

Neuro-Dil in Vegetable Oil

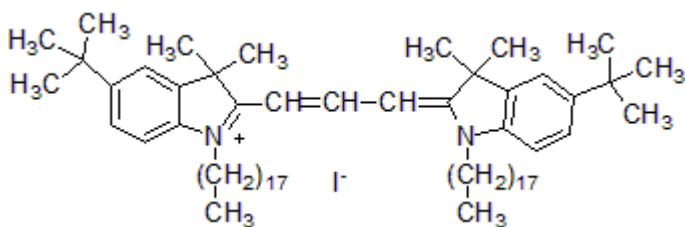
Neuro-Dil uniformly dissolved in vegetable oil in a formulation optimized for microinjection. Neuro-Dil was developed at Biotium as an alternative to the widely used fluorescent membrane probe Dil.



Product Description

Neuro-Dil uniformly dissolved in vegetable oil in a formulation optimized for microinjection. Neuro-Dil was developed at Biotium as an alternative to the widely used fluorescent membrane probe Dil. Like [Dilinoyley Dil](#), Neuro-Dil has structural features that may make the probe diffuse faster than Dil on cell membranes. However, Neuro-Dil dye has saturated carbon chains, making it more hydrophobic than Dilinoyley Dil, for potentially more stable labeling with less dye transfer between cells. Neuro-Dil has nearly identical absorption and emission wavelengths to those of Dil. We also offer [Neuro-Dil](#) in solid form.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 549/565 nm
- Store at room temperature and protect from light; do not freeze
- $C_{67}H_{115}IN_2$
- MW: 1072



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	Red
Excitation/Emission	549/565 nm

This datasheet was generated on January 9, 2026 at 11:55:54 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/neuro-dil-in-vegetable-oil/>