

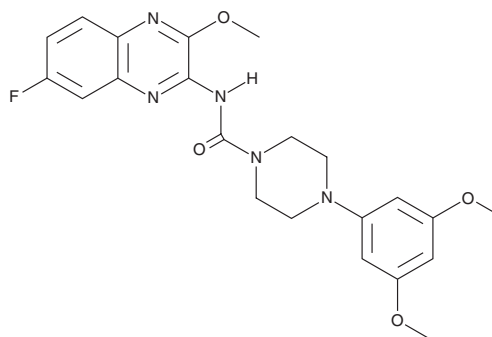
PRODUCT INFORMATION



Supinoxin

Item No. 36573

CAS Registry No.: 888478-45-3
Formal Name: 4-(3,5-dimethoxyphenyl)-N-(7-fluoro-3-methoxy-2-quinoxaliny)-1-piperazinecarboxamide
Synonym: RX-5902
MF: C₂₂H₂₄FN₅O₄
FW: 441.5
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 252 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

Supinoxin is supplied as a solid. A stock solution may be made by dissolving the Supinoxin in the solvent of choice, which should be purged with an inert gas. Supinoxin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Supinoxin in these solvents is approximately 10 and 25 mg/ml, respectively.

Description

Supinoxin is an inhibitor of Y593-phosphorylated DEAD-box helicase 5 (DDX5).¹ It binds to Y593-phosphorylated DDX5 (K_d = ~19 nM), but does not interact with unphosphorylated DDX5, and selectively inhibits β-catenin-dependent ATPase activity over RNA-dependent ATPase activity at 0.2 μM. Supinoxin induces apoptosis in MDA-MB-231 cells (IC₅₀ = 12 nM) and inhibits cell growth in SK-MEL-28 cells (IC₅₀ = 20 nM).

Reference

1. Kost, G.C., Yang, M.Y., Li, L., *et al.* A novel anti-cancer agent, 1-(3,5-dimethoxyphenyl)-4-[(6-fluoro-2-methoxyquinoxalin-3-yl)aminocarbonyl] piperazine (RX-5902), interferes with β-catenin function through Y593 phospho-p68 RNA helicase. *J. Cell. Biochem.* **116**(8), 1595-1601 (2015).

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, 05/31/2023

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