

# PRODUCT INFORMATION



## BCECF AM

Item No. 15922

**Formal Name:** 3',6'-bis[(acetyloxy)methoxy]-5(or 6)-[[[(acetyloxy)methoxy]carbonyl]-3-oxo-spiro[isobenzofuran-1(3H),9'-[9H]xanthene]-2',7'-dipropanoic acid, 2',7'-bis[(acetyloxy)methyl] ester

**Synonym:** BCECF Acetoxymethyl ester

**MF:** C<sub>42</sub>H<sub>40</sub>O<sub>21</sub>

**FW:** 880.7

**Purity:** ≥95%

**Ex./Em. Max:** 440, 490/535 nm

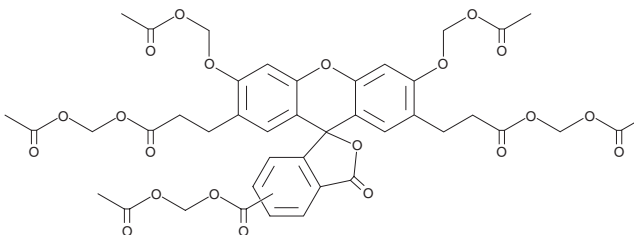
**Supplied as:** An oil

**Storage:** -20°C

**Stability:** ≥4 years

**Special Conditions:** Protect from light and moisture

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



### Laboratory Procedures

BCECF AM is supplied as an oil. A stock solution may be made by dissolving the BCECF AM in the solvent of choice. BCECF AM is soluble in organic solvents such as acetonitrile, DMSO, and dimethyl formamide.

### Description

BCECF AM is a widely-used fluorescent indicator for estimating intracellular pH.<sup>1,2</sup> This membrane-permeant compound is modified by intracellular esterases to produce BCECF, a polar fluorescein derivative that is retained by cells. Intracellular BCECF is stable and has an efflux half-life of greater than two hours.<sup>3</sup> Intracellular pH measurements are made by ratio imaging, which involves determining the pH-dependent ratio of emission intensity (detected at 535 nm) when the dye is excited at 490 nm vs. the emission intensity when excited at 440 nm.<sup>2,4</sup> This approach is amenable to either spectrofluorometry or flow cytometry.<sup>2,5</sup> BCECF AM can also be used to investigate intracellular changes in other ions, including potassium.<sup>6</sup>

### References

1. Rink, T.J., Tsien, R.Y., and Pozzan, T. Cytoplasmic pH and free Mg<sup>2+</sup> in lymphocytes. *J. Cell Biol.* **95**(1), 189-196 (1982).
2. Ozkan, P. and Mutharasan, R. A rapid method for measuring intracellular pH using BCECF-AM. *Biochim. Biophys. Acta* **1572**(1), 143-148 (2002).
3. Dive, C., Cox, H., Watson, J.V., et al. Polar fluorescein derivatives as improved substrate probes for flow cytometry assay of cellular esterases. *Mol. Cell Probe* **2**(2), 131-145 (1988).
4. O'Connor, N. and Silver, R.B. Ratio imaging: Practical considerations for measuring intracellular Ca<sup>2+</sup> and pH in living cells. *Methods Cell Biol.* **81**, 415-433 (2007).
5. Chow, S., Hedley, D., and Tannock, I. Flow cytometric calibration of intracellular pH measurements in viable cells using mixtures of weak acids and bases. *Cytometry* **24**(4), 360-367 (1996).
6. Balkay, L., Márián, T., Emri, M., et al. Flow cytometric determination of intracellular free potassium concentration. *Cytometry* **28**(1), 42-49 (1997).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/12/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM