

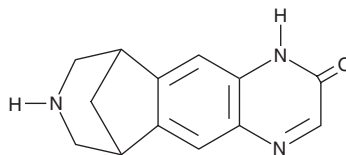
# PRODUCT INFORMATION



## Hydroxy Varenicline

Item No. 22388

**CAS Registry No.:** 357424-21-6  
**Formal Name:** 1,6,7,8,9,10-hexahydro-6,10-methano-2H-pyrazino[2,3-h][3]benzazepin-2-one  
**Synonym:** 2-Hydroxyvarenicline  
**MF:** C<sub>13</sub>H<sub>13</sub>N<sub>3</sub>O  
**FW:** 227.3  
**Purity:** ≥95%  
**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

Hydroxy varenicline is supplied as a solid. A stock solution may be made by dissolving the hydroxy varenicline in the solvent of choice, which should be purged with an inert gas. Hydroxy varenicline is soluble in a 1:1 solution of acetonitrile:methanol.

### Description

Hydroxy varenicline is a metabolite of the nicotinic acetylcholine receptor (nAChR) agonist varenicline (Item No. 15030).<sup>1</sup>

### Reference

1. Obach, R.S., Reed-Hagen, A.E., Krueger, S.S., *et al.* Metabolism and disposition of varenicline, a selective α4β2 acetylcholine receptor partial agonist, in vivo and in vitro. *Drug. Metab. Dispos.* **34**(1), 121-130 (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, 12/12/2022

#### 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](http://WWW.CAYMANCHEM.COM)