

**Biotin \*CAS 58-85-5\***

Catalog Number: 3001

Unit Size: 1 g

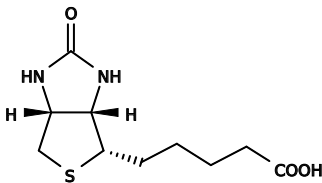
**Product Details**

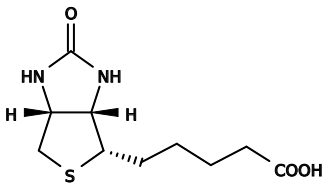
---

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

**Chemical Properties**

---

Appearance	Solid white
Molecular Weight	244.31
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

**Applications**

---

Biotin binds with high affinity to both avidins and streptavidins. Biotin and its derivatives can be conjugated to many biomolecules without significantly altering the biological activity of the target molecules since biotin is relatively a small molecule. A biopolymer (such as proteins) can react with several molecules of biotin that, in turn, can each bind one avidin. This characteristic greatly increases the sensitivity of many biological assays. Biotin derivatives are widely used for biological detections and purification. Our biotin is highly purified and tested for avidin-binding affinity.