

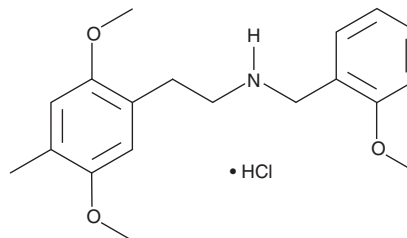
PRODUCT INFORMATION



25D-NBOMe (hydrochloride)

Item No. 11946

CAS Registry No.: 1539266-35-7
Formal Name: 2,5-dimethoxy-N-[(2-methoxyphenyl)methyl]-4-methyl-benzeneethanamine, monohydrochloride
Synonym: 2C-D-NBOMe
MF: C₁₉H₂₅NO₃ • HCl
FW: 351.9
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 220, 282 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

2C-D (Item No. 11888), a 2,5-dimethoxy phenethylamine, is designer drug with mild hallucinogenic activity, due in part to weak activity at serotonin receptors.^{1,2} 25D-NBOMe is a derivative of 2C-D, characterized by the addition of a benzyl-methoxy (BOMe) group to the amine. While the properties of this compound are not known, the N-(2-methoxybenzyl) addition significantly increases the affinity and activity of other disubstituted phenethylamines at the serotonin receptor 5-HT_{2A}.^{3,4} This product is intended for research and forensic applications.

References

1. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11**(5), 468-482 (2010).
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).
3. Braden, M.R., Parrish, J.C., Naylor, J.C., *et al.* Molecular interaction of serotonin 5-HT_{2A} receptor residues Phe339^{6.51} and Phe340^{6.52} with superpotent N-benzyl phenethylamine agonists. *Mol. Pharm.* **70**(6), 1956-1965 (2006).
4. Ettrup, A., Palner, M., Gillings, N., *et al.* Radiosynthesis and evaluation of ¹¹C-CIMBI-5 as a 5-HT_{2A} receptor agonist radioligand for PET. *J.Nuclear Med.* **51**(11), 1763-1770 (2010).

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, 07/24/2023

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