

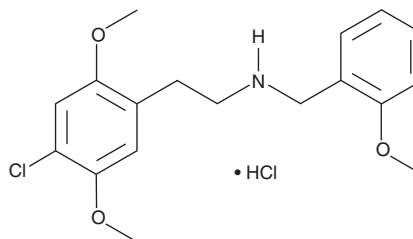
PRODUCT INFORMATION



25C-NBOMe (hydrochloride)

Item No. 9001096

CAS Registry No.: 1539266-19-7
Formal Name: 2-(4-chloro-2,5-dimethoxyphenyl)-
N-(2-methoxybenzyl)ethanamine,
monohydrochloride
Synonym: 2C-C-NBOMe
MF: C₁₈H₂₂ClNO₃ • HCl
FW: 372.3
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 295 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-NBOMe (Item No. 9001096) is an analytical reference standard that is structurally categorized as a phenethylamine. 25C-NBOMe is a derivative of 2C-C having an N-(2-methoxybenzyl) addition at the amine. 2C-C is a known hallucinogen that stimulates monoamine receptor activity and inhibits the re-uptake of serotonin and norepinephrine in rat brain synaptosomes (IC₅₀ = 31 and 63 μM, respectively).^{1,2} The N-(2-methoxybenzyl) addition to 2C-NBOMe increases the affinity and selectivity for the 5-HT_{2A} receptor over other serotonin receptors.³ Radiolabeled 25C-NBOMe has been used for positron emission tomography imaging of the 5-HT_{2A} receptor in porcine brain.⁴ This product is intended for forensic and research uses.

References

1. Nonaka, R., Nagai, F., Ogata, A., *et al.* In vitro screening of psychoactive drugs by [³⁵S]GTPγS binding in rat brain membranes. *Biol. Pharm. Bull.* **30(12)**, 2328-33 (2007).
2. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (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., 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, 06/23/2023

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