

PRODUCT INFORMATION



4-hydroxy Xylazine O-Glucuronide (lithium salt)

Item No. 38355

Formal Name: (2S,3S,4S,5R,6S)-6-((4-((5,6-dihydro-4H-1,3-thiazin-2-yl)amino)-3,5-dimethylphenoxy)-3,4,5-trihydroxytetrahydro-2H-pyran-2-carboxylate, monolithium salt

Synonyms: p-hydroxy Xylazine O-Glucuronide, para-hydroxy Xylazine O-Glucuronide

MF: $C_{18}H_{23}N_2O_7S \cdot Li$

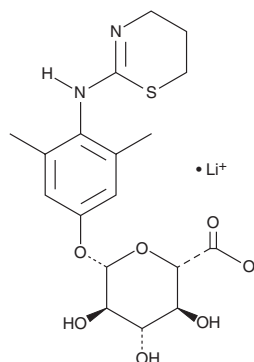
FW: 418.4

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

4-hydroxy Xylazine O-glucuronide (lithium salt) is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy xylazine O-glucuronide (lithium salt) in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Xylazine O-glucuronide (lithium salt) is soluble in DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

4-hydroxy Xylazine O-glucuronide is a metabolite of the α_2 -adrenergic receptor agonist xylazine (Item No. 14113).¹ It is formed from xylazine via the intermediate metabolite 4-hydroxy xylazine (Item No. 22920).

Reference

1. Matos, R.R., Martucci, M.E.P., de Anselmo, C.S., *et al.* Pharmacokinetic study of xylazine in a zebrafish water tank, a human-like surrogate, by liquid chromatography Q-Orbitrap mass spectrometry. *Forensic Toxicol.* **38**, 108-121 (2020).

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, 03/30/2023

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