

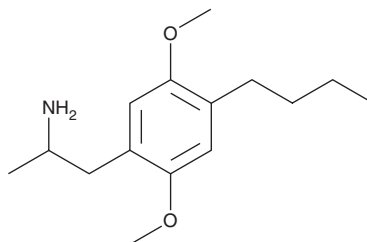
PRODUCT INFORMATION



DOBU

Item No. 17655

CAS Registry No.: 63779-89-5
Formal Name: 4-butyl-2,5-dimethoxy- α -methylbenzeneethanamine
Synonym: 2,5-Dimethoxy-4-butylamphetamine
MF: C₁₅H₂₅NO₂
FW: 251.4
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 225, 291 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

DOBU (Item No. 17655) is an analytical reference standard that is structurally categorized as an amphetamine. Experiments with isolated rat thoracic aorta suggested that this compound may act as an antagonist of the 5-HT₂ receptor.¹ This product is intended for research and forensic applications.

Reference

1. Seggel, M.R., Yousif, M.Y., Lyon, R.A., *et al.* A structure-affinity study of the binding of 4-substituted analogues of 1-(2,5-dimethoxyphenyl)-2-aminopropane at 5-HT₂ serotonin receptors. *J. Med. Chem.* **33**(3), 1032-1036 (1990).

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, 06/16/2023

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