invitrogen USER GUIDE

# Rainbow Calibration Particles

Catalog Numbers A34305

Pub. No. MAN0017318 Rev. B.0

<u>^</u>!\

**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

### **Product description**

Rainbow Calibration Particles contain a mixture of 3.0 µm particles in 8 different fluorescent intensities. The calibration particles can be used for routine calibration of flow cytometers using wavelengths ranging from 365–650 nm.

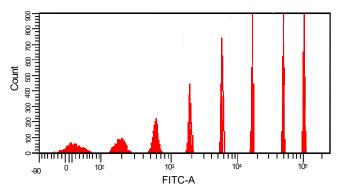


Fig. 1 Histogram analysis of particle fluorescence at 488-nm excitation and 530/30-nm emission.

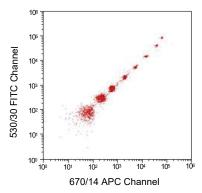


Fig. 2 Particle fluorescence plotted at 488-nm excitation and 530/30-nm emission versus 633-nm excitation and 670/14-nm emission.

## Contents and storage

Component	Amount	Composition	Storage
Rainbow Calibration Particles	5 mL	$1 \times 10^7$ particles/mL in deionized water with 0.02% sodium azide and 0.01% NP-40	2–8°C DO NOT FREEZE Protect from light

#### Product use

For routine flow cytometer calibration, dilute 3-5 drops of Rainbow Calibration Particles in 1 mL of buffer.

### 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 at <a href="https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html">www.thermofisher.com/us/en/home/global/terms-and-conditions.html</a>. If you have any questions, please contact Life Technologies at <a href="https://www.thermofisher.com/support">www.thermofisher.com/support</a>.



Life Technologies Corporation | 29851 Willow Creek | Eugene, Oregon 97402 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0017318

· · · · · · · · · · · · · · · · · · ·		
Revision	Date	Description
B.0	18 November 2021	Updating the Figure 1 graph.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

