

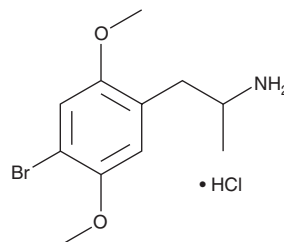
PRODUCT INFORMATION



4-bromo-2,5-DMA (hydrochloride)

Item No. 11142

CAS Registry No.: 29705-96-2
Formal Name: 4-bromo-2,5-dimethoxy- α -methylbenzeneethanamine, monohydrochloride
Synonyms: 4-bromo-2,5-Dimethoxyamphetamine, Dimethoxybromoamphetamine, DOB
MF: C₁₁H₁₆BrNO₂ • HCl
FW: 310.6
Purity: ≥98%
UV/Vis.: λ_{max} : 206, 297 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

4-bromo-2,5-DMA is a common hallucinogenic designer drug which avidly binds and activates serotonin receptors ($K_i = 2.2$ and 2.8 nM for 5-HT_{2A} and 5-HT_{2C}, respectively).^{1,2} Methods have been developed for the identification of this compound in serum and in urine.^{3,4} 4-bromo-2,5-DMA is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. de Boer, D. and Bosman, I. A new trend in drugs-of-abuse; the 2C-series of phenethylamine designer drugs. *Pharm. World Sci.* **26(2)**, 110-113 (2004).
2. Chambers, J.J., Kurrasch-Orbaugh, D.M., Parker, M.A., et al. Enantiospecific synthesis and pharmacological evaluation of a series of super-potent, conformationally restricted 5-HT_{2A/2C} receptor agonists. *J. Med. Chem.* **44(6)**, 1003-1010 (2001).
3. Wohlfarth, A., Weinmann, W., and Dresen, S. LC-MS/MS screening method for designer amphetamines, tryptamines, and piperazines in serum. *Anal. Bioanal. Chem.* **396(7)**, 2403-2414 (2010).
4. Kerrigan, S., Banuelos, S., Perrella, L., et al. Simultaneous detection of ten psychedelic phenethylamines in urine by gas chromatography-mass spectrometry. *J. Anal. Toxicol.* **35(7)**, 459-469 (2011).

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, 04/13/2022

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