

PRODUCT INFORMATION



SGC-AAK1-1

Item No. 41118

CAS Registry No.: 2247894-32-0

Formal Name: N-[6-[3-[[[(diethylamino)sulfonyl]amino]phenyl]-1H-indazol-3-yl]-cyclopropanecarboxamide]

MF: $C_{21}H_{25}N_5O_3S$

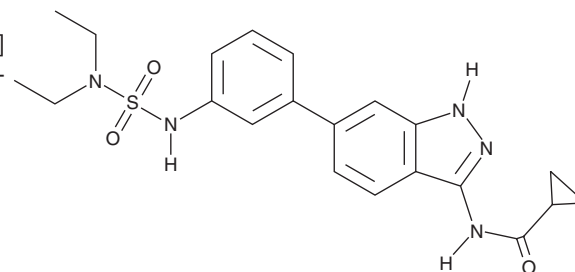
FW: 427.5

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-AAK1-1 is supplied as a solid. A stock solution may be made by dissolving the SGC-AAK1-1 in the solvent of choice, which should be purged with an inert gas. SGC-AAK1-1 is sparingly soluble (1-10 mg/ml) in the organic solvent DMSO.

Description

SGC-AAK1-1 is a dual inhibitor of AP2-associated protein kinase 1 (AAK1) and BMP2-inducible kinase (BMP2K; K_i s = 9.1 and 17 nM, respectively).¹ It is selective for AAK1 and BMP2K over cyclin G-associated kinase (GAK) and serine/threonine kinase 16 (STK16; K_i s = 1,700 and 270 nM, respectively). See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Wells, C., Couñago, R.M., Limas, J.C., *et al.* SGC-AAK1-1: A chemical probe targeting AAK1 and BMP2K. *ACS Med. Chem. Lett.* **11**(3), 340-345 (2019).

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, 06/26/2024

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