Dilinoleyl DiO

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



Product Description

Dilinoleyl DiO (1,1'-dilinoleyl-3,3'-oxacarbocynanine perchlorate), also known as FAST DiOTM, 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.

- λεx/λεm (MeOH) = 484/499 nm
- ε = 140,000
- Yellow orange solid soluble in ethanol, DMF or DMSO
- Store at -20°C
- C₅₃H₇₇CIN₂O₆
- MW: 873.65

Fast DiO is a trademark of Thermo Fisher Scientific.

References

1. Neuron 17, 91 (1996).

This datasheet was generated on November 4, 2025 at 10:32:20 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-oxacarbocynanine-perchlorate/

Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

| 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 |
| Fixation options | Fix before staining (formaldehyde), Fix after sta (formaldehyde), Permeabilize before staining |
| Colors | Green |
| Excitation/Emission | 484/499 nm |