

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 CellBriteTM Cell Membrane Labeling Solutions.



Product attributes

Call us: 800-304-5357

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

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

This datasheet was generated on November 23, 2025 at 03:08:31 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dioc143-hydroxyethanesulfonate/