

ActinGreen™ 488 ReadyProbes® Reagent

Cat. no. R37122

Pub. no. MAN0010749

Revision A.0

WARNING! For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

| Product | Amount | Storage |
|---|--------|--|
| ActinGreen™ 488 ReadyProbes® Reagent | 2.5 mL | <ul style="list-style-type: none">• Room temperature• Do not freeze |

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

Quick Protocol



1. Apply to fixed and permeabilized cells using 2 drops per mL of media.



2. Incubate 30 minutes.



3. Wash with PBS, HBSS, or LCIS.



4. Image cells.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate[s] warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE[S] DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE[S] BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit lifetechnologies.com/support
lifetechnologies.com