

# HIS Lite™ Cy5 Tris NTA Chelator

Catalog Number: 12659

Unit Size: 100 ug

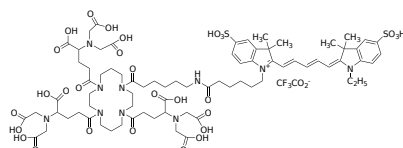
## Product Details

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

## Chemical Properties

Appearance	Solid blue
Molecular Weight	1801.87
Soluble In	Water

### Chemical Structure



## Spectral Properties

Excitation Wavelength	651 nm
Emission Wavelength	670 nm

## Applications

Cy5-Tris NTA compound is a sensitive fluorescent chelator for detecting polyhistidine-labeled proteins in cells, solution and solid surfaces with the addition of certain metal ions (such as Ni<sup>2+</sup>, Co<sup>2+</sup> etc.) as a mediator. It is the chelator of our HIS Lite™ Cy5 Tris NTA-Ni Complex (#12619). In combination with other color tris-NTA compounds (such as #12615 and #12617), it can be used for multicolor analysis of polyhistidine-tagged proteins. 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. The transition metal ions (e.g., Ni<sup>2+</sup>)-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.