

DiOC₁₄(3), Hydroxyethanesulfonate

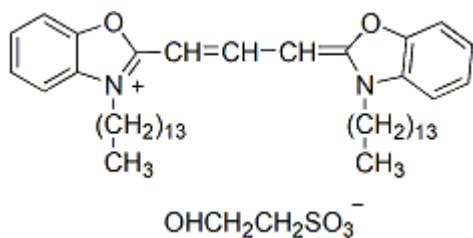
DiOC₁₄(3) is a DiO derivative, but is more soluble in aqueous media than the latter because it has shorter hydrocarbon chains. Also see our ready-to-use CellBrite™ Cell Membrane Labeling Solutions.



Product Description

DiOC₁₄(3) is a DiO derivative, but is more soluble in aqueous media than the latter because it has shorter hydrocarbon chains. Also see our ready-to-use CellBrite™ Cell Membrane Labeling Solutions.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 484/501 nm
- ϵ (MeOH) = 150,000
- Orange solid soluble in methanol, ethanol or DMSO
- Store desiccated at 4 °C
- C₄₇H₇₄N₂O₆S
- MW: 795



Product attributes

Probe cellular localization	Membrane/cell surface, Membrane/vesicular
For live or fixed cells	For fixed cells, For live/intact cells
Assay type/options	Co-cultures, Extended staining (several days to weeks)
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
Colors	Green
Excitation/Emission	484/501 nm