

## IDA maleimide

Catalog Number: 12631

Unit Size: 5 mg

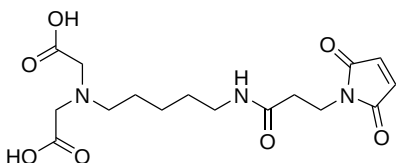
### Product Details

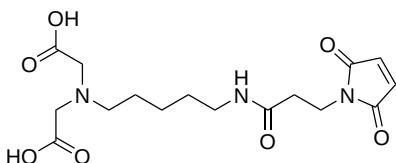
---

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

### Chemical Properties

---

Appearance	Solid white
Molecular Weight	369.37
Soluble In	DMSO
Chemical Structure	



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

Immobilized metal affinity chromatography (IMAC) is a popular method for protein purification, particularly for recombinant proteins fused to a polyhistidine-tag. Transition metal ions immobilized to a matrix through a chelating ligand interact with the polyhistidine-tag, effectively sequestering the fused protein from a sample. Nitrilotriacetic acid (NTA) and iminodiacetic acid (IDA) are two such ligands commonly used in commercially available resins. AAT Bioquest offer a variety of NTA building blocks for developing either NTA-based purification and detection. IDA is complimentary to NTA. The tridentate IDA ligand requires a lower imidazole concentration to elute protein than the tetradentate NTA. IDA is a smaller molecule which can be coupled to the matrix at a higher density resulting in a higher metal loading capacity. IDA maleimide is an excellent building block that can be used to develop a variety of IDA probes for detecting polyHis-containing proteins, and tools for purifying polyHis tagged proteins.