

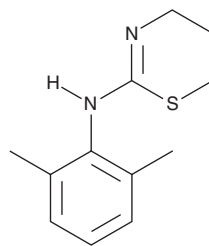
PRODUCT INFORMATION



Xylazine

Item No. 22641

CAS Registry No.: 7361-61-7
Formal Name: N-(2,6-dimethylphenyl)-5,6-dihydro-4H-1,3-thiazin-2-amine
MF: C₁₂H₁₆N₂S
FW: 220.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Xylazine (Item No. 22641) is an analytical reference standard categorized as an anesthetic.¹ It has been used as an adulterant in drugs of abuse and misused, leading to non-fatal and fatal intoxications.² This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 14113).

References

1. Richardson, C.A. and Flecknell, P.A. Anaesthesia and post-operative analgesia following experimental surgery in laboratory rodents: Are we making progress? *Altem. Lab. Anim.* **33(2)**, 119-127 (2005).
2. Ruiz-Colon, K., Chavez-Arias, C., Diaz-Alcala, J.E., *et al.* Xylazine intoxication in humans and its importance as an emerging adulterant in abused drugs: A comprehensive review of the literature. *Forensic Sci. Int.* **240**, 1-8 (2014).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/15/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM