

PRODUCT INFORMATION



SR 2595

Item No. 18291

CAS Registry No.: 1415252-61-7

Formal Name: 4'-[[[5-[[[(1S)-1-[4-(1,1-dimethylethyl)phenyl]ethyl]amino]carbonyl]-2,3-dimethyl-1H-indol-1-yl]methyl]-[1,1'-biphenyl]-2-carboxylic acid

MF: C₃₇H₃₈N₂O₃

FW: 558.7

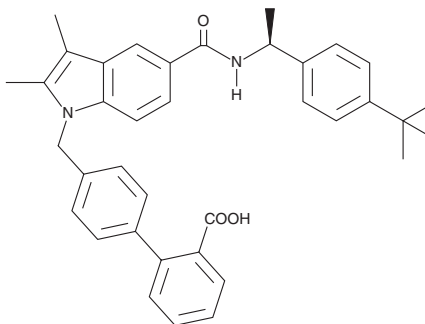
Purity: ≥98%

UV/Vis.: λ_{max}: 251 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

SR 2595 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 2595 in the solvent of choice. SR 2595 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of SR 2595 in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

SR 2595 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SR 2595 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SR 2595 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SR 2595 is an inverse agonist of PPAR γ (IC₅₀ = 30 nM) that represses both transactivation in a promoter:reporter assay and expression of the adipogenic marker fatty acid-binding protein 4 in differentiating murine preadipocytes.¹ Repression of PPAR γ with SR 2595 promotes osteogenesis, as measured by calcium phosphatase deposition, in cultured human mesenchymal stem cells (MSCs).¹ SR 2595 also increases expression of bone morphogenetic proteins BMP2 and BMP6 in MSCs.¹

Reference

1. Marciano, D.P., Kuruvilla, D.S., Boregowda, S.V., et al. Pharmacological repression of PPAR γ promotes osteogenesis. *Nature Communications* **6**, 1-7 (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, 09/23/2016

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