

Neuro-Dil in Vegetable Oil

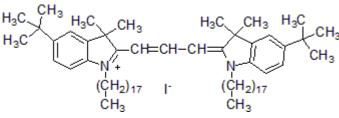
Neuro-Dil uniformly dissolved in vegetable oil in a formulation optimized for microinjection. Neuro-Dil was developed at Biotium as an alternative to the widely used fluorescent membrane probe Dil



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 Dilinoleyl 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 Dilinoleyl 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
- C₆₇H₁₁₅IN₂
- MW: 1072



Product link: https://biotium.com/product/neuro-dii-in-vegetable-oil/

This datasheet was generated on November 24, 2025 at 11:13:17 PM. Visit product page to check for updated information before use.

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	Red
Excitation/Emission	549/565 nm