

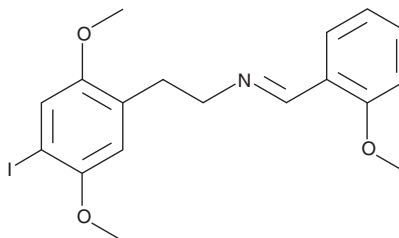
PRODUCT INFORMATION



25I-NBOMe imine analog

Item No. 11968

CAS Registry No.: 2177287-33-9
Formal Name: (E)-2-(4-iodo-2,5-dimethoxyphenyl)-N-(2-methoxybenzylidene)ethanamine
MF: C₁₈H₂₀INO₃
FW: 425.3
Purity: ≥95%
UV/Vis.: λ_{max}: 206, 236, 305 nm
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥5 years



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

Description

25I-NBOMe, a derivative of the serotonin 2A (5-HT_{2A}) receptor agonist 2C-I, is 16-fold more potent than 2C-I with a K_i value of 0.044 nM for human 5-HT_{2A} receptors.^{1,2} 25I-NBOMe imine analog differs from 25I-NBOMe (Item No. 9001128) by having a double bond linking the N-(2-methoxybenzyl) group to 25I. The properties of this compound have not been evaluated. This product is intended for research and forensic 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., Palner, M., Gillings, N., *et al.* Radiosynthesis and evaluation of ¹¹C-CIMBI-5 as a 5-HT_{2A} receptor agonist radioligand for PET. *J. Nucl. Med.* **51**(11), 1763-1770 (2010).

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, 11/21/2025

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