

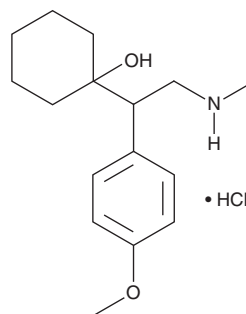
PRODUCT INFORMATION



(±)-N-desmethyl Venlafaxine (hydrochloride)

Item No. 35403

CAS Registry No.: 93413-90-2
Formal Name: 1-[1-(4-methoxyphenyl)-2-(methylamino)ethyl]-cyclohexanol, monohydrochloride
Synonym: Wy 45494
MF: C₁₆H₂₅NO₂ • HCl
FW: 299.8
Purity: ≥95%
UV/Vis.: λ_{max}: 226 nm
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

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

Description

(±)-N-desmethyl Venlafaxine is a minor active metabolite of the selective norepinephrine and serotonin reuptake inhibitor (SNRI) venlafaxine.¹ It is formed from venlafaxine by the cytochrome P450 (CYP) isoform CYP3A4.² (±)-N-desmethyl Venlafaxine inhibits norepinephrine and serotonin reuptake in rat synaptosome preparations (IC₅₀s = 4.7 and 1.6 μM, respectively).¹ *In vivo*, (±)-N-desmethyl venlafaxine reverses reserpine-induced hypothermia in mice with a minimum effective dose (MED) of 10 mg/kg.

References

1. Muth, E.A., Moyer, J.A., Haskins, J.T., *et al.* Biochemical, neurophysiological, and behavioral effects of Wy-45,233 and other identified metabolites of the antidepressant venlafaxine. *Drug Dev. Res.* **23(2)**, 191-199 (1991).
2. Otton, S.V., Ball, S.E., Cheung, S.W., *et al.* Venlafaxine oxidation *in vitro* is catalysed by CYP2D6. *Br. J. Clin. Pharmacol.* **41(2)**, 149-156 (1996).

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, 11/07/2022

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