

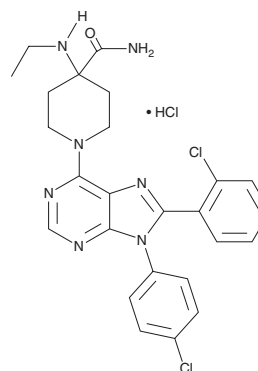
PRODUCT INFORMATION



CP 945,598 (hydrochloride)

Item No. 11955

CAS Registry No.: 686347-12-6
Formal Name: 1-[8-(2-chlorophenyl)-9-(4-chlorophenyl)-9H-purine-6-yl]-4-ethylamino-4-piperidinecarboxamide, monohydrochloride
Synonym: Otenabant
MF: C₂₅H₂₅Cl₂N₇O • HCl
FW: 546.9
Purity: ≥98%
UV/Vis.: λ_{max}: 281 nm
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

CP 945,598 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 945,598 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CP 945,598 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 0.30 mg/ml.

Description

CP 945,598 (hydrochloride) is a selective antagonist of the cannabinoid receptor CB₁ with a K_i value of 0.7 nM and demonstrates relatively low affinity for CB₂ receptors (K_i = 7.6 μM).^{1,2} It is reported to reverse cannabinoid agonist-mediated responses (i.e., hypo-locomotion, hypothermia, analgesia, and catalepsy), to reduce food intake, and to increase energy expenditure as well as fat oxidation in rats.^{1,2}

References

1. Griffith, D.A., Hadcock, J.R., Black, S.C., *et al.* Discovery of 1-[9-(4-chlorophenyl)-8-(2-chlorophenyl)-9H-purin-6-yl]-4-ethylaminopiperidine-4-carboxylic acid amide hydrochloride (CP-945,598), a novel, potent, and selective cannabinoid type 1 receptor antagonist. *J. Med. Chem.* **52**(2), 234-237 (2009).
2. Hadcock, J.R., Griffith, D.A., Iredale, P.A., *et al.* *In vitro* and *in vivo* pharmacology of CP-945,598, a potent and selective cannabinoid CB₁ receptor antagonist for the management of obesity. *Biochem. Biophys. Res. Commun.* **394**(2), 366-371 (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, 11/08/2022

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