

# PRODUCT INFORMATION



**SR10221**

Item No. 18292

**CAS Registry No.:** 1415321-64-0

**Formal Name:** 2S-[2-chloro-5-[[[5-[[[(1S)-1-[4-(1,1-dimethylethyl)phenyl]ethyl]amino]carbonyl]-2,3-dimethyl-1H-indol-1-yl]methyl]phenoxy]-propanoic acid

**MF:** C<sub>33</sub>H<sub>37</sub>ClN<sub>2</sub>O<sub>4</sub>

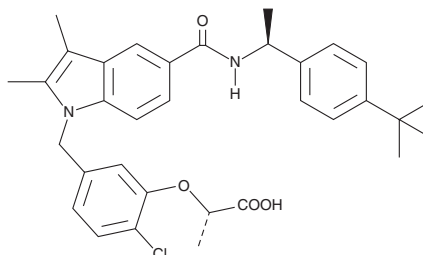
**FW:** 561.1

**Purity:** ≥98%

**Supplied as:** A 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

SR10221 is supplied as a solid. A stock solution may be made by dissolving the SR10221 in the solvent of choice, which should be purged with an inert gas. SR10221 is soluble (≥10 mg/ml) in DMSO. SR10221 is sparingly soluble (1-10 mg/ml) in ethanol.

## Description

SR10221 is an inverse agonist of peroxisome proliferator-activated receptor  $\gamma$  (PPAR $\gamma$ ).<sup>1</sup> It binds to PPAR $\gamma$  ( $K_d$  = 570 nM) and enhances PPAR $\gamma$  binding to corepressor peptides in a cell-free assay. SR10221 (1-1,000 nM) inhibits the growth of UM-UC-9 bladder cancer cells.

## Reference

1. Irwin, S., Karr, C., Furman, C., *et al.* Biochemical and structural basis for the pharmacological inhibition of nuclear hormone receptor PPAR $\gamma$  by inverse agonists. *J. Biol. Chem.* **298**(11), 102539 (2022).

### 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, 04/21/2025

## 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)