

Catalog Number C004
Product Name Hoechst 33258, pentahydrate (bis-benzimide), 10 mg/mL in water
CAS Number 23491-45-4
Molecular Formula C₂₅H₃₇Cl₃N₆O₆
Molecular Weight 623.96
Appearance yellow solution
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹		
Maximum	353 nm	352 ± 3 nm
Extinction	44000 cm ⁻¹ M ⁻¹	47000 ± 4000 cm ⁻¹ M ⁻¹
Fluorescence ¹		
Emission Maximum	460 nm	461 ± 4 nm
HPLC		
Purity	96 % at 254 nm	≥ 95 % at 254 nm
NMR		
Result	meets specification	consistent with structure

1. Solvent: 0.1 M HCl/Methanol.