

CERTIFICATE OF ANALYSIS

Catalog Number C010
Product Name ethidium homodimer-1 (EthD-1)
CAS Number 61926-22-5
Molecular Formula C₄₆H₅₀Cl₄N₈
Molecular Weight 856.77
Appearance red solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹		
Maximum	496 nm	494 ± 4 nm
Extinction	9300 cm ⁻¹ M ⁻¹	9000 ± 700 cm ⁻¹ M ⁻¹
Fluorescence ¹		
Emission Maximum	630 nm	630 ± 4 nm
HPLC		
Purity	95 % at 254 nm	≥ 90 % at 254 nm
NMR		
Result	meets specification	consistent with structure

1. Solvent: H₂O.