

PRODUCT INFORMATION



25B-NBF (hydrochloride)

Item No. 15967

CAS Registry No.: 1539266-17-5

Formal Name: 4-bromo-N-[(2-fluorophenyl)methyl]-
2,5-dimethoxy-benzeneethanamine,
monohydrochloride

Synonym: 2C-B-NBF

MF: $C_{17}H_{19}BrFNO_2 \cdot HCl$

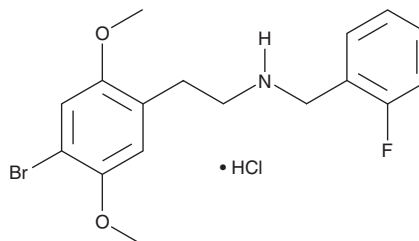
FW: 404.7

Purity: $\geq 98\%$

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

25B-NBF is an *N*-fluorobenzyl derivative of the phenethylamine hallucinogen 2C-B (Item No. 11734), which acts as a highly potent agonist for the 5-HT_{2C} receptor. 25B-NBF binds human 5-HT_{2A} receptors and rat 5-HT_{2C} receptors with pK_i values of 8.57 and 7.73, respectively.¹ This product is intended for forensic and research applications.

References

1. Hansen, M., Phonekeo, K., Paine, J.S., *et al.* Synthesis and structure-activity relationships of *N*-benzyl phenethylamines as 5-HT_{2A/2C} agonists. *ACS Chem. Neurosci.* **5**(3), 243-249 (2014).

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, 06/29/2023

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