

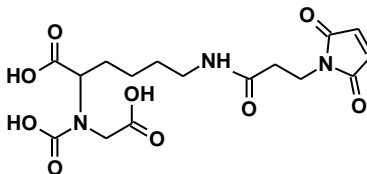
Product Information Sheet

Ordering Information

Product Number: 12609
Product Name: NTA maleimide
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 24 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 413.38
Appearance: Off-white powder



Chemical Structure:
Soluble in: DMSO
Excitation Wavelength: N/A
Emission Wavelength: N/A

Application Notes

NTA maleimide is used to attach NTA groups to a biological substrate or a surface that has thiol groups. Through the NTA moiety attached to the substrate or surface, genetically expressed proteins bearing the hexahistidine extension (HIS Tag) can be detected or immobilized via Ni-mediated His-Tag method. The NTA maleimide is an excellent building block for developing the HIS tag detection probes.