

## CERTIFICATE OF ANALYSIS

**Catalog Number** C018  
**Product Name** 'DiD'; DiI<sub>C<sub>18</sub>(5)</sub>  
(1,1'-dioctadecyl-3,3,3',3'-tetramethylindodicarbocyanine, perchlorate)  
**Appearance** blue solid  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup> Maximum Extinction	644 nm 240,000 cm <sup>-1</sup> M <sup>-1</sup>	644 ± 4 nm 245,000 ± 20,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup> Emission Maximum	664 nm	663 ± 4 nm
<b>HPLC</b> Purity	96 % at 640 nm	≥ 90 % at 640 nm
<b>NMR</b> Result	meets specification	consistent with structure

1. Solvent: Methanol.