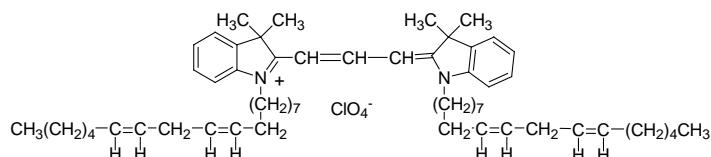


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** **Dilinoleyl DiI** (also called FAST DiI or 1,1'-dilinoleyl-3,3,3',3'-tetramethylindocarbocyanine perchlorate)**CATALOG #:** **60034****MOLECULAR INFORMATION:** C<sub>59</sub>H<sub>89</sub>ClN<sub>2</sub>O<sub>4</sub>  
Mwt: 925.82**PROPERTIES:**

**Color & Form** Dark red oily solid  
**Purity** ≥ 98% by TLC  
**Solubility** Soluble in DMSO, DMF or EtOH  
 $\lambda_{ex} \backslash \lambda_{em}$  549 nm/565 nm (MeOH)  
**Extinction Coefficient** 145,000 (MeOH)

**STORAGE AND HANDLING:**

Store at 4°C and protect from light especially when in solution.

**APPLICATION:**

Dilinoleyl DiI and DiI (#60010) have identical absorption and emission spectra. However, the unsaturated hydrocarbon chains of dilinoleyl DiI make the dye have a much faster lateral diffusion rate in cell membranes than regular DiI. This property makes dilinoleyl DiI particularly useful for tracing neurons in tissues.

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

**TOXICITY:**

Unknown.

**FIRST AID:**

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.