

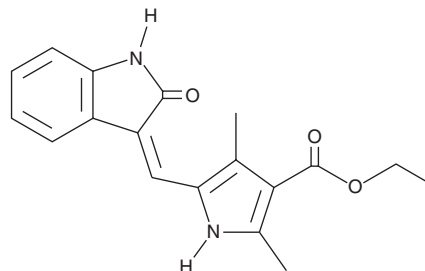
PRODUCT INFORMATION



VEGFR2 Kinase Inhibitor I

Item No. 13578

CAS Registry No.: 15966-93-5
Formal Name: 5-[(1,2-dihydro-2-oxo-3H-indol-3-ylidene)methyl]-2,4-dimethyl-1H-pyrrole-3-carboxylic acid, ethyl ester
Synonym: SU 5408
MF: C₁₈H₁₈N₂O₃
FW: 310.4
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 266, 418 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

VEGFR2 Kinase inhibitor I is supplied as a crystalline solid. A stock solution may be made by dissolving the VEGFR2 kinase inhibitor I in the solvent of choice, which should be purged with an inert gas. VEGFR2 Kinase inhibitor I is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of VEGFR2 kinase inhibitor I in these solvents is approximately 1 mg/ml.

Description

Vascular endothelial growth factor receptor 2 (VEGFR2, also known as KDR and FLK1) is a receptor tyrosine kinase that regulates angiogenesis, vascular development, and embryonic hematopoiesis in response to VEGF isoforms A, C, and D. VEGFR2 kinase inhibitor I is a potent, cell-permeable inhibitor of mouse VEGFR2 kinase (IC₅₀ = 70 nM).¹ It has little or no effect against receptors for platelet-derived growth factor, epidermal growth factor, or insulin-like growth factor (IC₅₀ > 100 μM).¹

Reference

1. Sun, L., Tran, N., Tang, F., *et al.* Synthesis and biological evaluations of 3-substituted indolin-2-ones: A novel class of tyrosine kinase inhibitors that exhibit selectivity toward particular receptor tyrosine kinases. *J. Med. Chem.* **41**(14), 2588-2603 (1998).

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, 10/03/2022

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