

PRODUCT INFORMATION



CCG-203971

Item No. 15075

CAS Registry No.: 1443437-74-8
Formal Name: N-(4-chlorophenyl)-1-[3-(2-furanyl)benzoyl]-3-piperidinecarboxamide

MF: C₂₃H₂₁ClN₂O₃

FW: 408.9

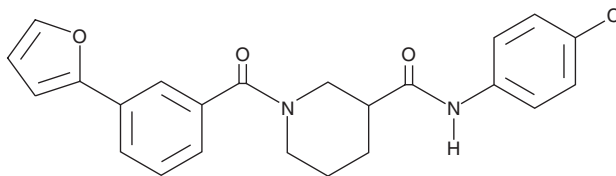
Purity: ≥90%

UV/Vis.: λ_{max}: 252, 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

CCG-203971 is supplied as a crystalline solid. A stock solution may be made by dissolving the CCG-203971 in the solvent of choice, which should be purged with an inert gas. CCG-203971 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CCG-203971 in these solvents is approximately 50, 25, and 30 mg/ml, respectively.

Description

Signaling through small G proteins of the RhoA subfamily, including RhoC, induces actin-regulated cytosolic-to-nuclear translocation of the oncogene product megakaryoblastic leukemia 1 (MKL1), which binds to serum response factor (SRF).¹ The MKL1/SRF complex, in turn, activates the transcription of serum response element (SRE) regulated genes, stimulating cell migration, a process that is central to metastasis.¹ CCG-203971 is an inhibitor of SRE activation in the prostate cancer cell line PC-3 (IC₅₀ = 6.4 μM), with 87% inhibition of SRE activation achieved at 100 μM.² This compound also inhibits PC-3 cell migration (IC₅₀ = 4.2 μM), as determined by a scratch wound assay.² CCG-203971 causes no cytotoxicity when evaluated by WST-1 assay.² It is well tolerated in normal mice up to doses of 100 mg/kg given intraperitoneally over five days.²

References

1. Evelyn, C.R., Wade, S.M., Wang, Q., *et al.* CCG-1423: A small-molecule inhibitor of RhoA transcriptional signaling. *Mol. Cancer Ther.* **6**(8), 2249-2260 (2007).
2. Bell, J.L., Haak, A.J., Wade, S.M., *et al.* Optimization of novel nipecotic bis(amide) inhibitors of the Rho/MKL1/SRF transcriptional pathway as potential anti-metastasis agents. *Bioorg. Med. Chem. Lett.* **23**(13), 3826-3832 (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, 10/06/2022

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