

# PRODUCT INFORMATION

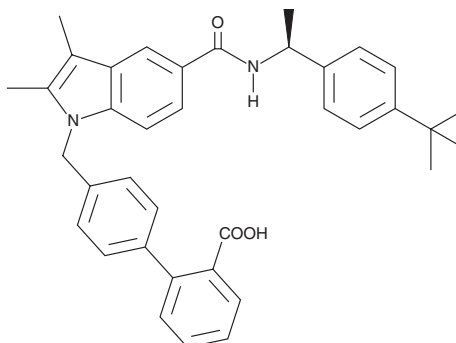


**SR 2595**

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:** C<sub>37</sub>H<sub>38</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 558.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251 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

SR 2595 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 2595 in the solvent of choice, which should be purged with an inert gas. SR 2595 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). 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 $\gamma$  (IC<sub>50</sub> = 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.<sup>1</sup> Repression of PPAR $\gamma$  with SR 2595 promotes osteogenesis, as measured by calcium phosphatase deposition, in cultured human mesenchymal stem cells (MSCs).<sup>1</sup> SR 2595 also increases expression of bone morphogenetic proteins BMP2 and BMP6 in MSCs.<sup>1</sup>

## Reference

1. Marciano, D.P., Kuruvilla, D.S., Boregowda, S.V., et al. Pharmacological repression of PPAR $\gamma$  promotes osteogenesis. *Nat. Commun.* **6**, 7443 (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, 02/05/2025

## 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](http://WWW.CAYMANCHEM.COM)