

CERTIFICATE OF ANALYSIS

Catalog Number C016
Product Name 'DiO'; DiOC₁₈(3) (3,3'-dioctadecyloxacarbocyanine perchlorate)
CAS Number 34215-57-1
Molecular Formula C₅₃H₈₅ClN₂O₆
Molecular Weight 881.72
Appearance orange solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹		
Maximum	485 nm	485 ± 4 nm
Extinction	154,000 cm ⁻¹ M ⁻¹	160,000 ± 12,000 cm ⁻¹ M ⁻¹
Fluorescence ¹		
Emission Maximum	500 nm	502 ± 4 nm
HPLC		
Purity	96 % at 490 nm	≥ 95 % at 490 nm
NMR		
Result	meets specification	consistent with structure

1. Solvent: Methanol.