

# NucBlue™ Fixed Cell Stain (R37606)

## quick start guide

life



1. Add 2 drops per mL of media



2. Incubate 5 mins at 20-25°C



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 intended for human or animal therapeutic or diagnostic use. © 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners. R37603\_QuickReferenceCard

Life Technologies Corporation  
29851 Willow Creek Road  
Eugene, OR USA 97405-8300

**molecular  
probes®**  
by *life technologies*