

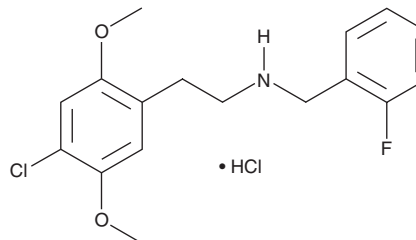
PRODUCT INFORMATION



25C-NBF (hydrochloride)

Item No. 15969

CAS Registry No.: 1539266-21-1
Formal Name: 4-chloro-N-[(2-fluorophenyl)methyl]-2,5-dimethoxy-benzeneethanamine, monohydrochloride
Synonym: 2C-C-NBF
MF: C₁₇H₁₉ClFNO₂ • HCl
FW: 360.3
Purity: ≥98%
UV/Vis.: λ_{max}: 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-C (Item No. 11735) is a 2,5-dimethoxyphenethylamine with a chlorine atom at the four position of the aromatic ring. A known hallucinogen, this compound stimulates monoamine receptor activity while inhibiting the re-uptake of serotonin (Item No. 14332) and norepinephrine.^{1,2} 25C-NBF (hydrochloride) is a derivative of 2C-C having an N-(2-fluorobenzyl) addition at the amine. It is an agonist of serotonin receptors that activates 5-HT_{2A} and 5-HT_{2C} similarly (EC₅₀s = ~0.3 μM for each).³ This product is intended for forensic and research applications.

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. Hansen, M., Phonekeo, K., Paine, J.S., *et al.* Synthesis and structure-activity relationships of N-benzyl phenethylamines as 5-HT_{2A/2C} agonists. *ACS Chem. Neurosci.* **5**(3), 243-249 (2014).

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, 06/29/2023

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