

HIS Lite™ iFluor™ 647 Tris NTA Chelator

Catalog Number: 12658

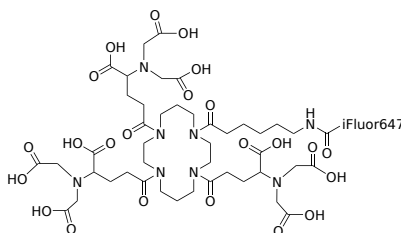
Unit Size: 100 ug

Product Details

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

Chemical Properties

Appearance	Solid blue
Molecular Weight	2006.23
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	656 nm
Emission Wavelength	670 nm

Applications

iFluor® 647 Tris-NTA Chelator is used as a sensitive red fluorescent probe for detecting polyhistidine-labeled proteins in combination with the addition of a heavy metal ion such as Ni²⁺ in cells, solution and solid surfaces. Fluorescent tris-NTA compounds 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 conjugated fluorophores form a much more stable complex with polyhistidine-tagged proteins. The high selectivity of tris-NTA compounds toward cumulated histidines enables the selective labeling of proteins in cell lysates and on the surface of live cells. Fluorescent tris-NTA conjugates can be applied for the analysis of a ternary protein complex in solution and on surfaces. In combination with OG488-tris-NTA compound iFluor® 647 Tris-NTA can be used for multicolor analysis of polyhistidine-tagged proteins. AAT Bioquest also offers HIS Lite™ iFluor™ 647 Tris NTA-Ni Complex that can be directly used for the specific and highly sensitive detection of His-tagged fusion proteins without the addition of a heavy metal ion. The transition metal ions (e.g., Ni ion)-mediated complexation of polyhistidine-labeled proteins with fluorescent tris-NTA conjugates provides a sensitive reporter for detecting and monitoring protein-protein interactions in real time.