

BCECF, AM Ester, 1 mg/mL in Anhydrous DMSO

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

CAS number	117464-70-7
Cell permeability	Membrane permeant
pKa	pKa 6.97
Indicator type	Ratiometric
Colors	Green

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](#).

- 1 mg/mL in anhydrous DMSO
- Store at -20°C and protect from light and moisture
- MW: approx. 615
- [117464-70-7]

