

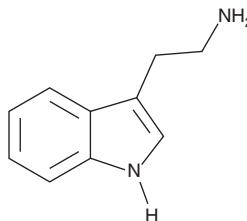
PRODUCT INFORMATION



Tryptamine

Item No. 20995

CAS Registry No.: 61-54-1
Formal Name: 1H-indole-3-ethanamine
Synonyms: 3-Indoleethylamine, NSC 165212
MF: C₁₀H₁₂N₂
FW: 160.2
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 282, 290 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

Tryptamine is supplied as a crystalline solid. A stock solution may be made by dissolving the tryptamine in the solvent of choice. Tryptamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of tryptamine in these solvents is approximately 10, 11, and 5 mg/ml, respectively.

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

Description

Tryptamine is an indole alkaloid and intermediate in the biosynthesis of serotonin and the phytohormone melatonin in plants.^{1,2} It increases the levels of the terpenoid indole alkaloids ajmalicine, strictosidine, and catharanthine in cultures of *C. roseus*.³ Tryptamine is also a product of tryptophan metabolism in mammals.⁴ Tryptamine derivatives have been synthetically produced as hallucinogenic drugs of abuse that act on the serotonergic system.⁵

References

1. Arnao, M.B. Phytomelatonin: Discovery, content, and role in plants. *Adv. Bot.* **815769**, (2014).
2. Fan, J., Xie, Y., Zhang, Z., *et al.* Melatonin: A multifunctional factor in plants. *Int. J. Mol. Sci.* **19(5)**, E1528 (2018).
3. Almagro, L., Fernández-Pérez, F., and Pedreño, M.A. Indole alkaloids from *Catharanthus roseus*: Bioproduction and their effect on human health. *Molecules* **20(2)**, 2973-3000 (2015).
4. Berumen, L.C., Rodríguez, A., Miledi, R., *et al.* Serotonin receptors in hippocampus. *ScientificWorldJournal* **823493**, (2012).
5. Araujo, A.M., Carvalho, F., Bastos, M.L., *et al.* The hallucinogenic world of tryptamines: An updated review. *Arch. Toxicol.* **89(8)**, 1151-1173 (2015).

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, 11/08/2022

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