

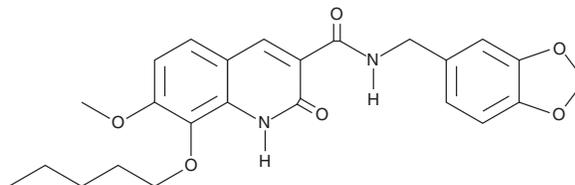
PRODUCT INFORMATION



JTE-907

Item No. 10009857

CAS Registry No.: 282089-49-0
Formal Name: N-(1,3-benzodioxol-5-ylmethyl)-1,2-dihydro-7-methoxy-2-oxo-8-(pentyloxy)-3-quinolinecarboxamide
MF: C₂₄H₂₆N₂O₆
FW: 438.5
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 264, 331 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

JTE-907 is supplied as a crystalline solid. A stock solution may be made by dissolving the JTE-907 in the solvent of choice, which should be purged with an inert gas. JTE-907 is soluble in organic solvents such as methanol, DMSO and dimethyl formamide (DMF). The solubility of JTE-907 in these solvents is approximately 1, 0.16 and 16.6 mg/ml, respectively.

JTE-907 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JTE-907 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JTE-907 has a solubility of approximately 0.2 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

JTE-907 is a cannabinoid 2 (CB₂) receptor inverse agonist.¹ It is selective for CB₂ over CB₁ receptors (K_s = 35.9 and 2,370 nM, respectively, for the human receptors). JTE-907 increases forskolin-induced cAMP production in CHO cells expressing human or mouse CB₂ receptors in a concentration-dependent manner. It promotes the differentiation of isolated mouse splenic CD4⁺ T cells into regulatory T cells (Tregs) when used at a concentration of 100 nM and decreases disease severity in a mouse model of inflammatory bowel disease (IBD) induced by dinitrobenzene sulfonic acid (DNBS).² JTE-907 (1 and 10 mg/kg) inhibits spontaneous scratching in a mouse model of chronic dermatitis, as well as reduces carrageenan-induced paw edema in mice (ED₅₀ = 0.05 mg/kg).^{1,3} Bilateral injection of JTE-907 (25 pmol/animal) into the anterior bed nucleus of the stria terminalis decreases the percentage of time spent in the open arms of the elevated plus maze in mice, indicating anxiety-like behavior.⁴

References

1. Iwamura, H., Suzuki, H., Ueda, Y., *et al.* In vitro and in vivo pharmacological characterization of JTE-907, a novel selection ligand for cannabinoid CB₂ receptor. *J. Pharmacol. Exp. Ther.* **296**(2), 420-425 (2001).
2. Gentili, M., Ronchetti, S., Ricci, E., *et al.* Selective CB2 inverse agonist JTE907 drives T cell differentiation towards a Treg cell phenotype and ameliorates inflammation in a mouse model of inflammatory bowel disease. *Pharmacol. Res.* **141**, 21-31 (2019).
3. Maekawa, T., Nojima, H., Kuraishi, Y., *et al.* The cannabinoid CB₂ receptor inverse agonist JTE-907 suppresses spontaneous itch-associated responses of NC mice, a model of atopic dermatitis. *Euro. J. Pharmacol.* **542**(1-3), 179-183 (2006).
4. Gomes-de-Souza, L., Bianchi, P.C., Costa-Ferreira, W., *et al.* CB₁ and CB₂ receptors in the bed nucleus of the stria terminalis differently modulate anxiety-like behaviors in rats. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **110**, 110284 (2021).

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, 04/04/2023

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