

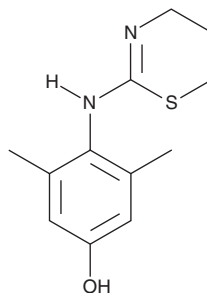
PRODUCT INFORMATION



4-hydroxy Xylazine

Item No. 22920

CAS Registry No.: 145356-32-7
Formal Name: 4-[(5,6-dihydro-4H-1,3-thiazin-2-yl)amino]-3,5-dimethyl-phenol
Synonyms: *p*-hydroxy Xylazine, *para*-hydroxy Xylazine
MF: C₁₂H₁₆N₂OS
FW: 236.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

4-hydroxy Xylazine is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy xylazine in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Xylazine is slightly soluble in methanol and DMSO.

Description

4-hydroxy Xylazine is a metabolite of xylazine (Item No. 14113).¹ Xylazine is an agonist of α_2 -adrenergic receptors used for sedation, anesthesia, and analgesia in non-human mammals.²⁻⁴ 4-hydroxy Xylazine has been used as a marker for xylazine doping in MS analysis of equine urine.¹

References

1. 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).
2. Virtanen, R., Savola, J.M., Saano, V., *et al.* Characterization of the selectivity, specificity and potency of medetomidine as an α_2 -adrenoceptor agonist. *Eur. J. Pharmacol.* **150(1-2)**, 9-14 (1988).
3. 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).
4. Vallverde, A. Alpha-2 agonists as pain therapy in horses. *Vet. Clin. North Am. Equine Pract.* **26(3)**, 515-532 (2010).

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, 09/29/2022

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