

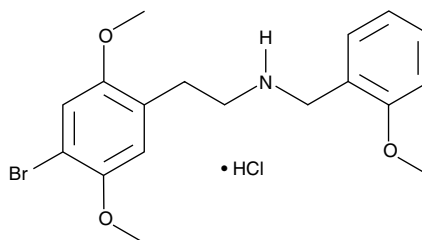
Product Information



25B-NBOMe (hydrochloride) (exempt preparation)

Item No. 15723

CAS Registry No.: 1539266-15-3
Formal Name: 4-bromo-2,5-dimethoxy-N-[(2-methoxyphenyl)methyl]-benzeneethanamine, monohydrochloride
Synonym: 2C-B-NBOMe
MF: C₁₈H₂₂BrNO₃ • HCl
FW: 416.7
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A solution in methanol
UV/Vis.: λ_{max}: 204, 282, 296 nm



Laboratory Procedures

For long term storage, we suggest that 25B-NBOMe (hydrochloride) (exempt preparation) be stored as supplied at -20°C. It should be stable for at least two years.

25B-NBOMe (hydrochloride) (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 25B-NBOMe (hydrochloride) (exempt preparation) in these solvents is approximately 20 mg/ml.

25B-NBOMe (hydrochloride) (exempt preparation) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 25B-NBOMe (hydrochloride) (exempt preparation) should be diluted with the aqueous buffer of choice. 25B-NBOMe (hydrochloride) (exempt preparation) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

2C-B is a hallucinogenic designer drug which potently activates the serotonin receptor 5-HT_{2C} (pEC₅₀ = 6.8 for arachidonic acid release).¹ 25B-NBOMe is a derivative of 2C-B, characterized by the addition of a benzyl-methoxy (BOMe) group to the amine. This N-(2-methoxybenzyl) addition significantly increases the affinity (K_i = 1.01 nM) and activity (ED₅₀ = 0.51 nM) at the serotonin receptor 5-HT_{2A}.² This product is intended for research and forensic applications.

References

1. 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**, 1054-1061 (2007).
2. Etrup, A., Hansen, M., Santini, M.A., *et al.* Radiosynthesis and *in vivo* evaluation of a series of substituted ¹¹C-phenethylamines as 5-HT_{2A} agonist PET tracers. *Eur. J. Nucl. Med. Mol. Imaging* **38**, 681-693 (2011).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/15723

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 04/09/2014

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com