

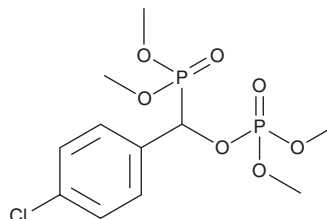
PRODUCT INFORMATION



SR 202

Item No. 21846

CAS Registry No.: 76541-72-5
Formal Name: phosphoric acid, (4-chlorophenyl)
(dimethoxyphosphinyl)methyl dimethyl ester
MF: $C_{11}H_{17}ClO_7P_2$
FW: 358.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 226 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

SR 202 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 202 in water. The solubility of SR 202 in water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

SR 202 is an antagonist of peroxisome proliferator-activated receptor γ (PPAR γ) transcriptional activity induced by troglitazone (Item No. 71750; IC_{50} = 140 μ M) but not of basal PPAR γ activity.¹ It is selective for PPAR γ , not affecting basal or agonist-induced transcriptional activity of PPAR α , PPAR β , or the farnesoid X receptor (FXR). It inhibits PPAR γ -dependent differentiation of preadipocyte 3T3-L1 cells in a dose-dependent manner. SR 202 (400 mg/kg) decreases the amount of weight gained and white adipose tissue mass accumulated by mice fed a standard or high-fat diet for ten weeks and is associated with lower PPAR γ mRNA levels. It protects against high-fat diet-induced insulin resistance in wild-type mice and improves insulin sensitivity in *ob/ob* mice.

Reference

1. Rieusset, J., Touri, F., Michalik, L., *et al.* A new selective peroxisome proliferator-activated receptor γ antagonist with antiobesity and antidiabetic activity. *Mol. Endocrinol.* **16(11)**, 2628-2644 (2002).

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, 11/02/2022

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