

PRODUCT INFORMATION



25H-NBOMe (hydrochloride)

Item No. 12061

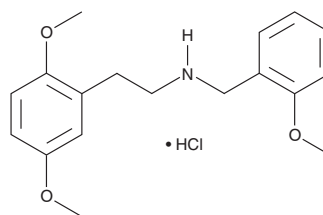
CAS Registry No.: 1566571-52-5
Formal Name: 2-(2,5-dimethoxyphenyl)-N-(2-methoxybenzyl)ethanamine, monohydrochloride

Synonyms: 2C-H-NBOMe
MF: C₁₈H₂₃NO₃ • HCl
FW: 337.8

Purity: ≥98%
Supplied as: A neat solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Description

25H-NBOMe, an N-benzyl analog of the phenethylamines mescaline and 2C-B, is an agonist for the rat and human serotonin 5-HT_{2A} receptors with K_i values of 1.19 and 2.83 nM, respectively.¹ In an *in vitro* activity assay, 25H-NBOMe demonstrates EC₅₀ values of 81.2 and 15.3 nM for the rat and human serotonin 5-HT_{2A} receptors, respectively.¹ This product is intended for research and forensic applications.

This product is a qualified Reference Material (RM) that has been manufactured and tested to meet ISO17025 and Guide 34 guidelines. These materials are tested using validated analytical methods on qualified instrumentation to ensure traceability of measurements. All traceable RMs may be distinguished by their CofAs and can be downloaded below using the batch number located on the product label. For a representative CofA please contact our technical support.

Reference

1. 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, 06/07/2016

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