

PRODUCT INFORMATION

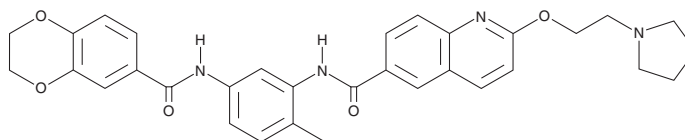


CCT251236

Item No. 27788

CAS Registry No.: 1693731-40-6
Formal Name: N-[5-[[[2,3-dihydro-1,4-benzodioxin-6-yl]carbonyl]amino]-2-methylphenyl]-2-[2-(1-pyrrolidinyl)ethoxy]-6-quinolinecarboxamide

MF: C₃₂H₃₂N₄O₅
FW: 552.6
Purity: ≥98%
UV/Vis.: λ_{max}: 244 nm
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

CCT251236 is supplied as a solid. A stock solution may be made by dissolving the CCT251236 in the solvent of choice, which should be purged with an inert gas. CCT251236 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CCT251236 in these solvents is approximately 3 mg/ml.

CCT251236 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CCT251236 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CCT251236 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CCT251236 is an inhibitor of transcription mediated by heat shock factor 1 (Hsf1) that binds to Pirin ($K_i = 28$ nM) and inhibits Hsf1-mediated induction of heat shock protein 27 (Hsp27) in U2OS cells ($IC_{50} = 19$ nM).¹ It inhibits growth of SKOV3 ovarian carcinoma cells ($GI_{50} = 1.1$ nM). CCT251236 (20 mg/kg) reduces tumor volume in an SKOV3 mouse xenograft model.

Reference

1. Cheeseman, M.D., Chessum, N.E., Rye, C.S., *et al.* Discovery of a chemical probe bisamide (CCT251236): An orally bioavailable efficacious pirin ligand from a heat shock transcription factor 1 (HSF1) phenotypic screen. *J. Med. Chem.* **60**(1), 180-201 (2017).

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, 11/01/2022

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