

PRODUCT INFORMATION



Sorafenib

Item No. 10009644

CAS Registry No.: 284461-73-0

Formal Name: 4-[4-[[[4-chloro-3-(trifluoromethyl)phenyl]amino]carbonyl]amino]phenoxy]-N-methyl-2-pyridinecarboxamide

Synonym: BAY 43-9006

MF: $C_{21}H_{16}ClF_3N_4O_3$

FW: 464.8

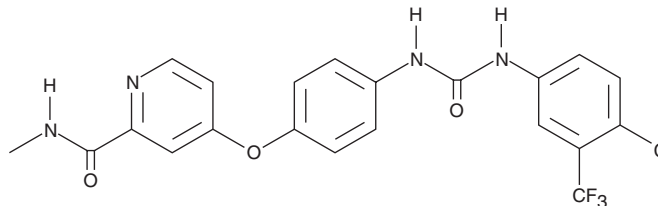
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 204, 266 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

Sorafenib is supplied as a crystalline solid. A stock solution may be made by dissolving the sorafenib in the solvent of choice, which should be purged with an inert gas. Sorafenib is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sorafenib in these solvents is approximately 20 mg/ml.

Sorafenib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sorafenib should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Sorafenib has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Sorafenib is a multi-kinase inhibitor that inhibits Raf-1 and B-RAF ($IC_{50}s = 6$ and $22 \mu M$, respectively), as well as the receptor tyrosine kinases VEGFR2, VEGFR3, PDGFR β , FLT3, and c-Kit ($IC_{50}s = 90, 15, 20, 57,$ and 58 nM, respectively).^{1,2} It is selective for these kinases over 12 other kinases, including ERK1, MEK1, EGFR, and HER2 ($IC_{50}s = >10 \mu M$ for all).² Sorafenib inhibits proliferation of PLC/PRF/5 and HepG2 cells ($IC_{50}s = 6.3$ and $4.5 \mu M$, respectively) and induces apoptosis in these cells.³ It completely inhibits tumor growth in a PLC/PRF/5 mouse xenograft model when administered at a dose of 30 mg/kg and reduces basic FGF-induced angiogenesis in a Matrigel™ assay *in vivo*.^{3,4} Sorafenib ($10 \mu M$) induces ferroptotic cell death in HT-1080 fibrosarcoma cells, an effect that can be blocked by the ferroptosis inhibitors ferrostatin-1 (Item No. 17729), deferoxamine (Item No. 14595), and β -mercaptoethanol, but does not induce ferroptosis in a variety of other cancer cell lines.^{5,7} It inhibits glutamate release by the system x_c^- cystine/glutamate transporter in HT-1080 cells when used at concentrations ranging from 2.5 to $10 \mu M$, decreases glutathione levels, and increases lipid peroxidation.⁵ Sorafenib also inhibits replication of hepatitis C virus (HCV) in Huh7.5 cells ($IC_{50} = 7.2 \mu M$).⁶ Formulations containing sorafenib have been used in the treatment of hepatocellular, renal cell, and thyroid carcinomas.

References

1. Lyons, J.F., Wilhelm, S., Hibner, B., *et al.* *Endocr. Relat. Cancer* **8**(3), 219-225 (2001).
2. Wilhelm, S.M., Carter, C., and Tang, L. *Cancer Res.* **64**(19), 7099-7109 (2004).
3. Liu, L., Cao, Y., Chen, C., *et al.* *Cancer Res.* **66**(24), 11851-11858 (2006).
4. Murphy, D.A., Makonnen, S., Lassoued, W., *et al.* *Am. J. Pathol.* **169**(5), 1875-1885 (2006).
5. Dixon, S.J., Patel, D.N., Welsch, M., *et al.* *Elife* **3**, e02523 (2014).
6. Himmelsbach, K., Sauter, D., Baumert, T.F., *et al.* *Gut* **58**(12), 1644-1653 (2009).
7. Zheng, J., Sato, M., Mishima, E., *et al.* *Cell Death Dis.* **12**(7), 698 (2021).

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, 09/29/2022

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