

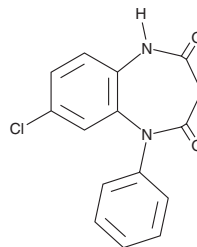
PRODUCT INFORMATION



N-Desmethylclobazam

Item No. 18564

CAS Registry No.: 22316-55-8
Formal Name: 8-chloro-1-phenyl-1H-1,5-benzodiazepine-2,4(3H,5H)-dione
Synonym: nor-Clobazam
MF: C₁₅H₁₁ClN₂O₂
FW: 286.7
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

N-Desmethylclobazam (Item No. 18564) is an analytical reference material that is structurally categorized as a benzodiazepine. It is a major active metabolite of clobazam, produced primarily by the action of cytochrome P450 isoform 3A4.¹ N-Desmethylclobazam positively modulates GABA_A receptors (EC₅₀s range from 98 to 151 μM, depending on subunit configuration).² This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. de Leon, J., Spina, E., and Diaz, F.J. Clobazam therapeutic drug monitoring: A comprehensive review of the literature with proposals to improve future studies. *Ther. Drug Monit.* **35(1)**, 30-47 (2013).
2. Hammer, H., Ebert, B., Jensen, H.S., *et al.* Functional characterization of the 1,5-benzodiazepine clobazam and its major active metabolite N-desmethylclobazam at human GABA_A receptors expressed in *Xenopus laevis* oocytes. *PLoS One* **10(3)**, (2015).

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, 08/13/2018

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