

PRODUCT INFORMATION



CCG-1423

Item No. 10010350

CAS Registry No.: 285986-88-1
Formal Name: N-[2-[4(4-chlorophenyl)amino]-1-methyl-2-oxoethoxy]-3,5-bis(trifluoromethyl)-benzamide

MF: C₁₈H₁₃ClF₆N₂O₃

FW: 454.8

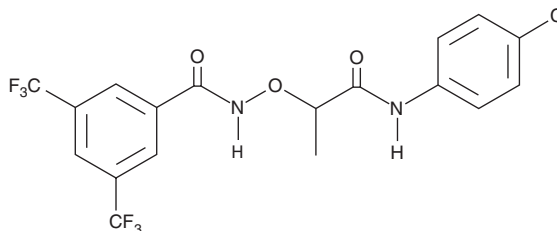
Purity: ≥98%

UV/Vis.: λ_{max}: 249 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-1423 is supplied as a crystalline solid. A stock solution may be made by dissolving the CCG-1423 in the solvent of choice, which should be purged with an inert gas. CCG-1423 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CCG-1423 in ethanol is approximately 0.25 mg/ml and approximately 10 mg/ml in DMSO and DMF.

CCG-1423 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CCG-1423 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CCG-1423 has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The Rho family of small GTPases play an important role in transduction of cell signaling events associated with several human cancers. CCG-1423 is a specific inhibitor of Rho pathway-mediated signaling and activation of serum response factor (SRF) transcription.¹ It inhibits SRF-driven luciferase expression in PC-3 cells following stimulation with constitutively active Gα₁₃ with an IC₅₀ value of 1 μM. The site of inhibition in the pathway is not precisely defined but CCG-1423 appears to act on some aspect of the interaction of SRF with its transcriptional cofactor megakaryoblastic leukemia 1 (MKL1) at a point upstream of DNA binding. CCG-1423 selectively inhibits DNA synthesis, proliferation and invasion of Rho-overexpressing cell lines at nanomolar to low micromolar concentrations.¹

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).

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