

CERTIFICATE OF ANALYSIS

Catalog Number C009
Product Name ethidium monoazide bromide (EMA)
CAS Number 58880-05-0
Molecular Formula C₂₁H₁₈BrN₅
Molecular Weight 420.31
Appearance red solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	464 nm 5900 cm ⁻¹ M ⁻¹	463 ± 4 nm 5400 ± 500 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	621 nm	625 ± 4 nm
HPLC Purity	95 % at 254 nm	≥ 95 % at 254 nm
NMR Result	meets specification	consistent with structure

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