### Glowing products for science

# **Dilinoleyl Dil**

Dilinoleyl Dil (1,1'-dilinoleyl-3,3,3',3'-tetramethylindocarbocynanine perchlorate), also known as FAST Dil $^{TM}$ , has identical absorption and emission spectra as Dil (#60010) .



## **Product Description**

Dilinoleyl Dil (1,1'-dilinoleyl-3,3,3',3'-tetramethylindocarbocynanine perchlorate), also known as FAST DilTM, has identical absorption and emission spectra as DilTM (#60010). Due to its unsaturated hydrocarbon chains, dilinoleyl Dil has a much faster lateral diffusion rate in cell membranes than Dil. This property makes dilinoleyl Dil particularly useful for tracing neurons in tissues. Also see Neuro-Dil.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 549/565 nm
- $\varepsilon = 148,000$
- Dark red oily paste soluble in ethanol, DMF or DMSO
- Store at -20°C and protect from light
- C<sub>59</sub>H<sub>89</sub>CIN<sub>2</sub>O<sub>4</sub>
- MW: 925.82

Chemical Structure:

Fast Dil is a trademark of Thermo Fisher Scientific.

### References

- 1. Neuroreport 8, 3965 (1997).
- 2. Biophys J 70, 988 (1996).
- 3. J Cell Biol 134, 1427 (1996).

This datasheet was generated on November 4, 2025 at 02:21:20 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dilinoleyl-dii-also-called-fast-dii1-or-11-dilinoleyl-3333-tetramethylindocarbocynanine-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	Red
Excitation/Emission	549/565 nm

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