

# Xylazine Amine

Catalog Number: 50611

Unit Size: 1 mg

## Product Details

---

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

Expiration Date 24 months upon receiving

## Chemical Properties

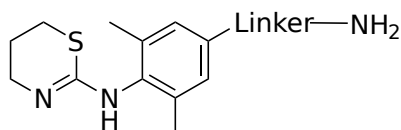
---

Appearance Solid off-white

Molecular Weight 534.48

Soluble In DMSO

Chemical Structure



## Spectral Properties

---

Excitation Wavelength N/A

Emission Wavelength N/A

## Applications

---

Xylazine amine is an excellent 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.