: See Table 2, page 2.						

Introduction

To help ensure accurate and reproducible results, flow cytometers should be checked daily for proper performance. Reference standards are indispensable for conducting these tests. AlignFlow[™] flow cytometry alignment beads are reliable references for aligning, focusing, and calibrating flow cytometers. These fluorescently stained, polystyrene microspheres are highly uniform with respect to size and fluorescence intensity. The AlignFlow[™] beads are available in three versions: for 350–370-nm excitation with UV lasers, for 488-nm excitation with blue lasers, and for 633-nm excitation with red lasers, and each version is available in two sizes: 2.5 µm–diameter and 6.0 µm–diameter (see Table 2, page 2).

The fluorescent dyes have been carefully selected for optimal excitation by laser sources commonly used in flow cytometry. The 488 nm–excitable beads for blue lasers emit broadly from 515 nm to 660 nm. The UV-excitable beads emit from 400 nm to 470 nm, and the 633 nm–excitable beads for red lasers, from 645 nm to 680 nm. Because the dyes are contained inside each microsphere, instead of merely on the bead surface, AlignFlow[™] beads exhibit superior signal stability.

Using AlignFlow[™], an operator can reproducibly adjust parameters crucial to flow cytometry before an experiment is started. The beads approximate the size, emission wavelength, and intensity of many biological samples and permit the calibration of the flow cytometer's laser source, optics, and stream flow without wasting valuable and sensitive experimental material. Because of their excellent stability, these beads can also serve as a daily reference standard by which an instrument's performance and reliability are evaluated.

Table 2 AlignFlow[™] flow cytometry alignment beads.

Catalog no.	Bead diameter	Ex *	Em range †	Laser type		
A16500	2.5 µm	488 nm	515–660 nm	Blue laser		
A16503	6.0 µm	466 000		Blue laser		
A16501	2.5 µm	633 nm	645–680 nm	Red laser		
A16504	6.0 µm	633 nm				
A16502	2.5 µm	350–370 nm	(00 (70	UV laser		
A16505	6.0 µm	350-370 nm	400–470 nm			
* Fluorescence excitation, in nm, for which the beads were designed. † Approximate useful emission range, in nm.						

Guidelines for Use

Life Technologies' AlignFlow[™] beads serve as a reference standard for calibrating flow cytometers. Experimental protocols depend somewhat on the flow cytometer and software used; please refer also to the reference materials applicable to your particular instrument. Before sampling, be sure that the polystyrene beads are uniformly suspended by vortex mixing and sonicating the suspension. Generally, one drop added to 1 mL of Haema-Line 2 sheath fluid or buffered saline solution provides an appropriate concentration for analysis; mix well before applying the sample.

During the alignment, set the approximate photomultiplier tube (PMT) voltage to place the sample in a convenient range for the flow cytometer, then maximize the signal and minimize the coefficient of variation (CV) by making the appropriate optical adjustments. Follow the instrument manufacturer's manual carefully for optimal performance.

When using the AlignFlow[™] beads as a daily reference source, maximize the peak heights and minimize the CVs, as outlined above. Once the settings are optimized, record all settings and print out the relevant data for your record. Each day, pre-run the flow cytometer using these reference beads and compare the output to the reference record. Deviations from the baseline readings may indicate instrument malfunction.

References

1. Flow Cytometry, A Practical Approach, 2nd Edition, M.G. Ormerod, Ed. IRL Press (1994); 2. Givan A.L., Flow Cytometry, First Principles, Wiley-Liss (1992).

Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
A16500	AlignFlow [™] Flow Cytometry Alignment Beads for Blue Lasers *2.5 µm*	2 × 1.5 mL
A16501	AlignFlow [™] Flow Cytometry Alignment Beads for Red Lasers *2.5 µm*	2 × 1.5 mL
A16502	AlignFlow [™] Flow Cytometry Alignment Beads for UV Lasers *2.5 μm*	2 × 1.5 mL
A16503	AlignFlow [™] Flow Cytometry Alignment Beads for Blue Lasers *6 µm*	2 × 1.5 mL
A16504	AlignFlow [™] Flow Cytometry Alignment Beads for Red Lasers *6 µm*	2 × 1.5 mL
A16505	AlignFlow [™] Flow Cytometry Alignment Beads for UV Lasers *6 µm*	2 × 1.5 mL

Purchaser Notification

Corporate Headquarters

5791 Van Allen Way Carlsbad, CA 92008 USA Phone: +1 760 603 7200 Fax: +1 760 602 6500 Email: techsupport@lifetech.com

European Headquarters

Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF UK Phone: +44 141 814 6100 Toll-Free Phone: 0800 269 210 Toll-Free Tech: 0800 838 380 Fax: +44 141 814 6260 Tech Fax: +44 141 814 6117 Email: euroinfo@invitrogen.com Email Tech: eurotech@invitrogen.com

Japanese Headquarters

LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Japan Phone: +81 3 5730 6509 Fax: +81 3 5730 6519 Email: jpinfo@invitrogen.com

Additional international offices are listed at www.lifetechnologies.com

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

Obtaining Support

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of
 analysis, citations, and other product support documents
- · Obtain information about customer training
- Download software updates and patches

SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

Certificate of Analysis

The Certificate of Ánalysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to **www.lifetechnologies.com/support** and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact **outlicensing@lifetech.com** or Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners. Tween is a registered trademark of Uniqema Americas, LLC.

©2013 Life Technologies Corporation. All rights reserved.

