

## 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.



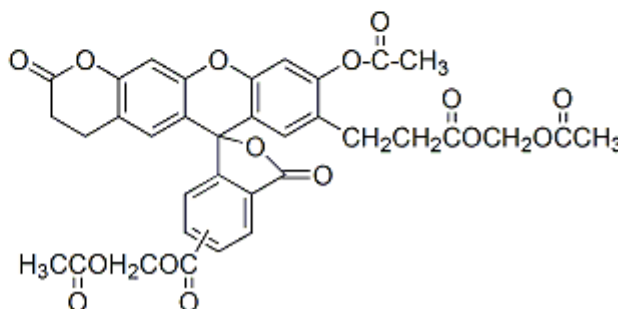
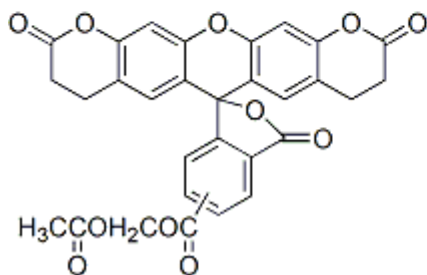
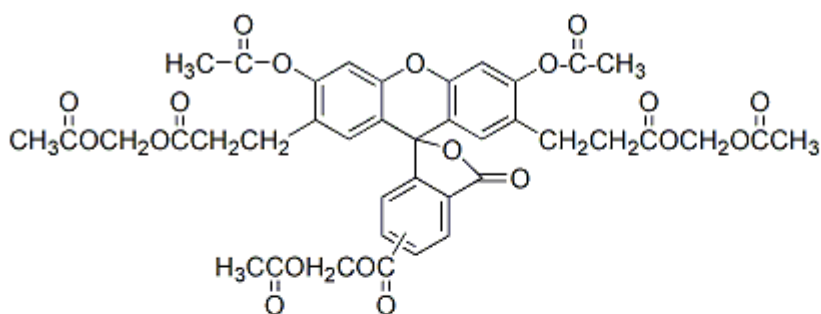
### Product attributes

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

## 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]



## 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