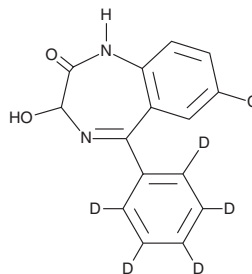


PRODUCT INFORMATION



Oxazepam-d₅ Item No. 15844

CAS Registry No.: 65854-78-6
Formal Name: 7-chloro-1,3-dihydro-3-hydroxy-5-(phenyl-d₅)-2H-1,4-benzodiazepin-2-one
Synonym: N-Desmethyletemazepam-d₅
MF: C₁₅H₆ClD₅N₂O₂
FW: 291.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

Oxazepam-d₅ is intended for use as an internal standard for the quantification of oxazepam (Item No. ISO60173) 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).

Oxazepam-d₅ is an analytical reference material that is structurally categorized as a benzodiazepine. Oxazepam is an active metabolite of diazepam (Item No. ISO60177), prazepam, and temazepam (Item No. ISO60182) that can positively modulate GABA_A receptors, which results in inhibitory effects in the central nervous system.¹ This product is intended for forensic and research purposes.

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.

Reference

1. Skerritt, J.H. and Johnston, G.A.R. Enhancement of GABA binding by benzodiazepines and related anxiolytics. *Eur. J. Pharmacol.* **89(3-4)**, 193-198 (1983).

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