

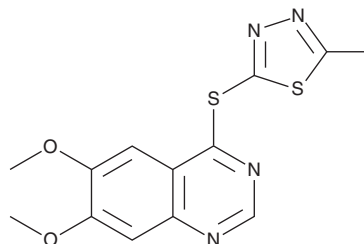
PRODUCT INFORMATION



SKLB1002

Item No. 34422

CAS Registry No.: 1225451-84-2
Formal Name: 6,7-dimethoxy-4-[(5-methyl-1,3,4-thiadiazol-2-yl)thio]-quinazoline
MF: C₁₃H₁₂N₄O₂S₂
FW: 320.4
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 249, 276, 342 nm
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

SKLB1002 is supplied as a solid. A stock solution may be made by dissolving the SKLB1002 in the solvent of choice, which should be purged with an inert gas. SKLB1002 is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

SKLB1002 is a VEGFR2 inhibitor (IC₅₀ = 32 nM).¹ It is selective for VEGFR2 over a panel of 16 additional kinases (IC₅₀s = 619->10,000 nM). SKLB1002 inhibits tube formation, migration, and VEGF-induced proliferation of human umbilical vein endothelial cells (HUVECs). It induces tumor cell apoptosis, decreases tumor microvessel density, and reduces tumor growth in an SW620 mouse xenograft model when administered at a dose of 50 or 100 mg/kg.

Reference

1. Zhang, S., Cao, Z., Tian, H., *et al.* SKLB1002, a novel potent inhibitor of VEGF receptor 2 signaling, inhibits angiogenesis and tumor growth *in vivo*. *Clin. Cancer Res.* **17(13)**, 4439-4450 (2011).

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, 11/22/2022

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