

**6-Aminohexanol \*CAS 4048-33-3\***

Catalog Number: 4001

Unit Size: 25 g

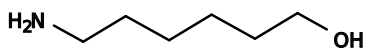
**Product Details**

---

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.