

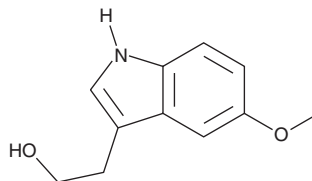
PRODUCT INFORMATION



5-methoxy Tryptophol

Item No. 21061

CAS Registry No.: 712-09-4
Formal Name: 5-methoxy-1H-indole-3-ethanol
Synonym: 5-MTOH
MF: C₁₁H₁₃NO₂
FW: 191.2
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 277 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

5-methoxy Tryptophol (5-MTOH) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 5-MTOH in these solvents is approximately 30 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. If an organic solvent-free solution of 5-MTOH is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 5-MTOH in PBS, pH 7.2, is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-MTOH is a natural indole that is produced by the pineal gland.^{1,2} It is a product of melatonin metabolism that may be biologically active.^{3,4} The levels of 5-MTOH in plasma vary in a diurnal pattern in rodents and humans.^{5,6}

References

1. Delvigs, P., Mclsaac, W.M., and Taborsky, R.G. The metabolism of 5-methoxytryptophol. *J. Biol. Chem.* **240**(1), 348-350 (1965).
2. Mullen, P.E., Leone, R.M., Hooper, J., et al. Pineal 5-methoxy tryptophol in man. *Psychoneuroendocrinology* **4**(2), 117-126 (1979).
3. Slominski, A., Semak, I., Pisarchik, A., et al. Conversion of L-tryptophan to serotonin and melatonin in human melanoma cells. *FEBS Lett.* **511**(1-3), 102-106 (2002).
4. Zisapel, N., Egozi, Y., and Laudon, M. Circadian variations in the inhibition of dopamine release from adult and newborn rat hypothalamus by melatonin. *Neuroendocrinology* **40**(2), 102-108 (1985).
5. Reddy, M.Y. and Jaqota, A. Melatonin has differential effects on age-induced stoichiometric changes in daily chronomics of serotonin metabolism in SCN of male Wistar rats. *Biogerontology* **16**(3), 285-302 (2015).
6. Wurtman, R.J. and Ozaki, Y. Physiological control of melatonin synthesis and secretion: Mechanisms, generating rhythms in melatonin, methoxytryptophol, and arginine vasotocin levels and effects on the pineal of endogenous catecholamines, the estrous cycle, and environmental lighting. *J. Neural Transm. Suppl.* **13**(13), 59-70 (1978).

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, 11/14/2022

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