

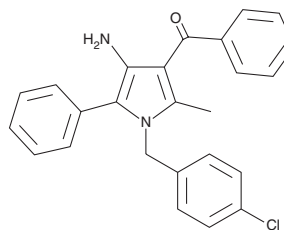
PRODUCT INFORMATION



URB447

Item No. 13261

CAS Registry No.: 1132922-57-6
Formal Name: [4-amino-1-[(4-chlorophenyl)methyl]-2-methyl-5-phenyl-1H-pyrrol-3-yl]phenyl-methanone
MF: C₂₅H₂₁ClN₂O
FW: 400.9
Purity: ≥98%
Supplied as: A crystalline 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

URB447 is supplied as a crystalline solid. A stock solution may be made by dissolving the URB447 in the solvent of choice, which should be purged with an inert gas. URB447 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of URB447 in these solvents is approximately 10 mg/ml.

Description

URB447 is a cannabinoid 1 (CB₁) receptor antagonist and CB₂ receptor agonist.¹ It inhibits the binding of the CB receptor agonist [³H]-WIN 55,212-2 to CHO-K1 cells expressing the human CB₂ receptor or rat cerebellar membranes (IC₅₀s = 41 and 313 nM, respectively). URB447 (20 mg/kg) reduces food intake and weight gain in *ob/ob* mice and reverses (R)-methanandamide-induced analgesia in mice. It decreases infarct volume and demyelination in a neonate rat model of hypoxia-ischemia-induced brain injury when administered at a dose of 1 mg/kg.²

References

- LoVerme, J., Duranti, A., Tontini, A., *et al.* Synthesis and characterization of a peripherally restricted CB₁ cannabinoid antagonist, URB447, that reduces feeding and body-weight gain in mice. *Bioorg. Med. Chem. Lett.* **19**(3), 639-643 (2009).
- Carloni, S., Crinelli, R., Palma, L., *et al.* The synthetic cannabinoid URB447 reduces brain injury and the associated white matter demyelination after hypoxia-ischemia in neonatal rats. *ACS Chem. Neurosci.* **11**(9), 1291-1299 (2020).

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, 03/11/2024

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