

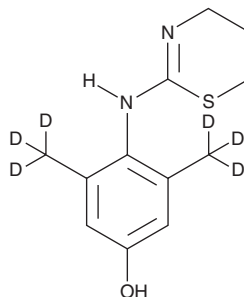
PRODUCT INFORMATION



4-hydroxy Xylazine-d₆ (CRM)

Item No. 39571

Formal Name: 4-((5,6-dihydro-4H-1,3-thiazin-2-yl)amino)-3,5-bis(methyl-d₃)phenol
Synonyms: p-hydroxy Xylazine-d₆, para-hydroxy Xylazine-d₆
MF: C₁₂H₁₀D₆N₂OS
FW: 242.4
Supplied as: A 100 µg/ml solution in methanol
Storage: -20°C
Stability: ≥2 years



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

Description

4-hydroxy Xylazine-d₆ (CRM) (Item No. 39571) is intended for use as an internal standard for the quantification of 4-hydroxy xylazine (Item No. 39607) 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).

4-hydroxy Xylazine is categorized as an anesthetic metabolite.¹ It is a metabolite of xylazine (Item Nos. 37854 | 22641) and has been used as a marker for xylazine doping in horses.^{2,3} This product is intended for research and forensic applications.

References

1. Mutlib, A.E., Chui, Y.C., Young, L.M., *et al.* Characterization of metabolites of xylazine produced *in vivo* and *in vitro* by LC/MS/MS and by GC/MS. *Drug Metab. Dispos.* **20(6)**, 840-848 (1992).
2. Spyrodaki, M.-H., Lyris, E., Georgoulakis, I., *et al.* Determination of xylazine and its metabolites by GC-MS in equine urine for doping analysis. *J. Pharm. Biomed. Anal.* **35(1)**, 107-116 (2004).
3. Kwok, W.H., Choi, T.L.S., Tsoi, Y.Y.K., *et al.* Screening of over 100 drugs in horse urine using automated on-line solid-phase extraction coupled to liquid chromatography-high resolution mass spectrometry for doping control. *J. Chromatogr. A* **1490**, 89-101 (2017).

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, 01/03/2025

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