

PRODUCT INFORMATION

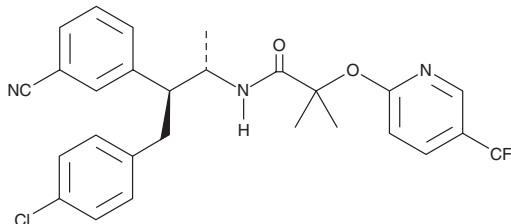


Taranabant

Item No. 10009795

CAS Registry No.: 701977-09-5
Formal Name: N-[(1S,2S)-3-(4-chlorophenyl)-2-(3-cyanophenyl)-1-methylpropyl]-2-methyl-2-[[5-(trifluoromethyl)-2-pyridinyl]oxy]-propanamide

Synonym: MK-0364
MF: C₂₇H₂₅ClF₃N₃O₂
FW: 516.0
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 270, 311 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

Taranabant is supplied as a crystalline solid. A stock solution may be made by dissolving the taranabant in the solvent of choice, which should be purged with an inert gas. Taranabant is soluble in the organic solvent DMSO at a concentration of approximately 50 mM.

Description

Taranabant is a cannabinoid (CB) receptor 1 inverse agonist ($K_i = 0.13$ nM).¹ It is selective for CB₁ over CB₂ receptors ($K_i = 170$ nM), as well as a panel of 170 receptors, enzymes, ion channels, and transporters at 1 μM. Taranabant increases forskolin-induced cAMP production in CHO cell membranes expressing human CB₁ ($EC_{50} = 2$ nM). *In vivo*, taranabant (3 mg/kg) reverses hypothermia induced by the CB receptor agonist CP 55,940 in rats. Taranabant (1 and 3 mg/kg) reduces two-hour and overnight food intake, as well as overnight weight gain in wild-type, but not CB₁ receptor-deficient, mice.

Reference

1. Fong, T.M., Guan, X.-M., Marsh, D.J., *et al.* Antiobesity efficacy of a novel cannabinoid-1 receptor inverse agonist, N-[(1S,2S)-3-(4-chlorophenyl)-2-(3-cyanophenyl)-1-methylpropyl]-2-methyl-2-[[5-(trifluoromethyl)pyridin-2-yl]oxy]propanamide (MK-0364), in rodents. *J. Pharmacol. Exp. Ther.* **321**(3), 1013-1022 (2007).

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, 12/14/2022

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