

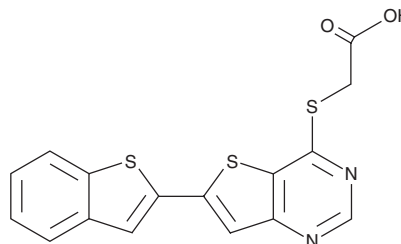
PRODUCT INFORMATION



SGC-STK17B-1

Item No. 41126

CAS Registry No.: 2650530-00-8
Formal Name: 2-[(6-benzo[b]thien-2-ylthieno[3,2-d]pyrimidin-4-yl)thio]-acetic acid
Synonym: SGC-Serine/Threonine Kinase 17b-1
MF: C₁₆H₁₀N₂O₂S₃
FW: 358.5
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-STK17B-1 is supplied as a solid. A stock solution may be made by dissolving the SGC-STK17B-1 in the solvent of choice, which should be purged with an inert gas. SGC-STK17B-1 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

SGC-STK17B-1 is an inhibitor of serine/threonine kinase 17b (STK17B), also known as DRAK2 (IC₅₀ = 43 nM).¹ It is selective for STK17B over STK17A (IC₅₀ = 4,700 nM), as well as eight additional kinases (IC₅₀s = 5,200->10,000 nM). See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Picado, A., Chaikuad, A., Wells, C.I., *et al.* A chemical probe for dark kinase STK17B derives its potency and high selectivity through a unique P-loop conformation. *J. Med. Chem.* **63**(23), 14626-14646 (2020).

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, 06/03/2024

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