



1. Add 1 drop per 10^5 cells in 0.5 mL of annexin-binding buffer or media with 2.5 mM calcium.



2. Incubate for 15 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.

Manufacturer: Life Technologies Corporation | 29851 Willow Creek Road | Eugene, OR 97402

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

MAN0017070 Rev. A