## BCECF AM (ultra pure)

pH indicator

BCECF AM (2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein acetoxymethyl ester) is a well established fluorescent probe that enables ratiometric monitoring of cellular pH. Intracellular pH plays an important role in modulating cellular events, including growth, calcium regulation, enzymatic activity, receptor-mediated signal transduction, ion transport, endocytosis, chemotaxis, adhesion and other cellular processes. Once inside, cellular esterases cleave the AM groups yielding a more hydrophilic BCECF trapped inside the cell. Wavelength Maxima: Excitation 505nm, Emission 520nm

Citations: 2

View Online »

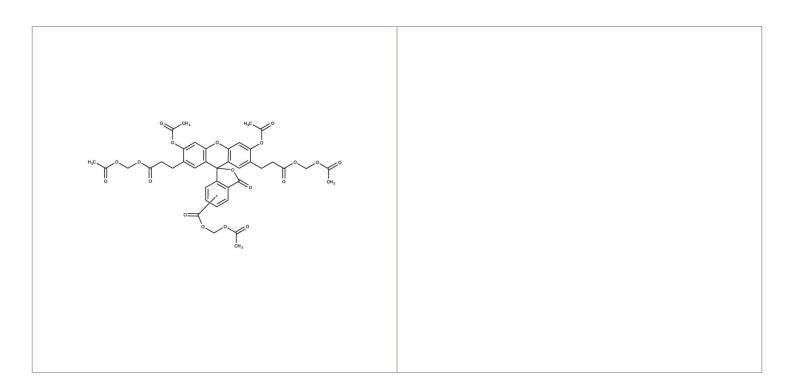
**Ordering Information** 

**Order Online** »

**ENZ-52102** 1mg

Manuals, SDS & CofA

**View Online** »



## **Handling & Storage**

**Use/Stability** Stable for at least one year after receipt when stored as recommended.

**Handling** Protect from light. Keep cool and dry.

Long Term Storage -20°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name 2',7'-bis-(2-Carboxyethyl)-5(6)-carboxyfluorescein

acetoxymethyl ester

**Appearance** Off white solid

CAS 117464-70-7

Formula  $C_{39}H_{36}O_{19}$ 

**MW** 808.7

Purity ≥95% (HPLC)

**Solubility** Soluble in DMSO.

**Technical Info / Product Notes**This product is a member of the CELLESTIAL<sup>®</sup> product

line, reagents and assay kits comprising fluorescent

molecular probes that have been extensively benchmarked for live cell analysis applications.

CELLESTIAL<sup>®</sup> reagents and kits are optimal for use in demanding imaging applications, such as confocal

microscopy, flow cytometry and HCS, where consistency

and reproducibility are required.

Last modified: May 29, 2024

