

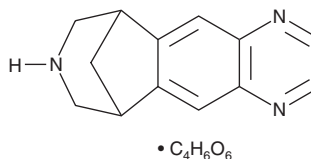
PRODUCT INFORMATION



Varenicline (tartrate)

Item No. 15030

CAS Registry No.: 375815-87-5
Formal Name: 7,8,9,10-tetrahydro-6,10-methano-6H-pyrazino[2,3-h][3]benzazepine-(2R,3R)-dihydroxybutanedioate
Synonym: CP 526,555-18
MF: C₁₃H₁₃N₃ • C₄H₆O₆
FW: 361.4
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 236, 320 nm
Supplied as: A crystalline 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 (tartrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the varenicline (tartrate) in the solvent of choice, which should be purged with an inert gas. Varenicline (tartrate) is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of varenicline (tartrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of varenicline (tartrate) in PBS (pH 7.2) is approximately 10 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, 12/22/2022

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