

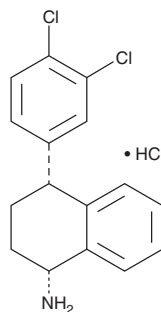
PRODUCT INFORMATION



(1R,4R)-N-desmethyl Sertraline (hydrochloride)

Item No. 35741

CAS Registry No.: 675126-09-7
Formal Name: (1R,4R)-4-(3,4-dichlorophenyl)-1,2,3,4-tetrahydro-1-naphthalenamine, monohydrochloride
Synonym: (1R,4R)-N-demethyl Sertraline
MF: C₁₆H₁₅Cl₂N • HCl
FW: 328.7
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

(1R,4R)-N-desmethyl Sertraline (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the (1R,4R)-N-desmethyl sertraline (hydrochloride) in the solvent of choice, which should be purged with an inert gas. (1R,4R)-N-desmethyl Sertraline (hydrochloride) is slightly soluble in methanol and DMSO.

Description

(1R,4R)-N-desmethyl Sertraline is a metabolite of the selective serotonin reuptake inhibitor (SSRI) sertraline (Item No. 14839).¹

Reference

1. Obach, R.S., Cox, L.M., and Tremaine, L.M. Sertraline is metabolized by multiple cytochrome P450 enzymes, monoamine oxidases, and glucuronyl transferases in human: An in vitro study. *Drug Metab. Dispos.* **33(2)**, 262-270 (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, 02/13/2024

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