

Diazepam BSA Conjugate

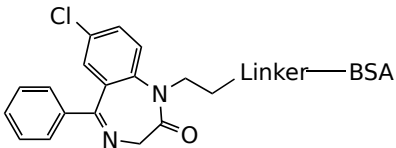
Catalog Number: 50665

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Diazepam BSA Conjugate is a bioconjugate in which diazepam is covalently linked to bovine serum albumin (BSA) to enhance its immunogenicity for use in immunoassay development. The conjugation is typically performed via active ester or linker-mediated chemistry to maintain the structural integrity of the diazepam pharmacophore while achieving efficient and stable attachment to the carrier protein. This conjugate is commonly employed as an immunogen for generating diazepam-specific polyclonal and monoclonal antibodies, as well as a coating antigen in competitive ELISA, lateral flow assays, and other immunodetection platforms. The high antigen density on BSA improves assay sensitivity and specificity, facilitating the reliable detection of diazepam and structurally related benzodiazepines in biological samples. This reagent is intended for research use only and is not approved for diagnostic or therapeutic applications.