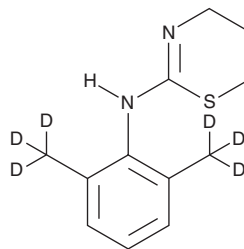


PRODUCT INFORMATION



Xylazine-d₆ Item No. 38353

CAS Registry No.: 1228182-53-3
Formal Name: N-(2,6-bis(methyl-d₃)phenyl)-5,6-dihydro-4H-1,3-thiazin-2-amine
MF: C₁₂H₁₀D₆N₂S
FW: 226.4
Chemical Purity: ≥98% (Xylazine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

Xylazine-d₆ (Item No. 38353) is intended for use as an internal standard for the quantification of xylazine (Item No. 22641) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Xylazine is categorized as an anesthetic.¹ Xylazine 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 research and forensic applications.

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, 02/26/2025

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