

PRODUCT INFORMATION

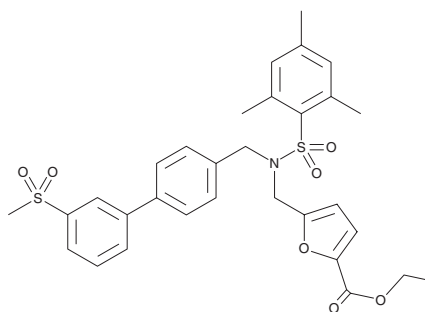


SR9238

Item No. 18771

CAS Registry No.: 1416153-62-2
Formal Name: 5-[[[3'-(methylsulfonyl)[1,1'-biphenyl]-4-yl]methyl][(2,4,6-trimethylphenyl)sulfonyl amino]methyl]-2-furancarboxylic acid, ethyl ester

MF: C₃₁H₃₃NO₇S₂
FW: 595.7
Purity: ≥98%
UV/Vis.: λ_{max}: 257 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

SR9238 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR9238 in the solvent of choice, which should be purged with an inert gas. SR9238 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SR9238 in these solvents is approximately 5 mg/ml.

SR9238 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SR9238 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SR9238 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The liver X receptors (LXRs) are nuclear receptors that act as ligand-dependent transcription factors.¹ They modulate lipid, cholesterol, and carbohydrate metabolism and homeostasis. SR9238 is an inverse agonist of the two LXR isoforms, LXRα and LXRβ (IC₅₀s = 214 and 43 nM, respectively).² It displays liver specificity *in vivo*, with little action at peripheral LXR. SR9238 suppresses hepatic lipogenesis, inflammation, lipid accumulation, and fibrosis in mouse models of non-alcoholic steatohepatitis.^{2,3} It also reduces plasma cholesterol levels in diet-induced obese mice.²

References

1. Gabbi, C., Warner, M., and Gustafsson, J.-Å. Action mechanisms of liver X receptors. *Biochem. Biophys. Res. Commun.* **446**(3), 647-650 (2013).
2. Griffett, K., Solt, L.A., El-Dien, B., *et al.* A liver-selective LXR inverse agonist that suppresses hepatic steatosis. *ACS Chem. Biol.* **8**(3), 559-567 (2013).
2. Griffett, K., Welch, R.D., Flaveny, C.A., *et al.* The LXR inverse agonist SR9238 suppresses fibrosis in a model of non-alcoholic steatohepatitis. *Mol. Metab.* **4**(4), 353-357 (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, 11/22/2022

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