

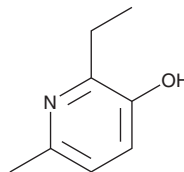
PRODUCT INFORMATION



Emoxypine

Item No. 45105

CAS Registry No.: 2364-75-2
Formal Name: 2-ethyl-6-methyl-3-pyridinol
Synonyms: Epigid, Epygid, Mexidol, Mexifin
MF: C₈H₁₁NO
FW: 137.2
Purity: ≥98%
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.

Description

Emoxypine (Item No. 45105) is an analytical reference standard categorized as a nootropic.¹ Emoxypine has been detected in dietary supplements marketed as nootropics.² This product is intended for research and forensic applications.

References

1. Volchegorskii, I.A., Miroshnichenko, I.Y., Rassokhina, L.M., *et al.* Anxiolytic and antidepressant effects of emoxipine, reamberin and mexidol in experimental diabetes mellitus. *Zh. Nevrol. Psikhiatr. Im. S.S. Korsakova* **117(5)**, 52-57 (2019).
2. Jędrejko, K., Catlin, O., Stewart, T., *et al.* Unauthorized ingredients in “nootropic” dietary supplements: A review of the history, pharmacology, prevalence, international regulations, and potential as doping agents. *Drug Test Anal.* **15(8)**, 803-839 (2023).

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, 03/20/2026

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