

CERTIFICATE OF ANALYSIS

Catalog Number C006
Product Name Hoechst 33342, 10 mg/mL in water
CAS Number 23491-52-3
Molecular Formula C₂₇H₃₇Cl₃N₆O₄
Molecular Weight 615.99
Appearance yellow solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	354 nm 43000 cm ⁻¹ M ⁻¹	352 ± 3 nm 43000 ± 4000 cm ⁻¹ M ⁻¹
Fluorescence ¹ Emission Maximum	463 nm	461 ± 4 nm
HPLC Purity	95 % at 254 nm	≥ 95 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: 0.1 M HCl/Methanol.