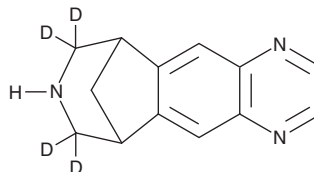


PRODUCT INFORMATION



Varenicline-d₄ Item No. 26102

CAS Registry No.: 2183239-01-0
Formal Name: 7,8,9,10-tetrahydro-6H-6,10-methanoazepino[4,5-g]quinoxaline-7,7,9,9-d₄
MF: C₁₃H₉D₄N₃
FW: 215.3
Chemical Purity: ≥98% (Varenicline)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Varenicline-d₄ is intended for use as an internal standard for the quantification of varenicline (Item No. 15030) 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).

Varenicline-d₄ is supplied as a solid. A stock solution may be made by dissolving the varenicline-d₄ in water. The solubility of varenicline-d₄ in water is approximately 0.0877 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Varenicline is a full agonist of α7 subunit-containing neuronal nicotinic acetylcholine receptors (nAChRs; EC₅₀ = 18 μM) and a partial agonist of α4β2 subunit-containing nAChRs (EC₅₀ = 2.3 μM).¹ Formulations containing varenicline have been used as aids in smoking cessation.

Reference

1. Mihalak, K.B., Carroll, F.I., and Luetje, C.W. Varenicline is a partial agonist at α4β2 and a full agonist at α7 neuronal nicotinic receptors. *Mol. Pharm.* **70**(3), 801-805 (2006).

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, 12/16/2022

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