

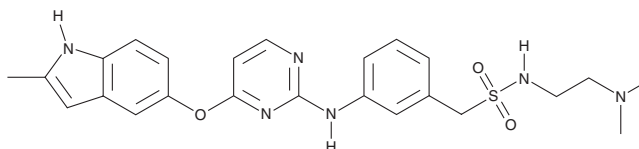
PRODUCT INFORMATION



Sulfatinib

Item No. 34206

CAS Registry No.: 1308672-74-3
Formal Name: N-[2-(dimethylamino)ethyl]-3-[[4-[(2-methyl-1H-indol-5-yl)oxy]-2-pyrimidinyl]amino]-benzenesulfonamide
Synonym: HMPL-012
MF: C₂₄H₂₈N₆O₃S
FW: 480.6
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 273 nm
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

Sulfatinib is supplied as a solid. A stock solution may be made by dissolving the sulfatinib in the solvent of choice, which should be purged with an inert gas. Sulfatinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of sulfatinib in these solvents is approximately 1, 10, and 30 mg/ml, respectively.

Sulfatinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sulfatinib should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Sulfatinib has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfatinib is a multi-kinase inhibitor.¹ It inhibits the receptor tyrosine kinases VEGFR1, VEGFR2, VEGFR3, FGFR1, colony stimulating factor 1 receptor (CSF1R), tropomyosin-related kinase B (TrkB), and FMS-related tyrosine kinase 3 (FLT3; IC₅₀s = 2, 24, 1, 15, 4, 41, and 67 nM, respectively). It is selective for these kinases over a panel of 278 kinases (IC₅₀s = >150 nM for all).

Reference

1. Qi, C. and Xu, J. Method of treating solid tumors. *Hutchison Medipharma Limited*. **WO2018090324A1** (2018).

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, 11/22/2022

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