

# Xylazine-Fluorescein Conjugate

Catalog Number: 50616

Unit Size: 1 mg

## Product Details

---

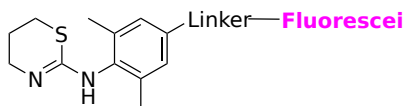
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

## Chemical Properties

---

Appearance	Solid orange
Molecular Weight	814.74
Soluble In	DMSO

Chemical Structure



## 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.