

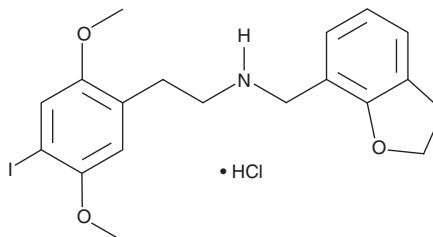
PRODUCT INFORMATION



25I-NBMD (hydrochloride)

Item No. 11947

CAS Registry No.: 1539266-14-2
Formal Name: N-(benzo[d][1,3]dioxol-4-ylmethyl)-2-(4-iodo-2,5-dimethoxyphenyl)ethanamine, monohydrochloride
Synonym: Cimbi-29
MF: C₁₈H₂₀INO₄ • HCl
FW: 477.7
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 235, 294 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

25I-NBMD is a derivative of the phenethylamine hallucinogen 2C-I, characterized by the addition of a methylenedioxy-benzyl group to the amine. This addition increases the affinity ($K_i = 0.049\text{--}0.21\text{ nM}$) and activity ($EC_{50} = 8.2\text{ nM}$) at the serotonin receptor 5-HT_{2A}.^{1,2} This product is intended for research and forensic applications.

References

1. Ettrup, A., Hansen, M., Santini, M.A., *et al.* Radiosynthesis and *in vivo* evaluation of a series of substituted ¹¹C-phenethylamines as 5-HT_{2A} agonist PET tracers. *Eur. J. Nucl. Med. Mol. Imaging* **38(4)**, 681-693 (2011).
2. 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).

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