

PRODUCT INFORMATION

4-hydroxy (S)-Duloxetine β-D-Glucuronide (sodium salt)

Item No. 27823

Formal Name: 4-[(1S)-3-(methylamino)-1-(2-thienyl)propoxy]-1-naphthalenyl, β-D-glucopyranosiduronic acid, monosodium salt

Synonym: LY550408

MF: C₂₄H₂₆NO₈S • Na

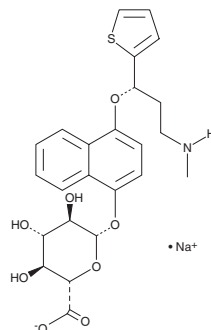
FW: 511.5

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 (S)-Duloxetine β-D-glucuronide (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy (S)-duloxetine β-D-glucuronide (sodium salt) in the solvent of choice, which should be purged with an inert gas. 4-hydroxy (S)-Duloxetine β-D-glucuronide (sodium salt) is slightly soluble in methanol and DMSO.

Description

4-hydroxy (S)-Duloxetine β-D-glucuronide is an inactive metabolite of the serotonin and norepinephrine reuptake inhibitor (S)-duloxetine (Item No. 14317).¹ It is formed via metabolism of (S)-duloxetine by the cytochrome P450 (CYP) isoforms CYP1A2 and CYP2D6 followed by glucuronidation.

Reference

1. 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, 12/05/2022

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