

Tyramide Sulfo-NHS Ester [Sulpho SHPP]

Catalog Number: 11044

Unit Size: 25 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	365.29
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Tyramide Sulfo-NHS Ester is an excellent building block to develop TSA (tyramide) probes. It is also called as Water-Soluble Bolton-Hunter Reagent (Sulfo-SHPP) used for preparing tyrosine-like conjugates for increasing the yield of subsequent iodination. The reagent is ideal for iodinating proteins that do not contain tyrosine residues or tyrosine residues with limited accessibility to iodination. Labeling with Sulfo-SHPP also can help to preserve tyrosine residues that are essential for protein function or immunogenicity. It contains an N-hydroxysuccinimide (NHS) ester reactive group that react efficiently with primary amino groups (-NH₂) in pH 7-9 buffers to form stable amide bonds. The sulfonated NHS ester of Sulfo-SHPP provides increased reagent solubility in comparison to the non-sulfonated form of Bolton-Hunter Reagent.