

Diazepam Biotin Conjugate

Catalog Number: 50664

Unit Size: 1 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

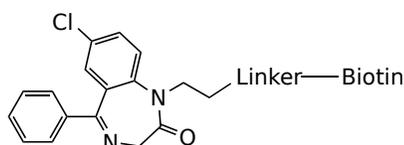
Chemical Properties

Appearance Solid

Molecular Weight 1 mg

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

Applications

Diazepam Biotin Conjugate is a structurally defined, biotinylated analog of diazepam engineered for affinity-based applications in neuropharmacological and biochemical research. The molecule features a stable covalent linkage between diazepam and biotin via a suitable spacer, designed to preserve the pharmacologically active benzodiazepine core while enabling high-affinity interaction with streptavidin, avidin, or neutravidin matrices. This dual functionality allows for the selective capture, enrichment, or immobilization of GABA_A receptor subunits or diazepam-binding proteins in complex biological samples. The conjugate is particularly useful in receptor-ligand binding assays, affinity purification, and target deconvolution workflows, where it facilitates sensitive and specific interrogation of GABAergic signaling pathways. Its compatibility with biotin-binding platforms also supports its use in biosensor development and high-throughput screening for drug discovery.