

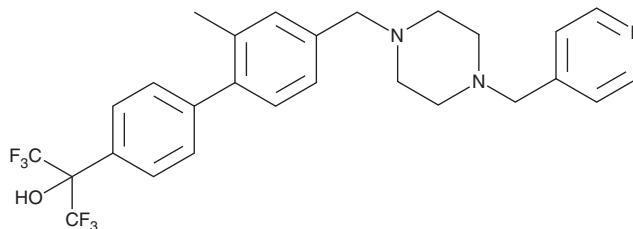
PRODUCT INFORMATION



SR 1903

Item No. 28253

CAS Registry No.: 1414248-06-8
Formal Name: 2'-methyl-4'-[[4-(4-pyridinylmethyl)-1-piperazinyl]methyl]- α,α -bis(trifluoromethyl)-[1,1'-biphenyl]-4-methanol
MF: C₂₇H₂₇F₆N₃O
FW: 523.5
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 249 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SR 1903 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 1903 in the solvent of choice, which should be purged with an inert gas. SR 1903 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SR 1903 in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

SR 1903 is a modulator of retinoic acid receptor-related orphan receptor γ (ROR γ) and liver X receptor (LXR).¹ It is an inverse agonist of ROR γ (IC₅₀ = ~100 nM in a cell-based reporter assay) and an agonist of LXR. It also binds to peroxisome proliferator-activated receptor γ (PPAR γ ; IC₅₀ = 209 nM) but does not activate it. SR 1903 (10 μ M) inhibits LPS-induced expression of triggering receptor expressed on myeloid cells 1 (TREM-1) in RAW 264.7 cells. It also inhibits LPS-induced expression of the LXR target genes IL-6 and IL-33 and increases expression of ABCG1, FASN, and SCD-1 in RAW 264.7 cells. SR 1903 (20 mg/kg twice per day) reduces severity score in a mouse model of collagen-induced arthritis. It reduces blood glucose levels in a glucose tolerance test, serum levels of total cholesterol and LDL, body weight, and fat mass in a mouse model of high-fat diet-induced obesity.

Reference

1. Chang, M.R., Ciesla, A., Strutzenberg, T.S., *et al.* Unique polypharmacology nuclear receptor modulator blocks inflammatory signaling pathways. *ACS Chem. Biol.* **14**(5), 1051-1062 (2019).

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, 12/05/2019

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