

RF002



1. Add 2 drops per 10^6 fixed cells in 1 mL.



2. Incubate for 30 minutes at 25°C.



3. Proceed with flow cytometry.

For Research Use Only. Not for use in diagnostic procedures.

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

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

MAN0016216 Rev. A.0

ThermoFisher
SCIENTIFIC