

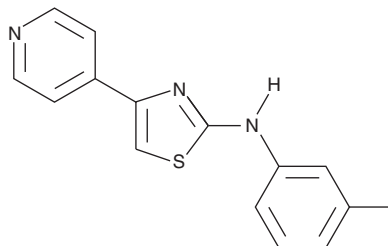
PRODUCT INFORMATION



STF-62247

Item No. 13084

CAS Registry No.: 315702-99-9
Formal Name: N-(3-methylphenyl)-4-(4-pyridinyl)-2-thiazolamine
MF: C₁₅H₁₃N₃S
FW: 267.3
Purity: ≥98%
UV/Vis.: λ_{max}: 246, 296 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

STF-62247 is supplied as a crystalline solid. A stock solution may be made by dissolving the STF-62247 in the solvent of choice, which should be purged with an inert gas. STF-62247 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of STF-62247 in these solvents is approximately 25 mg/ml.

STF-62247 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, STF-62247 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. STF-62247 has a solubility of approximately 0.2 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

STF-62247 is a small molecule agonist that induces autophagy and selectively causes lethality in renal cell carcinoma (RCC) cells that have lost the von Hippel-Lindau (VHL) tumor suppressor activity (IC₅₀ = 625 nM).¹ It significantly reduces the growth rate of tumors formed from VHL-deficient cells in mice.¹⁻³ STF-62247 induces cytotoxicity in VHL-positive RCC cells only at higher concentrations (IC₅₀ = 16 μM).¹

References

1. Turcotte, S., Sutphin, P.D., and Giaccia, A.J. Targeting therapy for the loss of von Hippel-Lindau in renal cell carcinoma; a novel molecule that induces autophagic cell death. *Autophagy* **4**(7), 944-946 (2008).
2. Turcotte, S., Chan, D.A., Sutphin, P.D., et al. A molecule targeting VHL-deficient renal cell carcinoma that induces autophagy. *Cancer Cell* **14**(1), 90-102 (2008).
3. Chan, D.A. and Giaccia, A.J. Targeting cancer cells by synthetic lethality; autophagy and VHL in cancer therapeutics. *Cell Cycle* **7**(19), 2987-2990 (2008).

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, 11/04/2022

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