

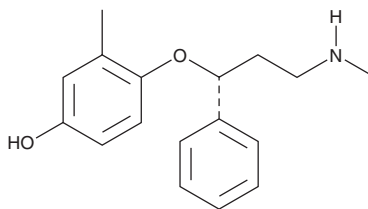
PRODUCT INFORMATION



4-hydroxy Atomoxetine

Item No. 34821

CAS Registry No.: 435293-66-6
Formal Name: 3-methyl-4-[(1R)-3-(methylamino)-1-phenylpropoxy]-phenol
MF: $C_{17}H_{21}NO_2$
FW: 271.4
Purity: $\geq 98\%$
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 is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy atomoxetine in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Atomoxetine is soluble in the organic solvent DMSO.

Description

4-hydroxy Atomoxetine is a metabolite of the norepinephrine transporter (NET) inhibitor atomoxetine (Item No. 22248).¹ It is formed from atomoxetine by the cytochrome P450 (CYP) isoform CYP2D6 in human liver microsomes.

Reference

1. 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, 10/17/2022

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