

Dilinoleyl DiO

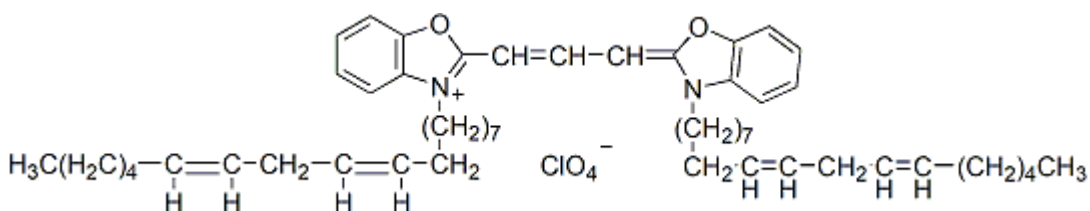
Dilinoleyl DiO (1,1'-dilinoyleyl-3,3'-oxacarbocyanine perchlorate), also known as FAST DiO™, has identical absorption and emission spectra as DiO (#[60011](#)).



Product Description

Dilinoleyl DiO (1,1'-dilinoyleyl-3,3'-oxacarbocyanine perchlorate), also known as FAST DiO™, has identical absorption and emission spectra as DiO (#[60011](#)). However, due to its unsaturated hydrocarbon chains, dilinoleyl DiO has a much faster lateral diffusion rate in cell membranes than DiO.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 484/499 nm
- ϵ = 140,000
- Yellow orange solid soluble in ethanol, DMF or DMSO
- Store at -20 °C
- $C_{53}H_{77}ClN_2O_6$
- MW: 873.65



Fast DiO is a trademark of Thermo Fisher Scientific.

References

1. Neuron 17, 91 (1996).

This datasheet was generated on June 14, 2026 at 07:08:48 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dilinoleyl-dio-also-called-fast-dio1-or-11-dilinoleyl-33-oxacarbocyanine-perchlorate/>

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	484/499 nm