

PRODUCT INFORMATION



SGC-CLK-1

Item No. 41311

CAS Registry No.: 748142-15-6

Formal Name: 4-(6-methoxy-1,2,3,4-tetrahydropyrazolo[1,5-b]pyridazin-3-yl)-N-[3-methoxy-5-(trifluoromethyl)phenyl]-2-pyrimidinamine

Synonyms: CAF-170, CDC-like Kinase 1

MF: $C_{19}H_{15}F_3N_6O_2$

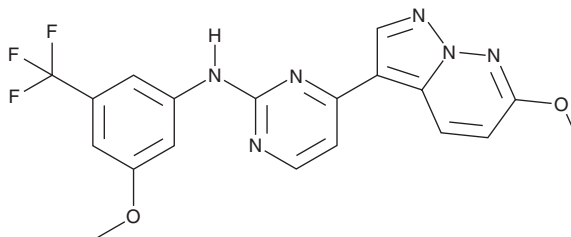
FW: 416.4

Purity: $\geq 98\%$

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-CLK-1 is supplied as a solid. A stock solution may be made by dissolving the SGC-CLK-1 in the solvent of choice, which should be purged with an inert gas. SGC-CLK-1 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SGC-CLK-1 is an inhibitor of CDC-like kinase 1 (Clk1) and Clk2 (IC_{50} s = 13 and 4 nM, respectively).¹ It is selective for Clk1 and Clk2 over Clk3, Clk4, homeodomain-interacting protein kinase 1 (HIPK1), HIPK2, and serine/threonine kinase 16 (STK16; IC_{50} s = 363, 46, 50, 42, and 49 nM, respectively). SGC-CLK-1 (1 μM) inhibits the proliferation of MDA-MB-435 breast cancer and U-118 MG glioma cells. It induces diffuse localization of Clk2 and phosphorylated serine/arginine-rich (pSR) proteins, targets of Clks, without reducing the levels of phosphorylated SR proteins in MDA-MB-435 and U-118 MG cells when used at a concentration of 500 nM. See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Tiek, D., Wells, C.I., Schröder, M., et al. SGC-CLK-1: A chemical probe for the Cdc2-like kinases CLK1, CLK2, and CLK4. *Curr. Res. Chem. Biol.* **3**, 100045 (2023).

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, 08/05/2024

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