

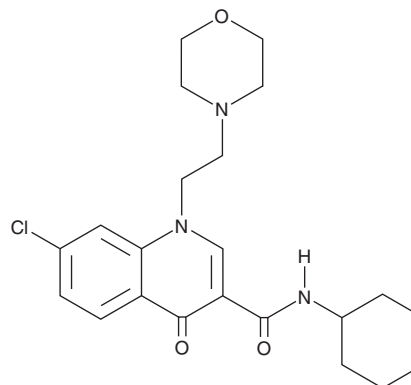
PRODUCT INFORMATION



CB-65

Item No. 43106

CAS Registry No.: 913534-05-1
Formal Name: 7-chloro-N-cyclohexyl-1,4-dihydro-1-[2-(4-morpholinyl)ethyl]-4-oxo-3-quinolinecarboxamide
MF: C₂₂H₂₈ClN₃O₃
FW: 417.9
Purity: ≥98%
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

CB-65 is supplied as a solid. A stock solution may be made by dissolving the CB-65 in the solvent of choice, which should be purged with an inert gas. CB-65 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

CB-65 is a cannabinoid 2 (CB₂) receptor inverse agonist.¹ It selectively binds to CB₂ over CB₁ receptors (K_is = 3.3 nM and >1,000 nM, respectively) and increases forskolin-induced cAMP production in CHO cells expressing the human CB₂ receptor (EC₅₀ = 1.2 nM).^{1,2} CB-65 (0.625, 2.5, and 20 nM) inhibits HepG2 hepatocellular carcinoma (HCC) cell invasion *in vitro*.³

References

1. Manera, C., Benetti, V., Castelli, M.P., *et al.* Design, synthesis, and biological evaluation of new 1,8-naphthyridin-4(1H)-on-3-carboxamide and quinolin-4(1H)-on-3-carboxamide derivatives as CB₂ selective agonists. *J. Med. Chem.* **49**(20), 5947-5957 (2006).
2. Ndong, C., O'Donnell, D., Ahmad, S., *et al.* Cloning and pharmacological characterization of the dog cannabinoid CB₂ receptor. *Eur. J. Pharmacol.* **669**(1-3), 24-31 (2011).
3. Pourkhalili, N., Ghahremani, M.H., Farsandaj, N., *et al.* Evaluation of anti-invasion effect of cannabinoids on human hepatocarcinoma cells. *Toxicol. Mech. Methods* **23**(2), 120-126 (2013).

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, 04/07/2025

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