

Xylazine-Biotin Conjugate

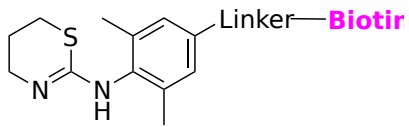
Catalog Number: 50615

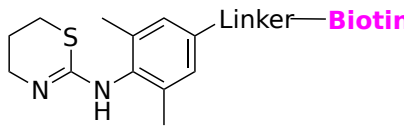
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid off-white
Molecular Weight	778.90
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Xylazine-Biotin conjugate is an excellent probe for developing assays and tests for detecting Xylazine. AAT Bioquest offers the largest collection of streptavidin conjugates that can be used with Xylazine-Biotin for detecting Xylazine. Xylazine is a non-opioid sedative or tranquilizer. It is not approved for use in people. It is commonly used in veterinary medicine for sedation, anesthesia, muscle relaxation, and analgesia in animals such as horses, cattle, and other mammals. In veterinary anesthesia, it is often combined with ketamine. Veterinarians also use xylazine as an emetic, especially in cats.