

## CERTIFICATE OF ANALYSIS

**Catalog Number** C003  
**Product Name** Hoechst 33258, pentahydrate (bis-benzimide)  
**CAS Number** 23491-45-4  
**Molecular Formula**  $C_{25}H_{37}Cl_3N_6O_6$   
**Molecular Weight** 623.96  
**Appearance** yellow solid  
**Lot Number**

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

1. Solvent: 0.1 M HCl/Methanol.