

## 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.



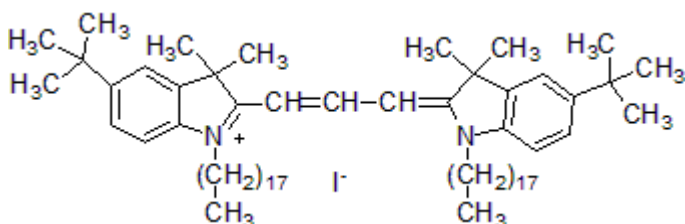
### Product attributes

<b>Probe cellular localization</b>	Membrane/cell surface, Membrane/vesicular
<b>For live or fixed cells</b>	For fixed cells, For live/intact cells
<b>Assay type/options</b>	Co-cultures, Extended staining (several days to weeks)
<b>Fixation options</b>	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
<b>Colors</b>	Red
<b>Excitation/Emission</b>	549/565 nm

## 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



This datasheet was generated on June 9, 2026 at 08:52:04 PM. Visit product page to check for updated information before use.

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