

CERTIFICATE OF ANALYSIS

Catalog Number C050
Product Name DiOC₆(3), (3,3'-dihexyloxacarbocyanine iodide)
CAS Number 53213-82-4
Molecular Formula C₂₉H₃₇IN₂O₂
Molecular Weight 572.53
Appearance red solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	484 nm 162,000 cm ⁻¹ M ⁻¹	484 ± 4 nm 155,000 ± 11,000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	500 nm	500 ± 4 nm
HPLC Purity	96 % at 484 nm	≥ 95 % at 484 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.