

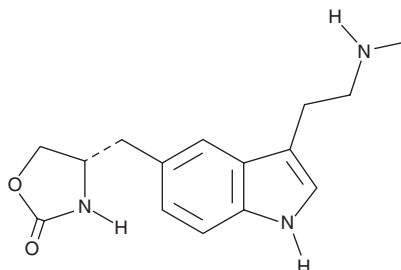
PRODUCT INFORMATION



N-desmethyl Zolmitriptan

Item No. 28058

CAS Registry No.: 139264-35-0
Formal Name: 4(S)-[[[3-[2-(methylamino)ethyl]-1H-indol-5-yl]methyl]-2-oxazolidinone
Synonyms: 183C91, DZT
MF: C₁₅H₁₉N₃O₂
FW: 273.3
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

N-desmethyl Zolmitriptan is supplied as a solid. A stock solution may be made by dissolving the N-desmethyl zolmitriptan in the solvent of choice, which should be purged with an inert gas. N-desmethyl Zolmitriptan is soluble in the organic solvent methanol.

Description

N-desmethyl Zolmitriptan is an active metabolite of the serotonin (5-HT) receptor subtype 5-HT_{1B} and 5-HT_{1D} agonist zolmitriptan (Item No. 15871).¹ N-desmethyl Zolmitriptan is an agonist of 5-HT_{1B} receptors that induces contraction of isolated human cerebral arteries (EC₅₀ = 100 nM).²

References

1. Patel, B., Suhagia, B.N., Jangid, A.G., *et al.* Systematic evaluation of matrix effect and cross-talk-free method for simultaneous determination of zolmitriptan and N-desmethyl zolmitriptan in human plasma: A sensitive LC-MS/MS method validation and its application to a clinical pharmacokinetic study. *Biomed. Chromatogr.* **30(3)**, 447-458 (2016).
2. Nilsson, T., Longmore, J., Shaw, D., *et al.* Contractile 5-HT_{1B} receptors in human cerebral arteries: Pharmacological characterization and localization with immunocytochemistry. *Br. J. Pharmacol.* **128(6)**, 1133-1140 (1999).

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, 11/01/2022

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