

PRODUCT INFORMATION

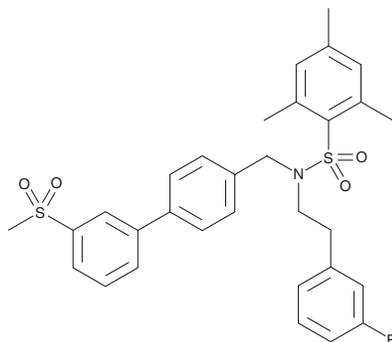


SR9243

Item No. 18420

CAS Registry No.: 1613028-81-1
Formal Name: N-[2-(3-bromophenyl)ethyl]-2,4,6-trimethyl-N-[[3'-(methylsulfonyl)[1,1'-biphenyl]-4-yl]methyl]benzenesulfonamide

MF: C₃₁H₃₂BrNO₄S₂
FW: 626.6
Purity: ≥95%
UV/Vis.: λ_{max}: 256 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

SR9243 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR9243 in the solvent of choice, which should be purged with an inert gas. SR9243 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SR9243 in these solvents is approximately 1 and 10 mg/ml, respectively.

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. SR9243 is a cell-permeable LXR inverse agonist that induces LXR-corepressor interaction at nanomolar concentrations.² It reduces cancer cell viability (IC₅₀ values range from 15 to 104 nM) without cytotoxicity against non-malignant cells.² SR9243 disrupts the Warburg effect in cancer cells, suppressing the expression of glycolytic and lipogenic genes and reducing glycolytic metabolites and lipid production. It is effective *in vivo*, blocking glycolytic and lipogenic gene expression and inducing apoptosis in colon cancer xenografts without inducing weight loss in mice.²

References

1. Gabbi, C., Warner, M., and Gustafsson, J.-Å. Action mechanisms of liver X receptors. *Biochem. Biophys. Res. Commun.* **446**(3), (2014).
2. Flaveny, C.A., Griffett, K., El-Gendy, B.E.D.M., et al. Broad anti-tumor activity of a small molecule that selectively targets the Warburg effect and lipogenesis. *Cancer Cell* **28**(1), 42-56 (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, 11/01/2022

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