

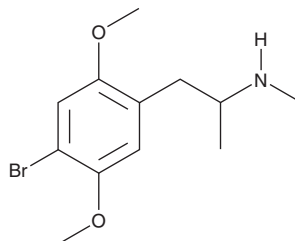
PRODUCT INFORMATION



4-bromo-2,5-DMMA

Item No. 14788

CAS Registry No.: 155638-80-5
Formal Name: 4-bromo-2,5-dimethoxy-N,α-dimethylbenzeneethanamine
Synonym: 4-bromo-2,5-Dimethoxymethamphetamine
MF: C₁₂H₁₈BrNO₂
FW: 288.2
Purity: ≥98%
UV/Vis.: λ_{max}: 296 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

4-bromo-2,5-DMMA (Item No. 14788) is a designer drug patterned after 2,5-dimethylamphetamine compounds on the illicit drug market.^{1,2} The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research purposes.

References

1. Maurer, H.H. Chemistry, pharmacology, and metabolism of emerging drugs of abuse. *Ther. Drug Monit.* **32**(5), 544-549 (2010).
2. 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**(1-3), 52-57 (2008).

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, 01/13/2026

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