

## Neuro-Dil

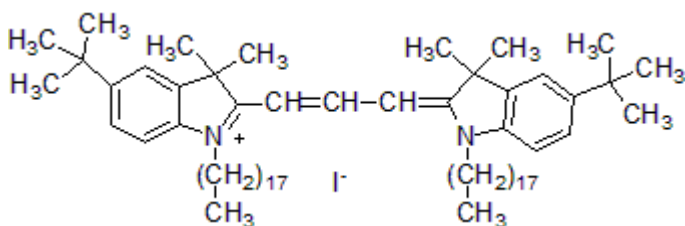
Neuro-Dil was developed at Biotium as an alternative to the widely used fluorescent membrane probe Dil. Neuro-Dil has structural features that may make the probe diffuse faster than Dil on cell membranes and also may result in a more stable labeling.



## Product Description

Neuro-Dil was developed at Biotium as an alternative to the widely used fluorescent membrane probe Dil. Like [Dilinoyleyl 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 Dilinoyleyl 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 Vegetable Oil](#) for microinjection studies. Also see our [CellBrite™ Orange Cytoplasmic Membrane Dye](#), a ready-to-use dye solution for cell labeling.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 549/565 nm
- $\epsilon$  = 148,000
- Red solid soluble at 1-2 mM (with heating) in DMF, DMSO, or ethanol
- Soluble at 1-2 mM in vegetable oil with heating and sonication
- Store at 4 °C and protect from light
- $C_{67}H_{115}IN_2$
- MW: 1072



## References

1. Proc Natl Acad Sci (2006) Feb 21;103(8):2938-42 [doi: 10.1073/pnas.0511159103](#)
2. J Neurosci. (2013) Oct 23;33(43):16853-64 [doi: 10.1523/JNEUROSCI.1844-13.2013](#)
3. PLoS One (2013) May 31;8(5):e64907 [doi: 10.1371/journal.pone.0064907](#)
4. J Neurosci (2013) Apr 10;33(15):6454-9 [doi: 10.1523/JNEUROSCI.0178-13.2013](#)
5. Bioconjug Chem (2014) Dec 17;25(12):2134-43 [doi: 10.1021/bc500465j](#)
6. Front Syst Neurosci (2016) Nov 15;10:95 [doi: 10.3389/fnsys.2016.00095](#)

This datasheet was generated on January 10, 2026 at 02:52:53 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/neuro-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  |