

CERTIFICATE OF ANALYSIS

Catalog Number C045
Product Name JC-1
(5,5',6,6'-tetrachloro-1,1',3,3'-tetraethylbenzimidazolylcarbocyanine iodide)
CAS Number 47729-63-5
Molecular Formula C₂₅H₂₇Cl₄I N₂
Molecular Weight 652.23
Appearance purple/red solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	514 nm 186,000 cm ⁻¹ M ⁻¹	514 ± 3 nm 197,000 ± 15,000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	533 nm	530 ± 4 nm
HPLC Purity	96 % at 514 nm	≥ 90 % at 514 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.