

1. Add 2 drops per mL of media.
2. Incubate for 30–60 minutes.
3. Image cells.

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

**Important Licensing Information:** This product 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

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

MAN0007449 Rev. A.0

**ThermoFisher**  
S C I E N T I F I C