

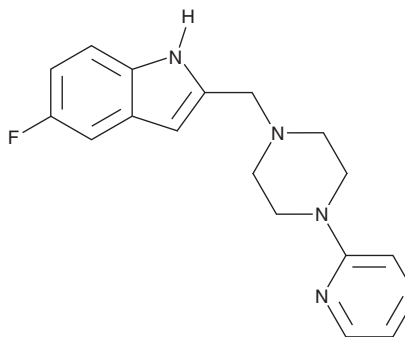
PRODUCT INFORMATION



CP 226,269

Item No. 36493

CAS Registry No.: 220941-93-5
Formal Name: 5-fluoro-2-[[4-(2-pyridinyl)-1-piperazinyl]methyl]-1H-indole
MF: C₁₈H₁₉FN₄
FW: 310.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

CP 226,269 is supplied as a solid. A stock solution may be made by dissolving the CP 226,269 in the solvent of choice, which should be purged with an inert gas. CP 226,269 is soluble in the organic solvent DMSO at a concentration of approximately 31 mg/ml.

Description

CP 226,269 is a dopamine D₄ receptor agonist.¹ It induces calcium flux in HEK293 cells co-expressing human dopamine D₄, but not D_{2L} or D₃, receptors and G_{αqo5} (EC₅₀s = 32, >10,000, and 985 nM, respectively). CP 226,269 induces phospholipid methylation in SK-N-MC neuroblastoma cells, which endogenously express dopamine D₁ and D₄ receptors (EC₅₀ = 2.9 nM).²

References

1. Moreland, R.B., Nakane, M., Donnelly-Roberts, D., *et al.* Comparative pharmacology of human dopamine D₂-like receptor stable cell lines coupled to calcium flux through G_{αqo5}. *Biochem. Pharmacol.* **68**(4), 761-772 (2004).
2. Sharma, A., Kramer, M.L., Wick, P.F., *et al.* D4 dopamine receptor-mediated phospholipid methylation and its implications for mental illnesses such as schizophrenia. *Mol. Psychiatry* **4**(3), 235-246 (1999).

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, 12/19/2022

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