

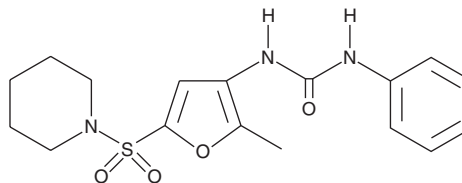
PRODUCT INFORMATION



GSK264220A

Item No. 13009

CAS Registry No.: 685506-42-7
Formal Name: N-[2-methyl-5-(1-piperidinylsulfonyl)-3-furanyl]-N'-phenyl-urea
MF: C₁₇H₂₁N₃O₄S
FW: 363.4
Purity: ≥98%
UV/Vis.: λ_{max}: 246 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

GSK264220A is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK264220A in the solvent of choice, which should be purged with an inert gas. GSK264220A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK264220A in ethanol is approximately 0.5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

GSK264220A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK264220A should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK264220A has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Endothelial lipase, a member of the triglyceride lipase gene family expressed in endothelial cells, prefers phospholipid-enriched high-density lipoprotein (HDL) over triglyceride as substrate. GSK264220A is a potent inhibitor of endothelial lipase (IC₅₀ = 16 nM).¹ Levels of HDL cholesterol are inversely correlated with endothelial lipase protein activity. Therefore, inhibition of endothelial lipase by compounds like GSK264220A may be expected to increase HDL levels and potentially decrease the risk of cardiovascular disease.

Reference

1. Keller, P.M., Rust, T., Murphy, D.J., *et al.* A high-throughput screen for endothelial lipase using HDL as substrate. *J. Biomol. Screen.* **13**(6), 468-475 (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, 01/04/2023

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