

## Flow Cytometer Calibration Beads Protocol-At-A-Glance

This protocol describes how to calibrate flow cytometers with 488 nm laser lines that excite the green fluorescent proteins AcGFP1 (*Aequorea coerulea* GFP) and EGFP, or flow cytometers with 561 nm laser lines that excite the red fluorescent protein, mCherry. Calibrate 488 nm laser lines using the **AcGFP Flow Cytometer Calibration Beads** (Cat. No. 632594). Calibrate 561 nm laser lines using the **mCherry Flow Cytometer Calibration Beads** (Cat. No. 632595).

### Protocol: Preparing Calibration Beads for Flow Cytometer Analysis

1. Add 1 ml of 1X Flow Cytometer Calibration Beads Dilution Buffer to a standard flow cytometer sample tube (BD Falcon 5 mL Polystyrene Round-Bottom Tubes, Disposable, VWR Cat. No. 60819-310).
2. Invert the stock tube of Calibration Beads 5 to 10 times to resuspend the beads.
3. After the beads are fully resuspended, transfer 20 µl of the bead suspension to the flow cytometer sample tube containing the dilution buffer.
4. Cap the sample tube and invert it 5 times to mix the diluted bead suspension.
5. Insert the sample tube into the sample port of the flow cytometer and run an analysis on the instrument.

Contact Us For Assistance	
Customer Service/Ordering	Technical Support
Telephone: 800.662.2566 (toll-free)	Telephone: 800.662.2566 (toll-free)
Fax: 800.424.1350 (toll-free)	Fax: 800.424.1350 (toll-free)
Web: <a href="http://www.clontech.com">www.clontech.com</a>	Web: <a href="http://www.clontech.com">www.clontech.com</a>
E-mail: <a href="mailto:orders@clontech.com">orders@clontech.com</a>	E-mail: <a href="mailto:tech@clontech.com">tech@clontech.com</a>

---

### Notice to Purchaser

Clontech® products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Clontech Laboratories, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Clontech and the Clontech logo are trademarks of Clontech Laboratories, Inc. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech is a Takara Bio Company. ©2013 Clontech Laboratories, Inc.

This document has been reviewed and approved by the Clontech Quality Assurance Department.