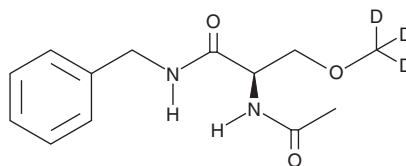


PRODUCT INFORMATION



Lacosamide-d₃ Item No. 25426

CAS Registry No.: 1217689-95-6
Formal Name: 2R-(acetylamino)-3-(methoxy-d₃)-N-(phenylmethyl)-propanamide
MF: C₁₃H₁₅D₃N₂O₃
FW: 253.3
Chemical Purity: ≥98% (Lacosamide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥5 years



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

Description

Lacosamide-d₃ (Item No. 25426) is an analytical reference standard intended for use as an internal standard for the quantification of lacosamide 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).

Lacosamide is categorized as an anticonvulsant.¹ Formulations containing lacosamide have induced euphoria, auditory and visual hallucinations, and psychosis in humans but have a low potential for abuse.^{1,2} Lacosamide-d₃ is regulated as a Schedule V compound in the United States. This product is intended for research and forensic applications.

References

1. Hawkins, K.L. and Gidal, B.E. When adverse effects are seen as desirable: Abuse potential of the newer generation antiepileptic drugs. *Epilepsy Behav.* **77**, 62-72 (2017).
2. Schoedel, K.A., Andreas, J.O., Doty, P., *et al.* Randomized, double-blind, placebo- and active comparator-controlled crossover study evaluating the abuse potential of the antiepileptic drug lacosamide in healthy recreational drug users. *J. Clin. Psychopharmacol.* **37(6)**, 675-683 (2017).

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, 07/06/2023

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