

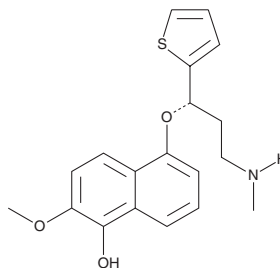
PRODUCT INFORMATION



5-hydroxy-6-methoxy (S)-Duloxetine

Item No. 28127

CAS Registry No.: 741693-79-8
Formal Name: 2-methoxy-5-[(1S)-3-(methylamino)-1-(2-thienyl)propoxy]-1-naphthalenol
MF: $C_{19}H_{21}NO_3S$
FW: 343.4
Purity: $\geq 95\%$
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

5-hydroxy-6-methoxy (S)-Duloxetine is supplied as a solid. A stock solution may be made by dissolving the 5-hydroxy-6-methoxy (S)-duloxetine in the solvent of choice, which should be purged with an inert gas. 5-hydroxy-6-methoxy (S)-Duloxetine is slightly soluble in chloroform and methanol.

Description

5-hydroxy-6-methoxy (S)-Duloxetine is a metabolite of (S)-duloxetine (Item No. 14317).^{1,2} It is formed from (S)-duloxetine via a 5- or 6-hydroxy duloxetine intermediate, which is formed by the cytochrome P450 (CYP) isoforms CYP1A2 and CYP2D6, and a catechol duloxetine intermediate.² 5-hydroxy-6-methoxy (S)-Duloxetine binds to the serotonin (5-HT), norepinephrine, and dopamine transporters with K_i values of 266, 920, and 2,814 nM, respectively.

References

1. Kuo, F., Gillespie, T.A., Kulanthaivel, P., *et al.* Synthesis and biological activity of some known and putative duloxetine metabolites. *Bioorg. Med. Chem.* **14**(13), 3481-3486 (2004).
2. Knadler, M.P., Lobo, E., Chappell, J., *et al.* Duloxetine: Clinical pharmacokinetics and drug interactions. *Clin. Pharmacokinet.* **50**(5), 281-294 (2011).

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, 10/24/2022

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