

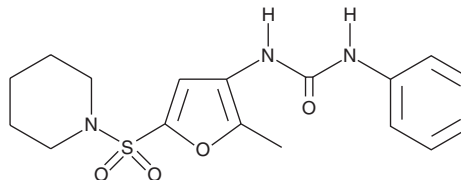
Product Information



GSK264220A

Catalog No. 13009

CAS Registry No.: 685506-42-7
Formal Name: N-[2-methyl-5-(1-piperidynylsulfonyl)-3-furanyl]-N'-phenyl-urea
MF: C₁₇H₂₁N₃O₄S
FW: 363.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that GSK264220A be stored as supplied at -20°C. It should be stable for at least two years.

GSK264220A is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK264220A in an organic solvent 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.

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. *Journal of Biomolecular Screening* **13**(6), 468-475 (2008).

Related Product

Endothelial Lipase Polyclonal Antibody - Cat. No. 100030

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 10/21/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com