

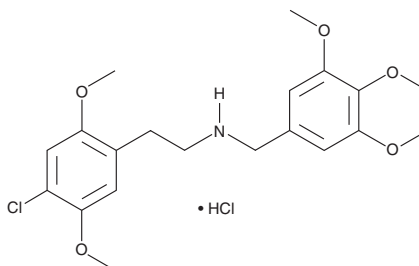
PRODUCT INFORMATION



30C-NBOMe (hydrochloride)

Item No. 14528

CAS No.: 2749391-65-7
Formal Name: 2-(4-chloro-2,5-dimethoxyphenyl)-
N-(3,4,5-trimethoxybenzyl)
ethanamine, monohydrochloride
MF: C₂₀H₂₆ClNO₅ • HCl
FW: 432.3
Purity: ≥95%
UV/Vis.: λ_{max}: 294, 351 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

25C-NBOMe (hydrochloride) (Item No. 9001096) is a derivate of the 2,5-dimethoxy phenethylamine hallucinogen 2C-C (Item No. 11735). The N-(2-methoxybenzyl) addition at the amine of 2C-C, to give 25C-NBOMe, increases affinity and selectivity for the 5-HT_{2A} receptor over other serotonin receptors.¹ 30C-NBOMe is an analog of 25C-NBOMe having methoxy groups at the three, four, and five positions, rather than the two position, on the benzene ring. While the 3,4,5-trimethoxybenzyl structure is like that found in the psychedelic alkaloid mescaline (Item No. 11950), this compound has a shortened chain length. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research purposes.

Reference

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).

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, 06/20/2023

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