

CERTIFICATE OF ANALYSIS

Catalog Number C008
Product Name propidium iodide, 1 mg/mL solution in water
CAS Number 25535-16-4
Molecular Formula C₂₇H₃₄I₂N₄
Molecular Weight 668.4
Appearance purple solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	494 nm 6100 cm ⁻¹ M ⁻¹	493 ± 3 nm 6000 ± 300 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	632 nm	635 ± 4 nm
HPLC Purity	95 % at 254 nm	≥ 92 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Buffer, 50 mM potassium phosphate, pH 7.