

Xylazine NHS Ester

Catalog Number: 50610

Unit Size: 1 mg

Product Details

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

Expiration Date 12 months upon receiving

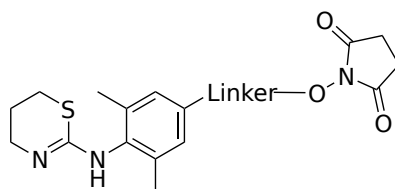
Chemical Properties

Appearance Solid off-white

Molecular Weight 361.42

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

Applications

Xylazine NHS ester is an excellent amino-reactive building block for developing 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.