



Product Data Sheet

Product Name: Hoechst 33342

Lot Number: See label on vial

Catalog Number: AS-83218

Size: 5 mL

Molecular Weight: 561.9

Formula: $C_{27}H_{37}Cl_3N_6O_4$

Appearance: Clear liquid

Purity:

HPLC: >95%

TLC: >95%

For Research Use Only