

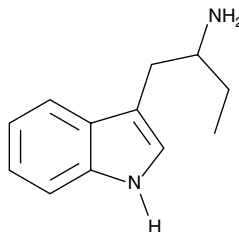
Product Information



α -Ethyltryptamine (exempt preparation)

Item No. 15671

CAS Registry No.: 2235-90-7
Formal Name: α -ethyl-1H-indole-3-ethanamine
Synonyms: α -ET, 3-Indolybutylamine,
NSC 88061, Ro 3-1932
MF: C₁₂H₁₆N₂
FW: 188.3
Purity: \geq 95%
Stability: \geq 2 years at -20°C
Supplied as: A solution in methanol
UV/Vis.: λ_{\max} : 222, 282 nm



Laboratory Procedures

For long term storage, we suggest that α -ethyltryptamine (α -ET) (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

α -ET (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol 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 α -ET (exempt preparation) in these solvents is approximately 20, 5, and 11 mg/ml, respectively.

α -ET (exempt preparation) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methanolic solution of α -ET (exempt preparation) should be diluted with the aqueous buffer of choice. α -ET (exempt preparation) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

α -ET is a tryptamine derivative with schedule I status in the United States.^{1,2} It has been reported to be a potent releaser of serotonin, norepinephrine, and dopamine as well as an inhibitor of monoamine oxidase.^{3,4} In tests of stimulus generalization in rats, α -ET generalized to DOM (Item No. 11145) and PMMA (Item No. 11562) with ED₅₀ values of 0.4 and 0.7 mg/kg, respectively.³ This product is intended for forensic and research applications.

References

1. Gey, K.F. and Pletscher, A. Effect of α -alkylated tryptamine derivatives on 5-hydroxytryptamine metabolism *in vivo*. *Brit. J. Pharmacol.* **19**, 161-167 (1962).
2. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).
3. Glennon, R.A., Bondareva, T., and Young, R. α -Ethyltryptamine (α -ET) as a discriminative stimulus in rats. *Pharmacol. Biochem. Behav.* **85(2)**, 448-453 (2006).
4. Martinez, D.L. and Geyer, M.A. Characterization of the disruptions of prepulse inhibition and habituation of startle induced by α -ethyltryptamine. *Neuropsychopharmacology* **16(3)**, 246-255 (1997).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/15671

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/31/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com