

PRODUCT INFORMATION



SU 16f

Item No. 19555

CAS Registry No.: 251356-45-3

Formal Name: 5-[(1,2-dihydro-2-oxo-6-phenyl-3H-indol-3-ylidene)methyl]-2,4-dimethyl-1H-pyrrole-3-propanoic acid

Synonym: PDGFR Tyrosine Kinase Inhibitor VII

MF: $C_{24}H_{22}N_2O_3$

FW: 386.4

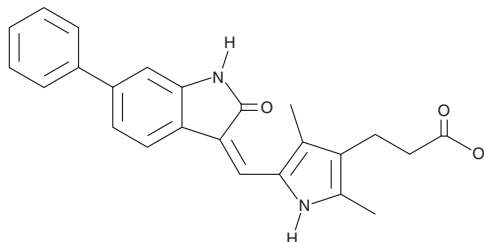
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 291, 459 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SU 16f is supplied as a crystalline solid. A stock solution may be made by dissolving the SU 16f in the solvent of choice. SU 16f is soluble in the organic solvent DMSO, which should be purged with an inert gas. The solubility of SU 16f in DMSO is 100 mM.

Description

SU 16f is a potent inhibitor of platelet-derived growth factor receptor β (PDGFR β) with an IC_{50} value of 10 nM.¹ It also inhibits VEGF receptor 2 (VEGFR2; IC_{50} = 140 nM). SU 16f selectively inhibits PDGF-over VEGF-, FGF-, and EGF-induced cell proliferation (IC_{50} s = 0.11, 10, 10, and 21.9 μ M, respectively). It also accelerates downregulation of fibroblast genes and increases the yield of beating clusters in human foreskin fibroblasts (HFFs) treated with 15 compounds to induce a cardiac myocyte-like phenotype.²

References

1. Sun, L., Tran, N., Liang, C., *et al.* Design, synthesis, and evaluations of substituted 3-[(3- or 4-carboxyethylpyrrol-2-yl)methylidenyl]indolin-2-ones as inhibitors of VEGF, FGF, and PDGF receptor tyrosine kinases. *J. Med. Chem.* **42**(25), 5120-5130 (1999).
2. Cao, N., Huang, Y., Zheng, J., *et al.* Conversion of human fibroblasts into functional cardiomyocytes by small molecules. *Science* **352**(6290), 1216-1220 (2016).

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, 12/22/2022

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