

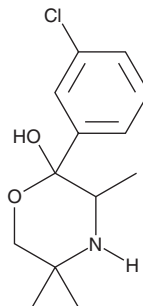
PRODUCT INFORMATION



hydroxy Bupropion

Item No. 9000522

CAS Registry No.: 357399-43-0
Formal Name: 2-(3-chlorophenyl)-3,5,5-trimethyl-2-morpholinol
Synonyms: BUPOH, OH-bupropion
MF: C₁₃H₁₈ClNO₂
FW: 255.7
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

hydroxy Bupropion is supplied as a crystalline solid. A stock solution may be made by dissolving the hydroxy bupropion in the solvent of choice, which should be purged with an inert gas. hydroxy Bupropion is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of hydroxy Bupropion in these solvents is approximately 20 mg/ml.

hydroxy Bupropion is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hydroxy bupropion should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. hydroxy Bupropion has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

hydroxy Bupropion (Item No. 9000522) is an analytical reference standard categorized as an antidepressant.¹ hydroxy Bupropion is a metabolite of bupropion (Item No. 10488).² It substitutes for (+)-amphetamine (D-amphetamine; Item No. 26204) in rats in a two-lever drug discrimination test.³ This product is intended for research and forensic applications.

References

1. *Antidepressants: From biogenic amines to new mechanisms of action.* Springer, Switzerland (2018).
2. Coles, R. and Kharasch, E.D. Stereoselective metabolism of bupropion by cytochrome P4502B6 (CYP2B6) and human liver microsomes. *Pharm. Res.* **25(6)**, 1405-1411 (2008).
3. Bondarev, M.L., Bondareva, T.S., Young, R., et al. Behavioral and biochemical investigations of bupropion metabolites. *Eur. J. Pharmacol.* **474(1)**, 85-93 (2003).

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, 02/01/2022

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