

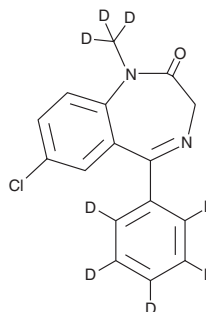
PRODUCT INFORMATION



Diazepam-d₈ (exempt preparation)

Item No. 23786

CAS Registry No.: 83056-50-2
Formal Name: 7-chloro-1,3-dihydro-1-(methyl-d₃)-5-(phenyl-d₅)-2H-1,4-benzodiazepin-2-one
MF: C₁₆H₅ClD₈N₂O
FW: 292.8
Chemical Purity: ≥98% (Diazepam)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

Diazepam-d₈ (exempt preparation) (Item No. 23786) is intended for use as an internal standard for the quantification of diazepam (Item No. ISO60177) 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).

Diazepam is categorized as a benzodiazepine.^{1,2} Diazepam is regulated as a Schedule IV compound in the United States. Diazepam-d₈ (exempt preparation) (Item No. 23786) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Rickels, K., Csanalosi, I., Greisman, P., *et al.* Ketazolam and diazepam in anxiety: A controlled study. *J. Clin. Pharmacol.* **20**(10), 581-589 (1980).
2. Birnir, B., Eghbali, M., Everitt, A.B., *et al.* Bicuculline, pentobarbital and diazepam modulate spontaneous GABA_A channels in rat hippocampal neurons. *Br. J. Pharmacol.* **131**(4), 695-704 (2000).

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, 02/13/2025

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