

PRODUCT INFORMATION



CCG-50014

Item No. 10802

CAS Registry No.: 883050-24-6
Formal Name: 4-[(4-fluorophenyl)methyl]-2-(4-methylphenyl)-1,2,4-thiadiazolidine-3,5-dione

MF: C₁₆H₁₃FN₂O₂S

FW: 316.3

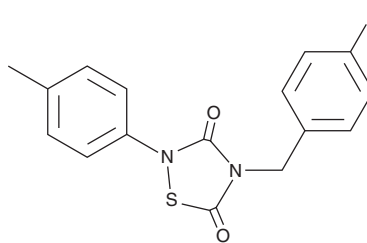
Purity: ≥98%

UV/Vis.: λ_{max}: 251 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

CCG-50014 is supplied as a crystalline solid. A stock solution may be made by dissolving the CCG-50014 in the solvent of choice. CCG-50014 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CCG-50014 in ethanol is approximately 1 mg/ml and 10 mg/ml in DMSO and DMF.

CCG-50014 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CCG-50014 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CCG-50014 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CCG-50014 is an inhibitor of regulator of G protein signaling 4 (RGS4) (IC₅₀ = 30 nM) and is >20-fold selective for RGS4 over other RGS proteins.¹ RGS proteins bind to the Gα subunits of activated heterotrimeric G proteins and accelerate the rate of GTP hydrolysis, acting as GTPase-activating proteins.¹ CCG-50014 irreversibly binds and destabilizes RGS4 with no effect on Gα_o, while blocking GTPase activity and Gα_o-dependent translocation of RGS4 to the membrane.¹

Reference

1. Blazer, L. L., Zhang, H., Casey, E. M., *et al.* A nanomolar-potency small molecule inhibitor of regulator of G protein signaling (RGS) proteins. *Biochemistry* (2011).

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, 10/04/2016

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