

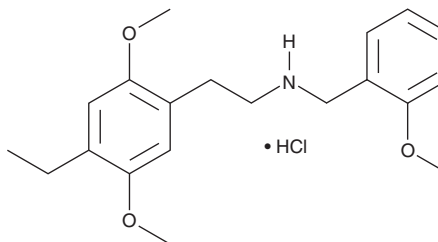
PRODUCT INFORMATION



25E-NBOMe (hydrochloride)

Item No. 14727

CAS Registry No.: 1539266-39-1
Formal Name: 4-ethyl-2,5-dimethoxy-N-
[(2-methoxyphenyl)methyl]-
benzeneethanamine, monohydrochloride
Synonym: 2C-E-NBOMe
MF: C₂₀H₂₇NO₃ • HCl
FW: 365.9
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 282 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-E, known formally as 2,5-dimethoxy-4-ethyl phenethylamine, is a psychoactive compound that is regulated as Schedule I in the United States.^{1,2} 25E-NBOMe is a derivative of 2C-E having an N-(2-methoxybenzyl) addition at the amine. This addition on related phenethylamines increases the affinity and selectivity for the 5-HT_{2A} receptor over other serotonin receptors.³ This product is intended for forensic and research purposes.

References

1. Blaazer, A.R., Smid, P., and Kruse, C.G. Structure-activity relationships of phenylalkylamines as agonist ligands for 5-HT_{2A} receptors. *ChemMedChem* **3**(9), 1299-1309 (2008).
2. Maurer, H.H. Chemistry, pharmacology, and metabolism of emerging drugs of abuse. *Ther. Drug Monit.* **32**(5), 544-549 (2010).
3. 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).

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, 06/21/2023

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