

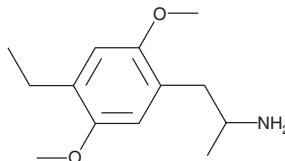
PRODUCT INFORMATION



DOET (exempt preparation)

Item No. 15704

CAS Registry No.: 22004-32-6
Formal Name: 4-ethyl-2,5-dimethoxy- α -methyl-
benzeneethanamine
MF: $C_{13}H_{21}NO_2$
FW: 223.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 225, 291 nm
Supplied as: A solution in methanol
Storage: $-20^{\circ}C$
Stability: ≥ 3 years



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

Description

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).

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, 07/02/2025

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