

CERTIFICATE OF ANALYSIS

Catalog Number C001
Product Name 4',6-diamidino-2-phenylindole, dihydrochloride (DAPI)
CAS Number 28718-90-3
Molecular Formula C₁₆H₁₇Cl₂N₅
Molecular Weight 350.25
Appearance yellow solid
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.