

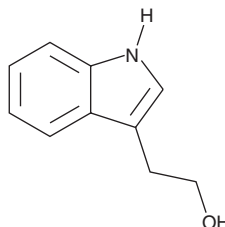
PRODUCT INFORMATION



Tryptophol

Item No. 28404

CAS Registry No.: 526-55-6
Formal Name: 1H-indole-3-ethanol
Synonyms: IEA, NSC 3884
MF: $C_{10}H_{11}NO$
FW: 161.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 223, 284 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Tryptophol is supplied as a crystalline solid. A stock solution may be made by dissolving the tryptophol in the solvent of choice, which should be purged with an inert gas. Tryptophol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of tryptophol in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Tryptophol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tryptophol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Tryptophol 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

Tryptophol is an indole that has been found in plants, bacteria, fungi, and sponges and has diverse biological activities, including antifungal, sleep-promoting, and quorum-sensing-related properties.¹⁻³ Tryptophol (100 μ M) inhibits the growth of the chytrid fungal skin pathogens *B. dendrobatidis* and *B. salamandrivorans* *in vitro*.¹ As a quorum-sensing molecule, it also induces autostimulation and increases tryptophol production in *B. dendrobatidis* and *B. salamandrivorans* cultures when used at a concentration of 1 μ M. Tryptophol (400 mg/kg) induces sleep in mice as assessed by loss of the righting reflex.³

References

1. Verbrugghe, E., Adriaensen, C., Martel, A., *et al.* Growth regulation in amphibian pathogenic chytrid fungi by the quorum sensing metabolite tryptophol. *Front. Microbiol.* **9**:3277 (2019).
2. Palmieri, A. and Petrini, M. Tryptophol and derivatives: Natural occurrence and applications to the synthesis of bioactive compounds. *Nat. Prod. Rep.* **36**(3), 490-530 (2019).
3. Feldstein, A., Chang, F.H., and Kucharski, J.M. Tryptophol, 5-hydroxytryptophol and 5-methoxytryptophol induced sleep in mice. *Life Sci.* **9**(6), 323-329 (1970).

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, 12/01/2022

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