

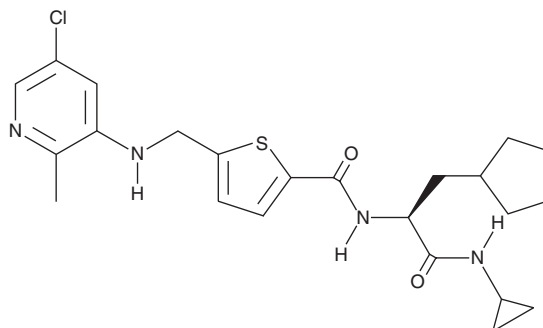
PRODUCT INFORMATION



GSK2830371

Item No. 16973

CAS Registry No.: 1404456-53-6
Formal Name: 5-[[[5-chloro-2-methyl-3-pyridinyl]amino]methyl]-N-[(1S)-1-(cyclopentylmethyl)-2-(cyclopropylamino)-2-oxoethyl]-2-thiophenecarboxamide
MF: C₂₃H₂₉ClN₄O₂S
FW: 461.0
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 255 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

GSK2830371 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2830371 in the solvent of choice, which should be purged with an inert gas. GSK2830371 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK2830371 in ethanol and DMSO is approximately 25 mg/ml and approximately 30 mg/ml in DMF.

GSK2830371 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK2830371 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GSK2830371 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK2830371 is a potent inhibitor of Wip1 (IC₅₀ = 6 nM).¹ It displays selectivity for Wip1 over 21 other phosphatases.¹ GSK2830371 increases phosphorylation of Wip1 substrates and blocks cell cycling in hematopoietic cancer cells and in Wip1-amplified cancer cells with wild-type p53.¹ This compound is orally bioavailable and causes inhibition of lymphoma xenograft growth in mice.¹

Reference

1. Gilmartin, A.G., Faitg, T.H., Richter, M., *et al.* Allosteric Wip1 phosphatase inhibition through flap-subdomain interaction. *Nat. Chem. Biol.* **10**(3), 181-187 (2014).

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, 12/19/2022

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