

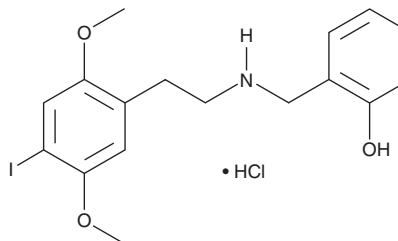
PRODUCT INFORMATION



25I-NBOH (hydrochloride)

Item No. 14909

CAS Registry No.: 1539266-12-0
Formal Name: 2-(((4-iodo-2,5-dimethoxyphenethyl)amino)methyl)phenol, monohydrochloride
Synonym: 2C-I-NBOH
MF: $C_{17}H_{20}INO_3 \cdot HCl$
FW: 449.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 299 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

25I-NBOH is a derivative of the phenethylamine hallucinogen 2C-I (Item No. 11113) with an N-(2-hydroxybenzyl) addition at the amine. It can act as a potent agonist for the human serotonin 5-HT_{2A} receptor ($K_i = 0.061$ nM) with 12-fold greater potency compared to 2C-I itself.¹⁻² This product is intended for research and forensic applications.

References

1. 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).
2. 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).

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, 06/22/2023

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