

## CERTIFICATE OF ANALYSIS

**Catalog Number** C019  
**Product Name** 'DiR'; DiIC<sub>18</sub>(7) (1,1'-dioctadecyl-3,3,3',3'-tetramethylindotricarbocyanine, iodide)  
**Appearance** blue/green solid  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup> Maximum Extinction	748 nm 263,000 cm <sup>-1</sup> M <sup>-1</sup>	748 ± 4 nm 270,000 ± 20,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup> Emission Maximum	775 nm	776 ± 4 nm
<b>HPLC</b> Purity	96 % at 750 nm	≥ 95 % at 750 nm
<b>NMR</b> Result	meets specification	consistent with structure

1. Solvent: Methanol.