BCECF, AM Ester

BCECF, AM ester is membrane-permeant and thus can be loaded into cells via incubation. This product is a mixture of three molecular species, all of which are readily hydrolyzed by endogenous esterases into a single BCECF free acid form once they are in the cells.



Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

CAS number	117464-70-7
Cell permeability	Membrane permeant
рКа	pKa 6.97
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	508/530 nm (high pH) (after hydro

Product Description

BCECF, AM ester is membrane-permeant and thus can be loaded into cells via incubation. This product is a mixture of three molecular species, all of which are readily hydrolyzed by endogenous esterases into a single BCECF free acid form once they are in the cells. For more information on BCECF, see 51010.

- White solid soluble in DMSO and DMF
- Store at -20°C and protect from light, especially in solution
- MW: approx. 688
- [117464-70-7]

$$\begin{array}{c|c} O & O & O & O \\ \hline O & O & O & O \\ \hline O & O & O & O \\ O & O & O & O \\ \hline O & O$$

References

1) DOI: 10.1152/ajpcell.1991.261.4.C642

2) DOI: 10.1139/m93-025

3) DOI: 10.1152/ajpcell.1992.262.1.C182 4) DOI: 10.1152/ajpcell.1992.262.3.C628 5) DOI: 10.1152/ajprenal.1989.256.5.F957

This datasheet was generated on November 5, 2025 at 02:55:29 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bcecf-am-ester/