

PRODUCT INFORMATION



25I-NBF (hydrochloride)

Item No. 14066

CAS Registry No.: 1539266-13-1

Formal Name: N-(2-fluorobenzyl)-2-(4-iodo-2,5-dimethoxyphenyl)ethanamine, monohydrochloride

Synonym: 2C-I-NBF

MF: $C_{17}H_{19}FINO_2 \cdot HCl$

FW: 451.7

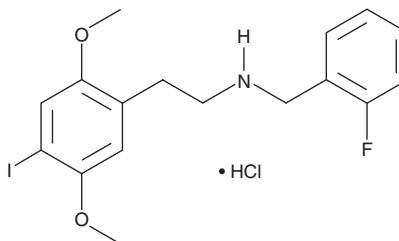
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 234, 300 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

25I-NBF is an N-benzyl derivative of the phenethylamine hallucinogen 2C-I which acts as a highly potent partial agonist for the 5-HT_{2A} receptor. 25I-NBF acts as a partial agonist of human 5-HT_{2A} receptors demonstrating a K_i value of 0.26 nM and an EC₅₀ value of 1.6 nM.¹ A radiolabelled form of 25I-NBF has been used for mapping the distribution of 5-HT_{2A} receptors in the brain, using positron emission tomography.² This product is intended for forensic and research applications.

References

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).
2. Ettrup, 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).

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, 07/25/2023

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