

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: BCECF, AM ester (2',7'-bis-(2-carboxyethyl)-5-(and 6-)carboxyfluorescein,

acetoxymethyl ester)

CATALOG #: 51011/51011-1/51012

MOLECULARBCECF AM ester is a mixture of three components, the structures of which are shown below. MWt: N/A [117464-70-7]

PROPERTIES:

Color & FormWhite solidPurity $\geq 95\%$ by HPLC

Solubility Soluble in DMSO and DMF

Absorption/Emission 508 nm/530 nm (for BCECF acid)

Extinction N/A

Coefficient

STORAGE AND Store desiccated at -20 °C. Protect from light, especially when in solution.

HANDLING:

APPLICATION: Membrane-permeant form of BCECF that can be loaded into cells via incubation.

Once in the cells, BCECF/AM is readily hydrolyzed by cytosolic esterases to BCECF

acid. See BCECF acid (51010) for more information on BCECF.

Ref: 1) Ho, A.K., et al. Am. J. Physiol. 261, C642(1991); 2) Futsaetner, C.M., et al. Can. J. Microbiol. 39, 180(1993); 3) Shepherd, R.M., et al. Am. J. Physiol. 262, C182(1992); 4) Hayasi, H., et al. Am. J. Physiol. 262, C628(1992); 5) Weiner,

I.D. and Hamm, L.L. *Am. J. Physiol.* **256**, F957(1989).

TOXICITY: Unknown

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on

clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.