

## CERTIFICATE OF ANALYSIS

**Catalog Number** C005  
**Product Name** Hoechst 33342, trihydrochloride, trihydrate  
**CAS Number** 23491-52-3  
**Molecular Formula**  $C_{27}H_{37}Cl_3N_6O_4$   
**Molecular Weight** 615.99  
**Appearance** yellow solid  
**Lot Number**

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

1. Solvent: 0.1 M HCl/Methanol.