

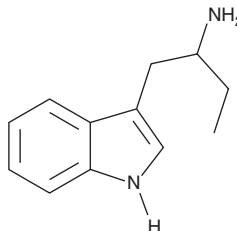
PRODUCT INFORMATION



α -Ethyltryptamine

Item No. 14653

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: $C_{12}H_{16}N_2$
FW: 188.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 222, 282 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 5 years



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

Description

α -Ethyltryptamine (α -ET) is a tryptamine derivative with schedule 1 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 ED_{50} 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. *Br. J. Pharmacol. Chemother.* **19(1)**, 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).

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, 06/21/2023

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