

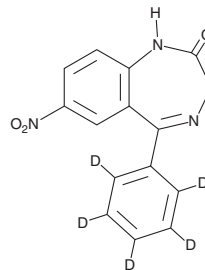
PRODUCT INFORMATION



Nitrazepam-d₅ (exempt preparation)

Item No. 38139

CAS Registry No.: 136765-45-2
Formal Name: 1,3-dihydro-7-nitro-5-(phenyl-d₅)-2H-1,4-benzodiazepin-2-one
MF: C₁₅H₆D₅N₃O₃
FW: 286.3
Purity: ≥98%
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥3 years



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

Description

Nitrazepam-d₅ (exempt preparation) (Item No. 38139) is intended for use as an internal standard for the quantification of nitrazepam (Item Nos. ISO60201 | 16196) by GC-or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Nitrazepam is categorized as a benzodiazepine. It is also categorized as a sedative and has been abused recreationally.^{1,2} Nitrazepam-d₅ is regulated as a Schedule IV compound in the United States. Nitrazepam-d₅ (exempt preparation) (Item No. 38139) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Yakushiji, T., Fukuda, T., Oyama, Y., *et al.* Effects of benzodiazepines and non-benzodiazepine compounds on the GABA-induced response in frog isolated sensory neurones. *Br. J. Pharmacol.* **98**(3), 735-740 (1989).
2. Lyphout, C., Yates, C., Margolin, Z.R., *et al.* Presentations to the emergency department with non-medical use of benzodiazepines and Z-drugs: Profiling and relation to sales data. *Eur. J. Clin. Pharmacol.* **75**(1), 77-85 (2019).

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, 01/10/2025

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