

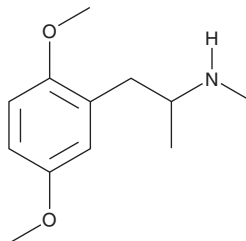
PRODUCT INFORMATION



2,5-DMMA

Item No. 15131

CAS Registry No.: 54687-43-3
Formal Name: 2,5-dimethoxy-N, α -dimethylbenzeneethanamine
Synonym: 2,5-Dimethoxymethamphetamine
MF: C₁₂H₁₉NO₂
FW: 209.3
Purity: $\geq$ 98%
UV/Vis.: λ_{max} : 291 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: $\geq$ 4 years



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

Description

2,5-DMMA (Item No. 15131) is a designer drug that combines the 2,5 disubstitution pattern of 2,5-dimethoxyphenethylamines with the backbone of methamphetamine. While the physiological and toxicological properties of this compound are poorly studied, the absence of an additional third substitution at either the three or four position on similar phenethylamines and amphetamines typically reduces effects on monoamine action.^{1,2} This product is intended for forensic and research applications.

References

1. Blaazer, A.R., Smid, P., and Kruse, C.G. Structure-activity relationships of phenylalkylamines as agonist ligands for 5-HT_{2A} receptors. *ChemMedChem* **3**(9), 1299-1309 (2008).
2. Moya, P.R., Berg, K.A., Gutiérrez-Hernandez, M.A., et al. Functional selectivity of hallucinogenic phenethylamine and phenylisopropylamine derivatives at human 5-hydroxytryptamine (5-HT)_{2A} and 5-HT_{2C} receptors. *J. Pharmacol. Exp. Ther.* **321**(3), 1054-1061 (2007).

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, 01/14/2026

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