

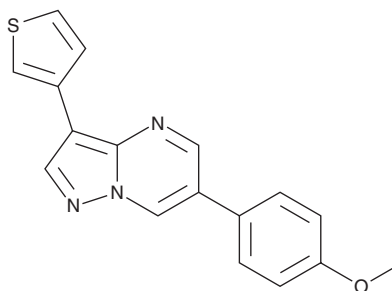
PRODUCT INFORMATION



VEGFR2 Kinase Inhibitor IV

Item No. 45474

CAS Registry No.: 216661-57-3
Formal Name: 6-(4-methoxyphenyl)-3-(3-thienyl)-pyrazolo[1,5-a]pyrimidine
Synonym: Vascular Endothelial Growth Factor Receptor 2 Kinase Inhibitor IV
MF: C₁₇H₁₃N₃OS
FW: 307.4
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

VEGFR2 kinase inhibitor IV is supplied as a solid. A stock solution may be made by dissolving the VEGFR2 kinase inhibitor IV in the solvent of choice, which should be purged with an inert gas. VEGFR2 kinase inhibitor IV is sparingly soluble (1-10 mg/ml) in 1:1 solution of methanol:chloroform and slightly soluble (0.1-1 mg/ml) in DMSO.

Description

VEGFR2 kinase inhibitor IV is an inhibitor of VEGFR2 (IC₅₀ = 19 nM).¹ It is greater than 100-fold selective for VEGFR2 over FGFR1 and Src but only 1.8-, 10-, and 10-fold selective over PDGFRβ, FLT1, and FLT2, respectively. VEGFR2 kinase inhibitor IV inhibits mitogenesis in human umbilical vein endothelial cells (HUVECs; IC₅₀ = 387 nM). It also reduces VEGF-induced increases in neuroglobin levels in primary mouse cerebrocortical neurons when used at a concentration of 20 nM.²

References

1. Fraley, M.E., Hoffman, W.F., Rubino, R.S., *et al.* Synthesis and initial SAR studies of 3,6-disubstituted pyrazolo[1,5-a]pyrimidines: A new class of KDR kinase inhibitors. *Bioorg. Med. Chem. Lett.* **12(19)**, 2767-2770 (2002).
2. Jin, K., Mao, X., Xie, L., *et al.* Interactions between vascular endothelial growth factor and neuroglobin. *Neurosci. Lett.* **519(1)**, 47-50 (2012).

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, 04/24/2026

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