

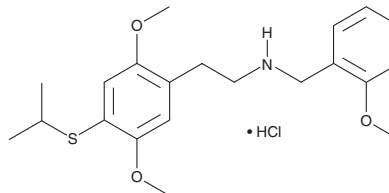
PRODUCT INFORMATION



25T4-NBOMe (hydrochloride)

Item No. 9001989

CAS Registry No.: 1566571-73-0
Formal Name: 2,5-dimethoxy-N-[(2-methoxyphenyl)methyl]-4-[(1-methylethyl)thio]benzeneethanamine, monohydrochloride
MF: C₂₁H₂₉NO₃S • HCl
FW: 412.0
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 306 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-T-4 (Item No. 11892) is described formally as 2,5-dimethoxy-4-(isopropylthio)phenethylamine. The physiological and toxicological effects of this designer drug have not been investigated, although there is a case report of 2C-T-4 causing acute psychosis.¹ 25T4-NBOMe is a derivative of 2C-T-4, characterized by the addition of a benzyl-methoxy (BOMe) group to the amine. While the properties of this compound are not known, the N-(2-methoxybenzyl) addition significantly increases the affinity and activity of other disubstituted phenethylamines at the serotonin receptor 5-HT_{2A}.^{2,3} This product is intended for forensic and research applications.

References

1. Miyajima, M., Matsumoto, T., and Ito, S. 2C-T-4 intoxication: Acute psychosis caused by a designer drug. *Psychiatry Clin. Neurosci.* **62**(2), 243 (2008).
2. 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).
3. 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. Nuclear 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, 07/14/2023

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