

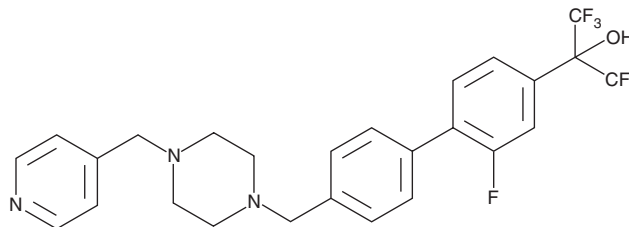
PRODUCT INFORMATION



SR 2211

Item No. 11972

CAS Registry No.: 1359164-11-6
Formal Name: 2-fluoro-4'-[[4-(4-pyridinylmethyl)-1-piperazinyl]methyl]- α,α -bis(trifluoromethyl)-[1,1'-biphenyl]-4-methanol
MF: C₂₆H₂₄F₇N₃O
FW: 527.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 251 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

SR 2211 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 2211 in the solvent of choice, which should be purged with an inert gas. SR 2211 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SR 2211 in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

Retinoic acid receptor-related nuclear receptor γ (ROR γ) plays a central role in T cell differentiation, particularly in the differentiation of pro-inflammatory T_H17 cells, which are implicated in autoimmune diseases like multiple sclerosis and rheumatoid arthritis.^{1,2} SR 2211 selectively binds ROR γ (K_i = 105 nM), acting as an inverse agonist of constitutive *in vitro* ROR γ activity (IC_{50} = 320 nM).³ It has minimal effects on ROR α , LXR α , and FXR activities. SR 2211 significantly inhibits gene expression of IL-17 and IL-23 receptor in activated EL-4 mouse T lymphocytes when given at 5 μ M.³

References

1. Ivanov, I.I., McKenzie, B.S., Zhou, L., *et al.* The orphan nuclear receptor ROR γ t directs the differentiation program of proinflammatory IL-17⁺ T helper cells. *Cell* **126(6)**, 1121-1133 (2006).
2. Solt, L.A., Kumar, N., Nuhant, P., *et al.* Suppression of T_H17 differentiation and autoimmunity by a synthetic ROR ligand. *Nature* **472(7344)**, 491-494 (2011).
3. Kumar, N., Lyda, B., Chang, M.R., *et al.* Identification of SR2211: A potent synthetic ROR γ -selective modulator. *ACS Chem. Biol.* **7(4)**, 672-677 (2012).

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, 06/21/2022

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