

XAV-939

Catalog #: 27100

Lot #: 111121

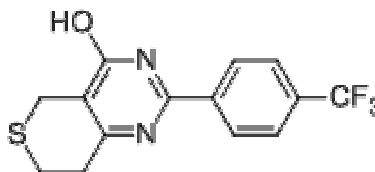
Size: 10 mg

Structure:

CAS Registry #: 284028-89-3

Purity: >99%

Chemical Formula: C₁₄H₁₁F₃N₂OS



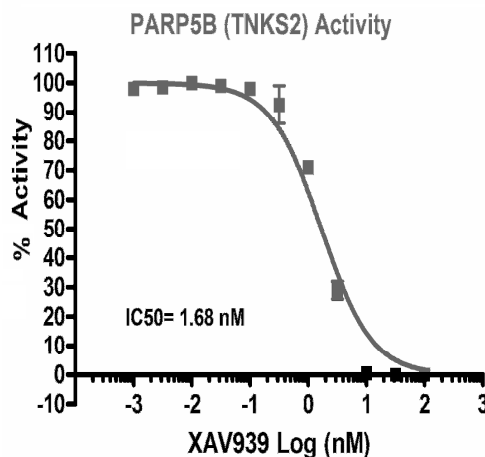
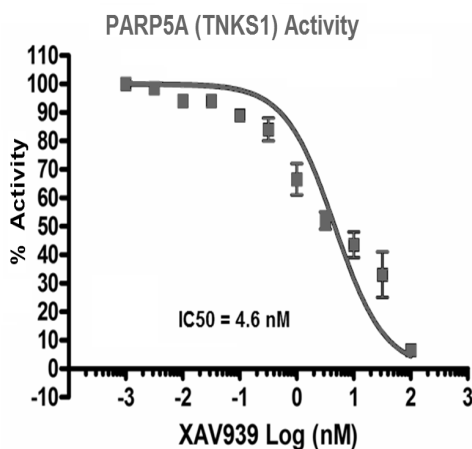
Molecular Weight: 312.31

Description: XAV 939 is a small molecule and selective Wnt pathway transcription factor β -catenin-mediated transcription inhibitor and axin stabilizing agent. It stimulates beta-catenin degradation by stabilizing axin, the concentration-limiting component of the destruction complex and stabilizes axin by inhibiting the poly-ADP-ribosylating enzymes tankyrase 1 (TNKS1) and tankyrase 2 (TNKS2).

Appearance: White to beige powder.

Solubility: Soluble in DMSO.

Biological Activity: XAV939 inhibited TNKS1 and TNKS2 with IC₅₀ of 4.6 nM and 1.68 nM, respectively using BPS Bioscience TNKS1 Assay Kit (Cat. # 80564) and TNKS2 Assay Kit (Cat. # 80566).



Storage/Stability: Store at or below -20 °C.

Quality Control: The purity was determined by HPLC analysis.

Reference: Huang, SM. *et al. Nature*. 2009 Oct 1; **461(7264)**:614-20.