

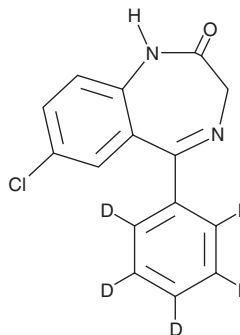
# PRODUCT INFORMATION



## Nordiazepam-d<sub>5</sub>

Item No. 15831

**CAS Registry No.:** 65891-80-7  
**Formal Name:** 7-chloro-1,3-dihydro-5-(phenyl-d<sub>5</sub>)-2H-1,4-benzodiazepin-2-one  
**Synonyms:** Dealkylprazepam-d<sub>5</sub>, Demethyldiazepam-d<sub>5</sub>, Desmethyldiazepam-d<sub>5</sub>, Nordazepam-d<sub>5</sub>, Norprazepam-d<sub>5</sub>  
**MF:** C<sub>15</sub>H<sub>6</sub>ClD<sub>5</sub>N<sub>2</sub>O  
**FW:** 275.8  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

Nordiazepam-d<sub>5</sub> (Item No. 15831) is an analytical reference material intended for use as an internal standard for the quantification of nordiazepam (Item No. ISO60174) 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).

Nordiazepam is categorized as a benzodiazepine. It is an active metabolite of diazepam (Item Nos. 15887 | ISO60177) and has sedative properties.<sup>1,2</sup> Nordiazepam-d<sub>5</sub> is regulated as a Schedule IV compound in the United States. 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 for reference materials.

### References

1. Calcaterra, N.E. and Barrow, J.C. Classics in chemical neuroscience: Diazepam (valium). *ACS Chem. Neurosci.* **5(4)**, 253-260 (2014).
2. Traversa, U., de Angelis, L., and Vertua, R. On the hypnogenic and anticonvulsant activities of demethyldiazepam and chlordemethyldiazepam: Time-effect relations. *J. Pharm. Pharmacol.* **29(8)**, 504-506 (1977).

#### 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, 08/27/2019

#### 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