

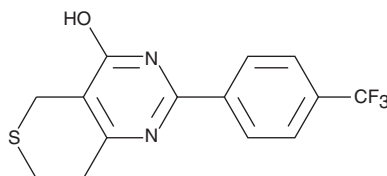
PRODUCT INFORMATION



XAV939

Item No. 13596

CAS Registry No.: 284028-89-3
Formal Name: 3,5,7,8-tetrahydro-2-[4-(trifluoromethyl)phenyl]-4H-thiopyrano[4,3-d]pyrimidin-4-one
MF: C₁₄H₁₁F₃N₂OS
FW: 312.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20C
Stability: ≥4 years



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

Laboratory Procedures

XAV939 is supplied as a crystalline solid. A stock solution may be made by dissolving the XAV939 in the solvent of choice, which should be purged with an inert gas. XAV939 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of XAV939 in DMSO is approximately 2 mg/ml and approximately 5 mg/ml in DMF.

XAV939 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, XAV939 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. XAV939 has a solubility of approximately 0.15 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The Wnt signaling pathway is integral to normal biological processes and inappropriately active in many cancers. It is regulated through degradation of the downstream effector β -catenin via a complex consisting of the tumor suppressor APC, axin, and glycogen synthase kinase 3 (GSK3). Axin is the concentration-limiting factor for this degradation complex. Tankyrases, initially identified as telomere-associated proteins¹, promote axin ubiquitination, possibly through poly-ADP-ribosylation (PARsylation).² XAV939 is a potent, small molecule inhibitor of tankyrase (TNKS) 1 and 2 with IC₅₀ values of 11 and 4 nM, respectively.² By inhibiting TNKS activity, XAV939 increases the protein levels of the axin-GSK3 β complex and promotes the degradation of β -catenin in SW480 cells.² At concentrations as low as 0.33 μ M, XAV939 inhibits colony formation of APC-deficient colorectal cancer cells.²

References

1. Smith, S., Gariat, I., Schmitt, A., *et al.* Tankyrase, a poly(ADP-ribose) polymerase at human telomeres. *Science* **282**, 1484-1487 (1998).
2. Huang, S.-M.A., Mishina, Y.M., Liu, S., *et al.* Tankyrase inhibition stabilizes axin and antagonizes Wnt signalling. *Nature* **461**, 614-619 (2009).

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, 11/17/2022

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