

CERTIFICATE OF ANALYSIS

Catalog Number C002
Product Name DAPI, 1 mg/mL solution in water
CAS Number 28718-90-3
Molecular Formula $C_{16}H_{17}Cl_2N_5$
Molecular Weight 350.25
Appearance yellow solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	347 nm 32000 cm ⁻¹ M ⁻¹	348 ± 3 nm 31000 ± 3000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	455 nm	457 ± 4 nm
HPLC Purity	96 % at 254 nm	• 95 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.