

Acridinium Biotin Conjugate

Catalog Number: 26050

Unit Size: 1 mg

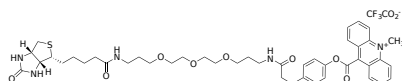
Product Details

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

Chemical Properties

Appearance	Solid off- white
Molecular Weight	928.03
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Acridinium biotin is a conjugate of acridinium ester and biotin. It is a chemiluminescent label that can be used to detect antibodies or antigens in a wide range of immunoassays. For example, it may be used to label antibodies or antigens that can be detected by streptavidin-coated magnetic beads in a chemiluminescent immunoassay system. Since biotin is a small molecule that can bind to streptavidin with high affinity and specificity, the acridinium biotin reagent can be used in a variety of immunological applications in combination of numerous commercially available streptavidin conjugates.