

# Rainbow Calibration Particles

Catalog nos. A34305

Pub. No. MAN0017318

Rev. A.0

## Product information

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 nm to 650 nm.

Table 1. Contents and storage

Material	Amount	Composition	Storage
Rainbow Calibration Particles	5 mL	1 × 10 <sup>7</sup> particles/mL in deionized water with 0.02% sodium azide and 0.01% NP-40	<ul style="list-style-type: none"> <li>• 2–8°C</li> <li>• DO NOT FREEZE</li> <li>• Protect from light</li> </ul>

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

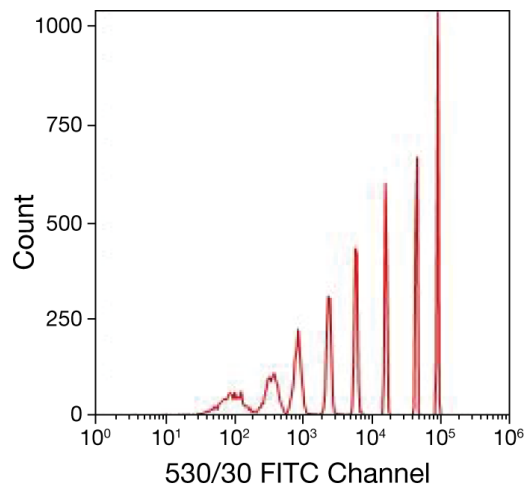
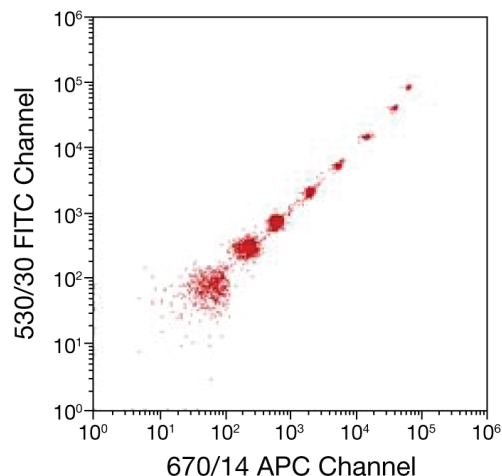


Figure 2. Particle fluorescence plotted at 488-nm excitation and 530/30-nm emission versus 633-nm excitation and 670/14-nm emission.



## Methods

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 found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

---



Manufacturer: Spherotech, Inc. | 27845 Irma Lee Circle | Lake Forest, IL 60045

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

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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
A.0	22 August 2017	New user guide.

**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.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.