

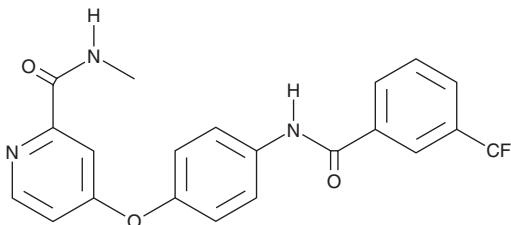
# PRODUCT INFORMATION



## SKLB610

Item No. 21561

**CAS Registry No.:** 1125780-41-7  
**Formal Name:** N-methyl-4-[4-[[3-(trifluoromethyl)benzoyl]amino]phenoxy]-2-pyridinecarboxamide  
**MF:** C<sub>21</sub>H<sub>16</sub>F<sub>3</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 415.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 272 nm  
**Supplied as:** A crystalline 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

SKLB610 is supplied as a crystalline solid. A stock solution may be made by dissolving the SKLB610 in the solvent of choice, which should be purged with an inert gas. SKLB610 is soluble in organic solvents such as ethanol, DMSO and dimethyl formamide (DMF). The solubility of SKLB610 in these solvents is approximately 5, 16, and 20 mg/ml, respectively.

SKLB610 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SKLB610 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SKLB610 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

SKLB610 is an inhibitor of VEGF receptor 2 (VEGFR2).<sup>1</sup> It inhibits VEGFR2 activity by 97% but also inhibits FGFR2 and PDGFRβ activity by 65 and 55%, respectively, when used at a concentration of 10 μM. It is selective for VEGFR2, FGFR2, and PDGFRβ over PI3K, EGFR, Aurora A, Cdk2/cyclin E, and Cdk6/cyclin D3 at 10 μM. SKLB610 inhibits phosphorylation of VEGFR2 induced by VEGF in human umbilical vein endothelial cells (HUVECs). It inhibits proliferation of HUVECs induced by VEGF and basic FGF (bFGF; IC<sub>50</sub>s = 2.2 and 4.7 μM, respectively). It also inhibits HUVEC capillary tube formation and migration when used at concentrations of 2.5 and 10 μM, respectively. SKLB610 inhibits proliferation of a variety of cancer cells, including A549 human lung cancer, HCT116 human colorectal carcinoma, MDA-MB-231 human mammary carcinoma, Raji human Burkitt's Lymphoma, and DU145 human prostate cancer cells (IC<sub>50</sub>s = 5.7, 5.3, 25.6, 6.4, and 6.3 μM, respectively). It reduces tumor growth in A549 and HCT116 mouse xenograft models when administered at a dose of 50 mg/kg per day.

### Reference

1. Cao, Z.X., Zheng, R.L., and Lin, H.J. SKLB610: A novel potential inhibitor of vascular endothelial growth factor receptor tyrosine kinases inhibits angiogenesis and tumor growth *in vivo*. *Cell Physiol. Biochem.* 27(5), 565-574 (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, 12/16/2022

#### 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