

Image-IT[®] Hypoxia Reagent

Pub. no. MAN0013415

Revision A.0

Detailed protocol is available online at www.lifetechnologies.com/manuals.

Product	Amount	Cat. no.	Storage*
Image-IT [®] Hypoxia Reagent	1 mg	H10498	<ul style="list-style-type: none">• $\leq -20^{\circ}\text{C}$• Protect from light• Dessicate

* When stored as directed, product is stable for at least 6 months.

MW: 711.87**Approximate excitation/emission maxima:** 490/610 nm

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

1. Plate cells at desired density on a glass bottom dish and incubate the cells overnight at 37°C in a CO₂ incubator.
2. Prepare 1 mM stock solution of Image-IT® Hypoxia Reagent in DMSO. Add the reagent to the cells at a final concentration of 1–10 µM and incubate for 15 to 30 minutes.
3. Place the cells in a hypoxia chamber attached to a fluorescence microscope and set to the desired oxygen concentration, and incubate for 1 hour.
4. Image the cells on the fluorescence microscope with excitation/emission of 490/610 nm.

Note: You can use the standard TRITC filter set for imaging; however, we recommend using FITC excitation and Texas Red® emission filters for best results.

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.

©2015 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