

Xylazine-Fluorescein Conjugate

Catalog Number: 50616

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	814.74
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Xylazine-Fluorescein conjugate is an excellent fluorescent probe for developing assays and tests for detecting Xylazine. AAT Bioquest offers the largest collection of dyes that can be used with Xylazine-Fluorescein conjugate for multicolor analysis of 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.