

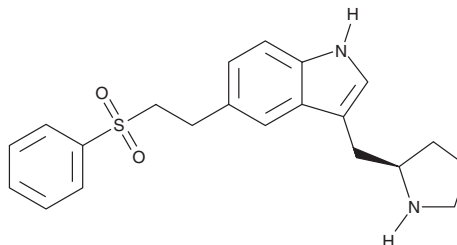
PRODUCT INFORMATION



N-desmethyl Eletriptan

Item No. 34302

CAS Registry No.: 153525-55-4
Formal Name: (R)-5-[2-(phenylsulfonyl)ethyl]-3-(2-pyrrolidinylmethyl)-1H-indole
MF: $C_{21}H_{24}N_2O_2S$
FW: 368.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

N-desmethyl Eletriptan is supplied as a solid. A stock solution may be made by dissolving the N-desmethyl eletriptan in the solvent of choice, which should be purged with an inert gas. N-desmethyl Eletriptan is soluble in organic solvents such as acetonitrile, DMSO, and methanol.

Description

N-desmethyl Eletriptan is a metabolite of eletriptan (Item No. 20048).¹ It is formed from eletriptan primarily by the cytochrome P450 (CYP) isoform CYP3A4 in human liver microsomes.

Reference

1. Evans, D.C., O'Connor, D., Lake, B.G., *et al.* Eletriptan metabolism by human hepatic CYP450 enzymes and transport by human P-glycoprotein. *Drug Metab. Dispos.* **31**(7), 861-869 (2003).

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, 11/18/2022

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