

## 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.



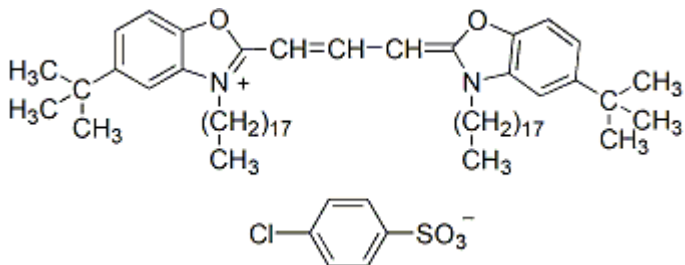
### Product attributes

<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>	Green
<b>Excitation/Emission</b>	484/501 nm

## 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 June 15, 2026 at 02:02:13 AM. Visit product page to check for updated information before use.

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