

CERTIFICATE OF ANALYSIS

Catalog Number C041
Product Name MitoBeacon™ Orange, 1 mg/mL in DMSO
Molecular Formula C₂₄H₂₄Cl₂N₂O
Molecular Weight 427.37
Appearance purple solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	552 nm 96,700 cm ⁻¹ M ⁻¹	551 ± 3 nm 102,000 ± 7,000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	576 nm	576 ± 4 nm
HPLC Purity	92 % at 254 nm	≥ 90 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.