

Glowing products for science

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 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 Vegetable Oil for microinjection studies. Also see our CellBriteTM Orange Cytoplasmic Membrane Dye, a ready-to-use dye solution for cell labeling.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 549/565 nm
- $\varepsilon = 148,000$
- 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

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

- C₆₇H₁₁₅IN₂
- 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 November 25, 2025 at 12:49:45 AM. 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