

Cy7 BCN

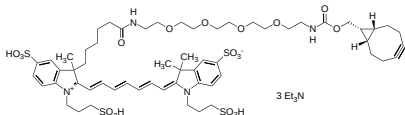
Catalog Number: 70544

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	1583.13
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	756 nm
Emission Wavelength	779 nm

Applications

Cy7 BCN is a clickable derivative of Cy7, a widely used near-infrared fluorescent dye for labeling peptides, oligonucleotides, and other biomolecules. Cy7 is characterized by a strong fluorescence intensity, high photostability, and pH-independent emission across a broad range (pH 4–10). To improve conjugation performance, Cy7-BCN incorporates a PEG spacer, which reduces steric hindrance and minimizes potential interference with target binding sites. This design maximizes conjugation efficiency while preserving the biological activity of the resulting conjugate. The bicyclononyne (BCN) moiety enables strain-promoted azide-alkyne cycloaddition (SPAAC) with azido groups, forming stable triazole linkages under catalyst-free conditions. In addition, unlike dibenzocyclooctyne (DBCO), BCN also reacts efficiently with tetrazines through an inverse electron-demand Diels–Alder (IEDDA) reaction. This reaction is rapid, selective, and bioorthogonal, allowing labeling of biomolecules under physiological conditions without the need for metal catalysts or disruption of native biological processes.