

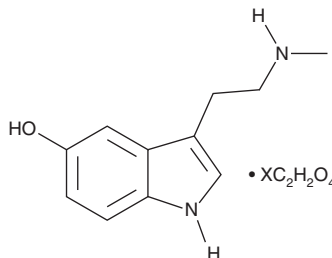
PRODUCT INFORMATION



5-hydroxy-N^ω-methyl Tryptamine (oxalate)

Item No. 20047

CAS Registry No.: 15558-50-6
Formal Name: 3-[2-(methylamino)ethyl]-1H-indol-5-ol, ethanedioate
Synonyms: N-methyl-5-HT, N-methyl Serotonin, N-methyl-5-hydroxy Tryptamine
MF: C₁₁H₁₄N₂O • XC₂H₂O₄
FW: 280.3
Purity: ≥95%
UV/Vis.: λ_{max}: 209, 276 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

5-hydroxy-N^ω-methyl Tryptamine (oxalate) is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-hydroxy-N^ω-methyl tryptamine (oxalate) in the solvent of choice, which should be purged with an inert gas. 5-hydroxy-N^ω-methyl Tryptamine (oxalate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 5-hydroxy-N^ω-methyl tryptamine (oxalate) in these solvents is approximately 3 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 5-hydroxy-N^ω-methyl tryptamine (oxalate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 5-hydroxy-N^ω-methyl tryptamine (oxalate) in PBS, pH 7.2, is approximately 2.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-hydroxy-N^ω-methyl Tryptamine is a metabolite of serotonin (5-HT; Item No. 14332) in humans that has also been found in plants.^{1,2,3} It is an agonist of the 5-HT receptor subtype 5-HT₇ (IC₅₀ = 23 pM in a radioligand binding assay).² 5-hydroxy-N^ω-methyl Tryptamine increases intracellular cyclic AMP production in HEK293 cells expressing 5-HT₇ (EC₅₀ = 22 nM). It inhibits serotonin uptake in HEK293 cells expressing human serotonin transporters (IC₅₀ = 490 nM). Elevated levels of urinary 5-hydroxy-N^ω-methyl tryptamine have been found in patients with schizophrenia, depression, and epilepsy.¹

References

1. Takeda, N., Ikeda, R., Ohba, K., *et al.* Bufotenine reconsidered as a diagnostic indicator of psychiatric disorders. *Neuroreport* **6**(17), 2378-2380 (1995).
2. Powell, S.L., Gödecke, T., Nikolic, D., *et al.* *In vitro* serotonergic activity of black cohosh and identification of N^ω-methylserotonin as a potential active constituent. *J. Agric. Food Chem.* **56**(24), 11718-11726 (2008).
3. 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, 10/04/2022

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