

HCS CellMask™ Stains

Cat. Nos. H32712, H32713, H32714, H32720, H32721

Rev. Date 1 October 2012

Pub. No. MAN0006875

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

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.

Chemical	CAS	Warning
DMSO	67-68-5	Components of the product may be absorbed into the body through the skin.

1. Prepare a 10-mg/mL HCS CellMask™ stock solution by dissolving the entire contents of Component A in 25 µL of Component B.
2. Prepare 4% paraformaldehyde and 0.1% Triton® X-100 solutions (available separately in the Image-iT® Fixation/Permeabilization Kit, Cat. no. R37602).
3. Prepare HCS CellMask™ staining solution by adding 2 µL of HCS CellMask™ stock solution (from step 1) to 10 mL of PBS.
4. Remove cell medium and add 4% paraformaldehyde solution. Incubate for 15 minutes at room temperature.

For research use only. Not for use in diagnostic procedures.

- Wash fixed cells 2–3 times with PBS and add 0.1% Triton® X-100 solution to each well. Incubate the cells for 15 minutes at room temperature.
- Wash the cells 2–3 times with PBS. Add 100 µL of HCS CellMask™ staining solution (from step 3) to each well and incubate for 30 minutes at room temperature.
- Wash each well with PBS 2–3 times. Image under appropriate filters.

Spectral Characteristics of the Fluorescent Dyes

HCS CellMask™ Stain	Ex/Em (maxima, in nm)
HCS CellMask™ Blue stain	346/442
HCS CellMask™ Green stain	493/516
HCS CellMask™ Orange stain	556/572
HCS CellMask™ Red stain	588/612
HCS CellMask™ Deep Red stain	650/655

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.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. Triton is a registered trademark of Dow Chemical Co.

For support visit www.lifetechnologies.com/support

www.lifetechnologies.com