



1. Add 2 drops per mL of media



2. Incubate 15–30 minutes



3. Image cells

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.

For Research Use Only. Not for use in diagnostic procedures. © 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. MAN0007236_Rev.1.0

