

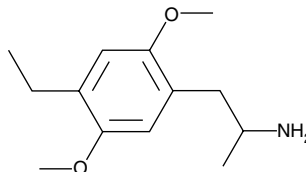
Product Information



DOET (exempt preparation)

Item No. 15704

CAS Registry No.: 22004-32-6
Formal Name: 4-ethyl-2,5-dimethoxy- α -methylbenzeneethanamine
MF: C₁₃H₂₁NO₂
FW: 223.3
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A solution in methanol
UV/Vis.: λ_{\max} : 225, 291 nm



Laboratory Procedures

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

DOET (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 and dimethyl formamide purged with an inert gas can be used. The solubility of DOET (exempt preparation) in ethanol is approximately 0.3 and 0.5 mg/ml, respectively.

DOET (exempt preparation) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

DOET is a potent, long-acting psychoactive drug of the phenethylamine and amphetamine chemical classes that, similar to other substituted amphetamines, likely acts as a serotonin 5-HT₂ receptor partial agonist.^{1,2} This compound is classified as a Schedule I substance in the United States, and is intended only for research and forensic applications.

References

1. Ewald, A.H. and Maurer, H.H. 2,5-dimethoxyamphetamine-derived designer drugs: Studies on the identification of cytochrome P450 (CYP) isoenzymes involved in formation of their main metabolites and on their capability to inhibit CYP2D6. *Toxicol. Lett.* **183**, 52-57 (2008).
2. Snyder, S.H., Weingartner, H., and Faillace, L.A. DOET (2,5-dimethoxy-4-ethylamphetamine), a new psychotropic drug. Effects of varying doses in man. *Arch. Gen. Psychiatry* **24(1)**, 50-55 (1971).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/15704

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, 04/07/2014

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