

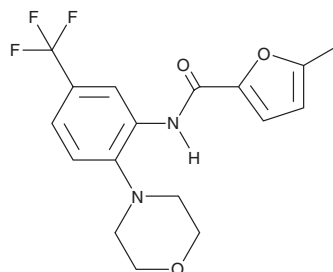
PRODUCT INFORMATION



SPHINX

Item No. 42958

CAS Registry No.: 848057-98-7
Formal Name: 5-methyl-N-[2-(4-morpholinyl)-5-(trifluoromethyl)phenyl]-2-furancarboxamide
MF: C₁₇H₁₇F₃N₂O₃
FW: 354.3
Purity: ≥98%
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

SPHINX is supplied as a solid. A stock solution may be made by dissolving the SPHINX in the solvent of choice, which should be purged with an inert gas. SPHINX is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

SPHINX is an inhibitor of serine/arginine-rich protein kinase 1 (SRPK1; IC₅₀ = 0.88 μM).¹ It is selective for SRPK1 over SRPK2 (IC₅₀ = >10 μM). SPHINX (10 μM) inhibits EGF-induced increases in phosphorylation of serine/arginine-rich splicing factor 1 (SRSF1) in ARPE-19 retinal epithelial cells. It decreases lesion area and choroidal neovascularization in mouse eyes in a model of laser-induced photocoagulation when administered at a dose of 25 ng/eye. SPHINX reduces tumor growth in a PC3 prostate cancer mouse xenograft model.²

References

1. Gammons, M.V., Fedorov, O., Ivson, D., *et al.* Topical antiangiogenic SRPK1 inhibitors reduce choroidal neovascularization in rodent models of exudative AMD. *Invest. Ophthalmol. Vis. Sci.* **54**(9), 6052-6062 (2013).
2. Mavrou, A., Brakspear, K., Hamdollah-Zadeh, M., *et al.* Serine-arginine protein kinase 1 (SRPK1) inhibition as a potential novel targeted therapeutic strategy in prostate cancer. *Oncogene* **34**(33), 4311-4319 (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, 03/07/2025

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