

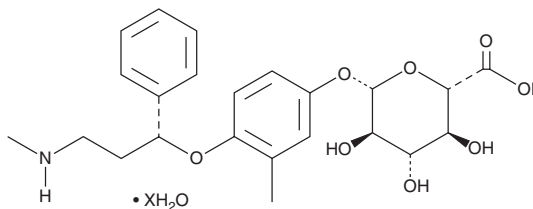
PRODUCT INFORMATION



4'-hydroxy Atomoxetine Glucuronide (hydrate)

Item No. 34822

Formal Name: 3-methyl-4-[(1R)-3-(methylamino)-1-phenylpropoxy]phenyl β-D-glucopyranosiduronic acid, hydrate
MF: C₂₃H₂₉NO₈ • XH₂O
FW: 447.5
Purity: ≥95%
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.

Laboratory Procedures

4'-hydroxy Atomoxetine glucuronide (hydrate) is supplied as a solid. A stock solution may be made by dissolving the 4'-hydroxy atomoxetine glucuronide (hydrate) in the solvent of choice, which should be purged with an inert gas. 4'-hydroxy Atomoxetine glucuronide (hydrate) is slightly soluble in methanol (sonicated).

4'-hydroxy Atomoxetine glucuronide (hydrate) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

4'-hydroxy Atomoxetine glucuronide is a metabolite of the norepinephrine transporter (NET) inhibitor atomoxetine (Item No. 22248).¹ It is formed from atomoxetine by glucuronidation of the intermediate metabolite 4-hydroxy atomoxetine.²

References

1. Todor, I., Popa, A., Neag, M., *et al.* Evaluation of a potential metabolism-mediated drug-drug interaction between atomoxetine and bupropion in healthy volunteers. *J. Pharm. Pharm. Sci.* **19(2)**, 198-207 (2016).
2. Sauer, J.-M., Ring, B.J., and Witcher, J.W. Clinical pharmacokinetics of atomoxetine. *Clin. Pharmacokinet.* **44(6)**, 571-590 (2005).

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, 12/06/2022

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