

CERTIFICATE OF ANALYSIS

Catalog Number C011
Product Name ethidium homodimer-2 (EthD-2), 1 mM solution in DMSO
Appearance purple solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption¹ Optical Density	0.23 at 497 nm	0.21 ± 0.02 at 500 ± 4 nm

1. Solvent: H₂O. Method: Optical Density, 51× dilution