

PRODUCT INFORMATION

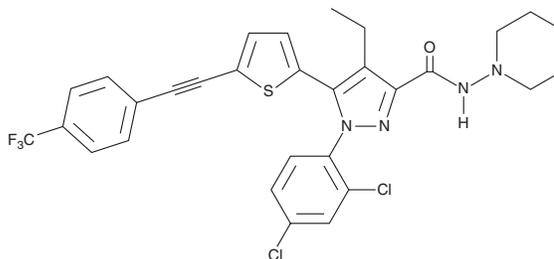


BPR0912

Item No. 10678

CAS Registry No.: 1253641-65-4
Formal Name: 1-(2,4-dichlorophenyl)-4-ethyl-N-1-piperidinyl-5-[5-[2-[4-(trifluoromethyl)phenyl]ethynyl]-2-thienyl]-1H-pyrazole-3-carboxamide

Synonym: TM38837
MF: C₃₀H₂₅Cl₂F₃N₄OS
FW: 617.5
Purity: ≥95%
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

BPR0912 is supplied as a solid. A stock solution may be made by dissolving the BPR0912 in the solvent of choice, which should be purged with an inert gas. BPR0912 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

BPR0912 is a cannabinoid 1 (CB₁) receptor inverse agonist (IC₅₀ = 8.5 nM for the human receptor).¹ It is selective for CB₁ over CB₂ receptors (IC₅₀ = 604.9 nM for the human receptor). *In vivo*, BPR0912 (10 mg/kg per day) inhibits diet-induced increases in body weight, epididymal and inguinal fat pad mass, and serum levels of insulin and triglycerides, as well as hepatic triglyceride levels, in a mouse model of high-fat diet-induced obesity.² It also increases the levels of mitochondrial uncoupling protein 1 (Ucp1) in brown adipose tissue from, as well as induces thermogenesis in, mice in the same mouse model.

References

1. Hung, M.-S., Chang, C.-P., Li, T.-C., *et al.* Discovery of 1-(2,4-dichlorophenyl)-4-ethyl-5-(5-(2-(4-(trifluoromethyl)phenyl)ethynyl)thiophen-2-yl)-N-(piperidin-1-yl)-1H-pyrazole-3-carboxamide as a potential peripheral cannabinoid-1 receptor inverse agonist. *ChemMedChem* **5**(9), 1439-1443 (2010).
2. Hsiao, W.C., Shia, K.S., Wang, Y.T., *et al.* A novel peripheral cannabinoid receptor 1 antagonist, BPR0912, reduces weight independently of food intake and modulates thermogenesis. *Diabetes Obes. Metab.* **17**(5), 495-504 (2015).

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, 04/24/2024

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