

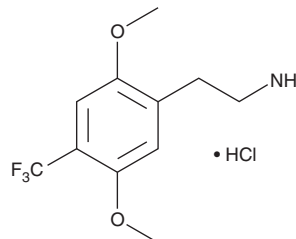
PRODUCT INFORMATION



2C-TFM (hydrochloride)

Item No. 14471

CAS Registry No.: 159277-13-1
Formal Name: 2,5-dimethoxy-4-(trifluoromethyl)-benzeneethanamine, monohydrochloride
Synonym: 2C-CF3
MF: C₁₁H₁₄F₃NO₂ • HCl
FW: 285.7
Purity: ≥95%
UV/Vis.: λ_{max}: 230, 299 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

A series of 2,5-dimethoxyphenethylamines, collectively referred to as 2Cs, have psychoactive effects.^{1,2} The most effective 2C compounds are substituted at the 4 position of the aromatic ring; many are scheduled as illegal substances.^{3,4} 2C-TFM is a 2C compound with a *para*-trifluoromethyl substitution. It is a potent agonist of the serotonin (5-HT) receptor subtypes 5-HT_{2A} and 5-HT_{2C} (K_{aff} = 165 and 32 nM).⁵ This product is intended for forensic and research purposes.

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(3)**, 1054-1061 (2007).
3. Meyer, M.R. and Maurer, H.H. Metabolism of designer drugs of abuse: An updated review. *Curr. Drug Metab.* **11(5)**, 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).
5. Nichols, D.E., Frescas, S., Marona-Lewicka, D., *et al.* 1-(2,5-Dimethoxy-4-(trifluoromethyl)phenyl)-2-aminopropane: A potent serotonin 5-HT_{2A/2C} agonist. *J. Med. Chem.* **37(25)**, 4346-4351 (1994).

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, 06/19/2023

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