

Dilinoleyl Dil

Dilinoleyl Dil (1,1'-dilinoyleyl-3,3,3',3'-tetramethylindocarbocyanine perchlorate), also known as FAST Dil™, has identical absorption and emission spectra as Dil ([#60010](#)) .



Product attributes

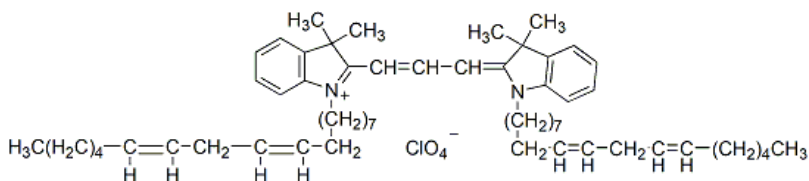
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
For live or fixed cells	For fixed cells, For live/intact cells
Assay type/options	Co-cultures, Extended staining (several days to weeks)
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Permeabilize before staining
Colors	Red
Excitation/Emission	549/565 nm

Product Description

Dilinoleyl Dil (1,1'-dilinoyleyl-3,3,3',3'-tetramethylindocarbocyanine perchlorate), also known as FAST Dil™, has identical absorption and emission spectra as Dil™ ([#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
- $\epsilon = 148,000$
- Dark red oily paste soluble in ethanol, DMF or DMSO
- Store at -20°C and protect from light
- $C_{59}H_{89}ClN_2O_4$
- 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 June 19, 2026 at 02:08:29 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/dilinoleyl-dii-also-called-fast-dii1-or-1-1-dilinoleyl-3333-tetramethylindocarbocyanine-perchlorate/>