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 Acid

CATALOG #: 51010

MOLECULAR $C_{27}H_{20}O_{11}$ **INFORMATION:** MWt: 520.5

[85138-49-4]

HO₂CH₂CH₂CO₂H
CO₂H
CO₂H

PROPERTIES:

Color & FormYellow orange solidPurity $\geq 95\%$ by HPLCSoluble in water (pH>6)

Absorption/Emission 508 nm/530 nm

Extinction 90,000

Coefficient

STORAGE AND HANDLING:

Store desiccated at 4 °C upon receipt. Protect from light, especially when in solution.

APPLICATION:

BCECF (full chemical name: 2',7'-bis-(carboxyethyl)-5-(and-6)-carboxyfluorescein) is the most widely used fluorescent pH sensor. Because its pKa (6.97) is close to physiological pH, BCECF can detect cytosolic pH change with high sensitivity. At low pH, the dye is weakly fluorescent but becomes more fluorescent with increasing pH. The excitation spectrum of the dye undergoes a slight shift during pH change, while the wavelength of the emission maximum remains unchanged. The pH is determined ratiometrically by the relative fluorescent intensities at 535 nm when the dye is excited at 439 nm and 505 nm respectively. BCECF acid is membrane-impermeant and therefore has to be loaded into cells via microinjection or scrape loading (See **51011**, **BCECF/AM**).

Ref: 1) Kolber, M.A., et al. *J. Immunol. Meth.* **108**, 255(1988); 2) Carpenter, L., et al. *Biochem. J.* **304**, 751(1994); 3) Sellers, J.R. et al. *J. Immunol. Meth.* **172**, 255(1994).

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.