

CERTIFICATE OF ANALYSIS

Catalog Number C017
Product Name 'DiI'; DiIC₁₈(3) (1,1'-dioctadecyl-3,3,3',3'-tetramethylindocarbocyanine perchlorate)
CAS Number 41085-99-8
Molecular Formula C₅₉H₉₇ClN₂O₄
Molecular Weight 933.88
Appearance red/purple solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹		
Maximum	549 nm	549 ± 4 nm
Extinction	149,000 cm ⁻¹ M ⁻¹	142,000 ± 12,000 cm ⁻¹ M ⁻¹
Fluorescence ¹		
Emission Maximum	565 nm	564 ± 4 nm
HPLC		
Purity	96 % at 550 nm	≥ 95 % at 550 nm
NMR		
Result	meets specification	consistent with structure

1. Solvent: Methanol.