

Xylazine-HRP Conjugate

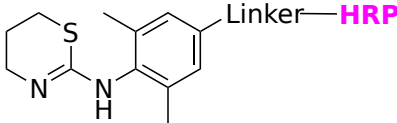
Catalog Number: 50622

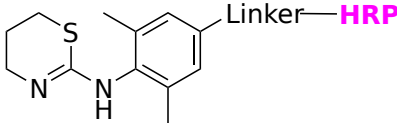
Unit Size: 1 mg

Product Details

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

Chemical Properties

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



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Xylazine-HRP conjugate is an excellent probe for developing ELISA assays and tests 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.