

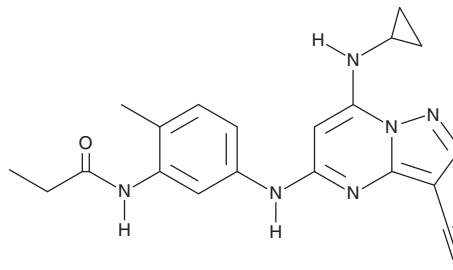
PRODUCT INFORMATION



SGC-CK2-1

Item No. 34103

CAS Registry No.: 2470424-39-4
Formal Name: N-[5-[[3-cyano-7-(cyclopropylamino)pyrazolo[1,5-a]pyrimidin-5-yl]amino]-2-methylphenyl]-propanamide
Synonym: SGC-Casein Kinase 2-1
MF: C₂₀H₂₁N₇O
FW: 375.4
Purity: ≥98%
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

SGC-CK2-1 is supplied as a solid. A stock solution may be made by dissolving the SGC-CK2-1 in the solvent of choice, which should be purged with an inert gas. SGC-CK2-1 is soluble in DMSO. SGC-CK2-1 is slightly soluble in acetonitrile.

Description

SGC-CK2-1 is an inhibitor of casein kinase 2 α (CK2 α ; IC₅₀s = 4.2 and 2.3 nM for the catalytic subunits CK2 α and CK2 α' , respectively).¹ It is selective for CK2 α over DYRK2, Plk4, MEK5, and HIPK1, -2, and -3 (IC₅₀s = 440, >10,000, 1,000, 3,700, 3,400, and 8,100 nM, respectively). SGC-CK2-1 (0.01-10 μ M) decreases phosphorylation of Akt in HCT116 cells. See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Wells, C.I., Drewry, D.H., Pickett, J.E., *et al.* Development of a potent and selective chemical probe for the pleiotropic kinase CK2. *Cell Chem. Biol.* **28(4)**, 546-558 (2021).

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, 11/06/2023

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