

Neuro-DiO in Vegetable Oil

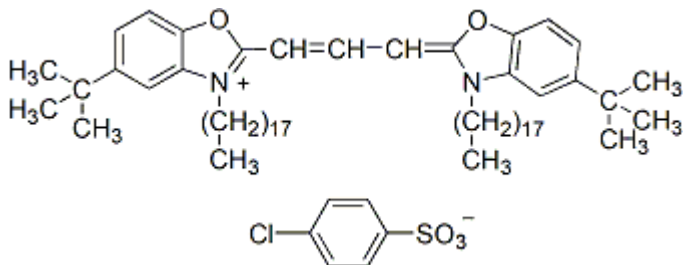
Neuro-DiO is a derivative of DiO and has spectral properties nearly identical to that of the latter. However, unlike DiO, Neuro-DiO is structurally more similar to DiI, which is known to be a much better lipophilic neuronal tracer than regular DiO.



Product Description

Neuro-DiO is a derivative of DiO and has spectral properties nearly identical to that of the latter. However, unlike DiO, Neuro-DiO is structurally more similar to DiI, which is known to be a much better lipophilic neuronal tracer than regular DiO. Like DiI in oil, Neuro-DiO is uniformly dissolved in vegetable oil in a formulation optimized for microinjection. The green fluorescent Neuro-DiO and red fluorescent DiI in oil can be used in combination for two color imaging. We also offer [Neuro-DiO](#) in solid form, and in our [CellBrite™ Green Cytoplasmic Membrane Dye](#), a ready-to-use dye solution for cell labeling.

- $\lambda_{Ex}/\lambda_{Em}$ = 484/501 nm
- Yellow-orange oil
- Store at room temperature and protect from light (do not freeze)
- $C_{67}H_{105}ClN_2O_5S$
- MW: 1086



This datasheet was generated on December 24, 2025 at 07:42:02 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/neuro-dio-in-vegetable-oil/>

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	Green
Excitation/Emission	484/501 nm