

## Tris NTA amine

Catalog Number: 12637

Unit Size: 1 mg

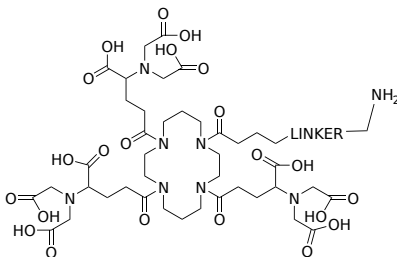
### Product Details

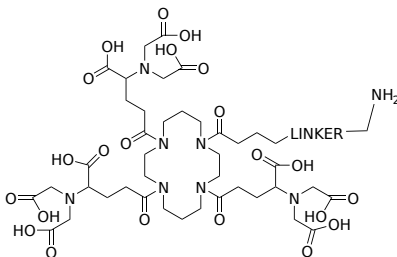
---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	1049.05
Soluble In	DMSO
Chemical Structure	



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

Tris NTA amine is an excellent building block to develop probes for detecting polyHIS tag proteins. Tris NTA compounds probably have the highest affinity for polyhistidine-labeled proteins. They can be used for the purification and analysis of polyhistidine-tagged proteins. AAT Bioquest offers a large family of fluorescent tris-NTA compounds that provide an efficient method for site-specific and stable noncovalent fluorescence labeling of polyhistidine-tagged proteins. In contrast to the transient binding of conventional mono-NTA, the multivalent interaction of tris-NTA compounds form a much more stable complex with polyhistidine-tagged proteins. The high selectivity of tris-NTA compounds toward cumulated histidines enable the selective labeling of proteins in cell lysates and on the surface of live cells.