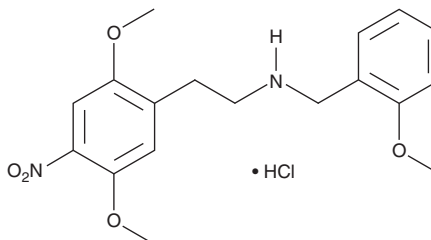


PRODUCT INFORMATION

25N-NBOMe (hydrochloride)

Item No. 14911

CAS Registry No.: 1566571-65-0
Formal Name: 2,5-dimethoxy-N-[(2-methoxyphenyl)methyl]-4-nitro-benzeneethanamine, monohydrochloride
Synonym: 2C-N-NBOMe
MF: C₁₈H₂₂N₂O₅ • HCl
FW: 382.8
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 274, 348, 356 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-N (Item No. 11890) is described formally as 2,5-dimethoxy-4-nitrophenethylamine and is structurally analogous to the psychedelic drug 2C-B (Item No. 11734). 2C-N potently stimulates arachidonic acid (Item No. 90010) release through the serotonin receptors 5-HT_{2C} (pEC₅₀ = 5.91) and 5-HT_{2A} (pEC₅₀ = 4.78).¹ 25N-NBOMe is a derivative of 2C-N having an N-(2-methoxybenzyl) addition at the amine. For similar dimethoxyphenethylamines including 2C-B, this N-(2-methoxybenzyl) addition significantly increases activity at the serotonin receptor.² This product is intended for forensic and research applications.

References

1. 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).
2. 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. Nucl. 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, 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