

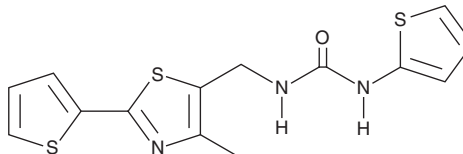
PRODUCT INFORMATION



VU0542270

Item No. 43074

CAS Registry No.: 1396814-79-1
Formal Name: N-[[4-methyl-2-(2-thienyl)-5-thiazolyl]methyl]-N'-2-thienyl-urea
MF: C₁₄H₁₃N₃OS₃
FW: 335.5
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

VU0542270 is supplied as a solid. A stock solution may be made by dissolving the VU0542270 in the solvent of choice, which should be purged with an inert gas. VU0542270 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

VU0542270 is an inhibitor of sulfonylurea receptor 2B (SUR2B) linked to ATP-sensitive potassium channel K_{ir}6.1 (SUR2B/K_{ir}6.1; IC₅₀ = 0.11 μM).¹ It is selective for SUR2B/K_{ir}6.1 over ten other inward-rectifier potassium channels (K_{ir}s; IC₅₀s = >30 μM for all). VU0542270 (10 or 100 μM) induces constriction in isolated mouse ductus arteriosus vessels.

Reference

1. Li, K., McClenahan, S.J., Han, C., *et al.* Discovery and characterization of VU0542270, the first selective inhibitor of vascular K_{ir}6.1/SUR2B KATP channels. *Mol. Pharmacol.* **105**(3), 202-212 (2024).

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, 04/21/2025

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