

BTC, acetoxymethyl (AM) ester

Cat. no. B6791

Pub. no. MAN0006890

Revision 1.0

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

WARNING! Before handling, read all applicable Safety Data Sheets (SDSs) at www.lifetechnologies.com/support.

Quick Protocol

1. Add 100 μ L DMSO to 100 μ g of dye to make a \sim 1 mM stock.
2. Dilute the stock 1:1,000 in a buffered physiological medium of choice (e.g., Phosphate-Buffered Saline).
3. Incubate cells in the working solution (from step 2) for 15–60 minutes.
4. Wash the cells and image.
5. Calibration is required; refer to the product manual for details at www.lifetechnologies.com/manuals.

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

Guidelines for using BTC AM Ester

- Empirically determine the exact loading concentration, time, and temperature. Use the minimum dye concentration required to yield fluorescence signals with adequate signal-to-noise level.
- Add non-ionic detergent Pluronic® F-127 to assist in dispersing the non-polar AM ester in aqueous media.
- You may add the organic anion transport inhibitors probenecid (1–2.5 mM) or sulfinpyrazone (0.1–0.25 mM) to the cell medium to reduce leakage of the dye.
- Properties of the BTC Indicator:

Indicator	Ion-free Ex/Em*	Ion-bound Ex/Em*
BTC	464/533	401/529

*Ex/Em: Fluorescence excitation and emission maxima in nm.

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.

© 2013 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. Pluronic is a registered trademark of BASF Corporation.

For support visit lifetechnologies.com/support
lifetechnologies.com