

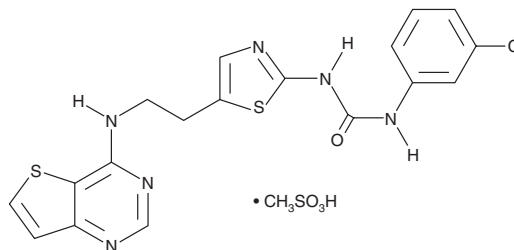
PRODUCT INFORMATION



SNS-314 (mesylate)

Item No. 14974

CAS Registry No.: 1146618-41-8
Formal Name: N-(3-chlorophenyl)-N'-[5-[2-(thieno[3,2-d]pyrimidin-4-ylamino)ethyl]-2-thiazoly]-urea, monomethanesulfonate
MF: C₁₈H₁₅ClN₆OS₂ • CH₃SO₃H
FW: 527.0
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 275 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

SNS-314 (mesylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the SNS-314 (mesylate) in the solvent of choice. SNS-314 (mesylate) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of SNS-314 (mesylate) in these solvents is approximately 20 mg/ml.

SNS-314 (mesylate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SNS-314 (mesylate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SNS-314 (mesylate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SNS-314 is a pan-Aurora kinase inhibitor (IC₅₀s = 9, 31, and 3.4 nM for Aurora A, B, and C, respectively).^{1,2} In a panel of 219 kinases, SNS-314 also inhibits TRKB, TRKA, FLT4, FMS, DDR2, AXL, and C-RAF (IC₅₀s = 5-100 nM).¹ SNS-314 inhibits Aurora B phosphorylation of histone H3 in HCT116 human colorectal carcinoma cells *in vitro* (EC₅₀ = <16 nM) and in a mouse xenograft model at a dose of 50 mg/kg. It also inhibits tumor growth in A2780 ovarian, A375 melanoma, H1299 lung, MDA-MB-231 breast, and PC3 prostate cancer mouse xenograft models when administered biweekly for six weeks at a dose of 170 mg/kg.²

References

- Oslob, J.D., Romanowski, M.J., Allen, D.A., *et al.* Discovery of a potent and selective aurora kinase inhibitor. *Bioorg. Med. Chem. Lett.* **18(17)**, 4880-4884 (2008).
- Arbitrario, J.P., Belmont, B.J., Evanchik, M.J., *et al.* SNS-314, a pan-Aurora kinase inhibitor, shows potent anti-tumor activity and dosing flexibility *in vivo*. *Cancer Chemother. Pharmacol.* **65(4)**, 707-717 (2010).

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, 11/23/2022

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