

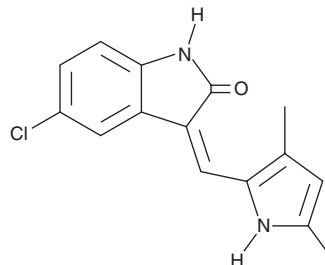
PRODUCT INFORMATION



SU 5614

Item No. 17410

CAS Registry No.: 1055412-47-9
Formal Name: (3Z)-5-chloro-3-[(3,5-dimethyl-1H-pyrrol-2-yl)methylene]-1,3-dihydro-2H-indol-2-one
MF: C₁₅H₁₃ClN₂O
FW: 272.7
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 280, 445 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

SU 5614 is supplied as a crystalline solid. A stock solution may be made by dissolving the SU 5614 in the solvent of choice, which should be purged with an inert gas. SU 5614 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SU 5614 in these solvents is approximately 0.25, 1, and 15 mg/ml, respectively.

Description

SU 5614 is an inhibitor of rearranged during transfection (RET) kinase with an IC₅₀ value of 170 nM.¹ It inhibits proliferation of, and RET/PTC2 autophosphorylation in, RET expressing Ba/F3 cells in a dose-dependent manner. It also inhibits growth of TPC-1 human papillary carcinoma cells with an IC₅₀ value of 2.7 μM. *In vivo*, SU 5614 inhibits RET autophosphorylation in a mouse NIH-PTC xenograft model. SU 5614 also inhibits the VEGF receptor (VEGFR; IC₅₀ = 1.23 μM) but has no effect on EGF or FGF kinase activities.² It inhibits VEGF-induced mitogenesis of human umbilical vein endothelial cells (HUVECs) *in vitro* and inhibits growth of A375 tumors in a mouse xenograft model.

References

1. Mologni, L., Sala, E., Cazzaniga, S., *et al.* Inhibition of RET tyrosine kinase by SU5416. *J. Mol. Endocrinol.* **37**(2), 199-212 (2006).
2. Fong, T.A.T., Shawver, L.K., Sun, L., *et al.* SU5416 is a potent and selective inhibitor of the vascular endothelial growth factor receptor (Flk-1/KDR) that inhibits tyrosine kinase catalysis, tumor vascularization, and growth of multiple tumor types. *Cancer Res.* **59**(1), 99-106 (1999).

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, 11/08/2022

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