

Diazepam NHS ester

Catalog Number: 50660

Unit Size: 1 mg

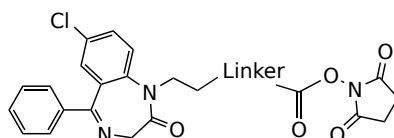
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	481.93
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Diazepam NHS ester is a chemically activated derivative of diazepam incorporating an N-hydroxysuccinimide (NHS) ester functional group for targeted bioconjugation. The NHS ester reacts selectively with primary amines under physiological to slightly basic conditions (pH 7.0–8.5), enabling the formation of stable amide bonds with proteins, peptides, and other amine-containing biomolecules. The diazepam scaffold remains intact, preserving its affinity for GABA_A receptors and related pharmacological targets. This reagent is well-suited for the development of diazepam-functionalized affinity ligands, targeted delivery systems, and molecular probes for receptor-binding assays and imaging applications. Its compatibility with aqueous systems and high coupling efficiency make Diazepam NHS ester a useful tool in neuropharmacological research, investigating benzodiazepine mechanisms, receptor-ligand interaction studies, and the development of functionalized delivery platforms or diagnostic probes.