

CERTIFICATE OF ANALYSIS

Catalog Number C043
Product Name MitoView Deep Red, 1 mg/mL in DMSO
Molecular Formula $C_{34}H_{36}Cl_2N_2$
Molecular Weight 543.58
Appearance purple solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	640 nm 186,000 cm ⁻¹ M ⁻¹	640 ± 3 nm 194,000 ± 22,000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	663 nm	662 ± 4 nm
HPLC Purity	95 % at 640 nm	≥ 90 % at 640 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.