

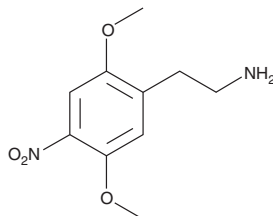
PRODUCT INFORMATION



2C-N (exempt preparation)

Item No. 15717

CAS Registry No.: 261789-00-8
Formal Name: 2,5-dimethoxy-4-nitrobenzeneethanamine
Synonym: 2,5-Dimethoxy-4-nitrophenethylamine
MF: C₁₀H₁₄N₂O₄
FW: 226.2
Purity: ≥98%
Storage: -20°C
Stability: ≥3 years
Supplied as: A solution in methanol
UV/Vis.: λ_{max}: 220, 275, 361 nm



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

Description

A series of 2,5-phenethylamines, collectively referred to as 2Cs, have psychoactive effects.^{1,2} The most effective 2C compounds are 2,5-dimethoxy-4-substituted phenethylamines; many are regulated as illegal substances.^{3,4} 2C-N is described formally as 2,5-dimethoxy-4-nitrophenethylamine and is structurally analogous to the psychedelic drug 2C-B. 2C-N potentially stimulates arachidonic acid release through the serotonin receptors 5-HT_{2C} (pEC₅₀ = 5.91) and 5-HT_{2A} (pEC₅₀ = 4.78).² This product is intended for forensic and research applications.

References

1. Bruno, R., Matthews, A.J., Dunn, M., *et al.* Emerging psychoactive substance use among regular ecstasy users in Australia. *Drug Alcohol Depend.* **124**(1-2), 19-25 (2012).
2. 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**, 1054-1061 (2007).
3. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11**, 468-482 (2010).
4. 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).

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, 04/08/2026

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