

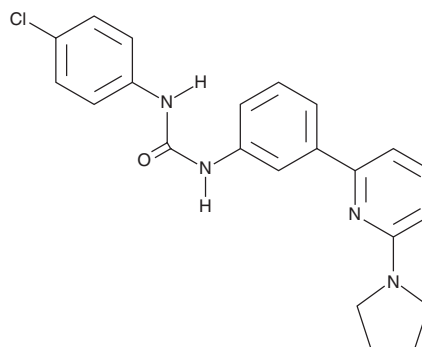
PRODUCT INFORMATION



PSNCBAM-1

Item No. 25855

CAS Registry No.: 877202-74-9
Formal Name: N-(4-chlorophenyl)-N'-[3-[6-(1-pyrrolidinyl)-2-pyridinyl]phenyl]-urea
MF: C₂₂H₂₁ClN₄O
FW: 392.9
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 260, 342 nm
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

PSNCBAM-1 is supplied as a solid. A stock solution may be made by dissolving the PSNCBAM-1 in the solvent of choice, which should be purged with an inert gas. PSNCBAM-1 is soluble in the organic solvent DMSO.

Description

PSNCBAM-1 is a cannabinoid 1 (CB₁) receptor negative allosteric modulator (NAM).¹ It is selective for CB₁ over CB₂ in HEK293 cells expressing the human receptors (IC₅₀s = 74.3 and >10,000 nM, respectively). PSNCBAM-1 inhibits activation of the CB₁ receptor induced by (±)-CP 55,940 (Item No. 13241), anandamide (AEA; Item No. 90050), or 2-arachidonoyl glycerol (2-AG; Item No. 62160) in yeast expressing human CB₁ (IC₅₀s = 45.2, 37 and 230.3 nM, respectively). It reduces food intake within one hour and decreases body weight in male rats when administered at a dose of 30 mg/kg. PSNCBAM-1 (10 mg/kg) reduces THC-induced increases in tail-flick latency in mice.²

References

1. Horswill, J.G., Bali, U., Shaaban, S., *et al.* PSNCBAM-1, a novel allosteric antagonist at cannabinoid CB₁ receptors with hypophagic effects in rats. *Br. J. Pharmacol.* **152(5)**, 805-814 (2007).
2. Gamage, T.F., Farquhar, C.E., Lefever, T.W., *et al.* The great divide: Separation between in vitro and in vivo effects of PSNCBAM-based CB₁ receptor allosteric modulators. *Neuropharmacology* **125**, 365-375 (2017).

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, 04/03/2023

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