

6-Aminohexanol

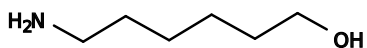
Catalog Number: 4002

Unit Size: 100 g

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid off white
Molecular Weight	117.19
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

6-Aminohexanol is a precursor that is widely used for preparing biotin and dye phosphoramidites.