

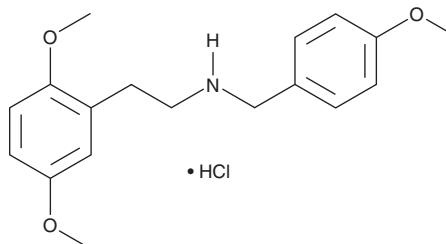
PRODUCT INFORMATION



25H-NB4OMe (hydrochloride)

Item No. 19365

CAS Registry No.: 1566571-54-7
Formal Name: 2,5-dimethoxy-N-[(4-methoxyphenyl)methyl]-benzeneethanamine, monohydrochloride
Synonym: 25H-NBOMe 4-methoxy isomer
MF: $C_{18}H_{23}NO_3 \cdot HCl$
FW: 337.8
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 282, 291 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

25H-NB4OMe (hydrochloride) (Item No. 19365) is an analytical reference standard that is structurally categorized as a phenethylamine. The biological properties of this compound have not been evaluated, although mass spectra characterization has been reported.¹ This product is intended for research and forensic applications.

Reference

1. Casale, J.F. and Hays, P.A. Characterization of eleven 2,5-dimethoxy-N-(2-methoxybenzyl)phenethylamine (NBOMe) derivatives and differentiation from their 3- and 4-methoxybenzyl analogues - Part I. *Microgram Journal* **9**(2), 84-109 (2012).

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