

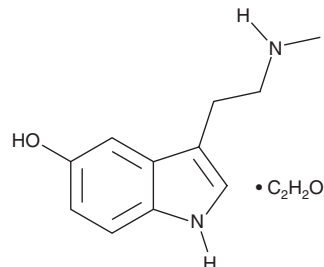
PRODUCT INFORMATION



5-hydroxy NMT (oxalate)

Item No. 26699

CAS Registry No.: 1975-81-1
Formal Name: 3-[2-(methylamino)ethyl]-1H-indol-5-ol, monoethanedioate
Synonyms: N-methyl-5-HT, N-methyl-5-Hydroxytryptamine, N-Methylserotonin, NSC 91541, 5-hydroxy-N-methyl Tryptamine, 5-hydroxy-N^ω-methyl Tryptamine
MF: C₁₁H₁₄N₂O • C₂H₂O₄
FW: 280.3
Purity: ≥95%
UV/Vis.: λ_{max}: 219, 277 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥3 years



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

Description

5-hydroxy NMT (oxalate) (Item No. 26699) is an analytical reference standard categorized as a tryptamine. 5-hydroxy NMT is a metabolite of serotonin (5-HT).^{1,2} Elevated plasma levels of 5-hydroxy NMT have been detected in individuals with cocaine dependence.¹ This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 20047).

References

1. Patkar, A.A., Rozen, S., Mannelli, P., *et al.* Alterations in tryptophan and purine metabolism in cocaine addiction: A metabolomic study. *Psychopharmacology (Berl)* **206(3)**, 479-489 (2009).
2. Heller, B. N-methylating enzyme in blood of schizophrenics. *Psychosomatics* **12(4)**, 273-274 (1971).

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, 07/03/2023

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