

Medetomidine BSA conjugate

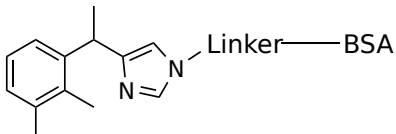
Catalog Number: 50655

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

The Medetomidine-BSA conjugate is a synthetic immunogen designed for the generation of antibodies against medetomidine, a selective α_2 -adrenergic receptor agonist commonly used in veterinary anesthesia and research. Due to its low molecular weight and limited immunogenicity, medetomidine must be conjugated to a carrier protein to elicit a robust immune response. In this construct, medetomidine is covalently linked to bovine serum albumin (BSA) via a chemically defined linker, ensuring stable attachment while preserving the pharmacologically relevant epitope. This conjugate is suitable for applications in antibody production and the development of immunoassays such as ELISA, competitive binding assays, and lateral flow devices for the qualitative or quantitative detection of medetomidine or structurally related compounds. This reagent is intended for research use only and is not approved for diagnostic or therapeutic applications.