

PRODUCT INFORMATION



S26948

Item No. 29722

CAS Registry No.: 353280-43-0

Formal Name: 2-[[4-[2-(6-benzoyl-2-oxo-3(2H)-benzothiazolyl)ethoxy]phenyl]methyl]-propanedioic acid, 1,3-dimethyl ester

MF: C₂₈H₂₅NO₇S

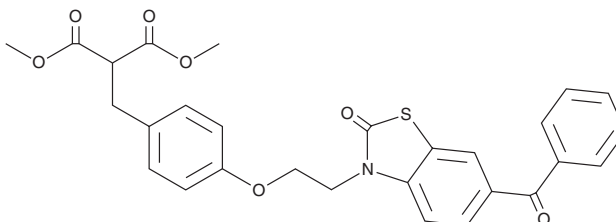
FW: 519.6

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

S26948 is supplied as a solid. A stock solution may be made by dissolving the S26948 in the solvent of choice, which should be purged with an inert gas. S26948 is soluble in the organic solvent DMSO.

Description

S26948 is a peroxisome proliferator-activated receptor γ (PPAR γ) agonist ($EC_{50} = 8.83$ nM in a transactivation assay).¹ It is selective for PPAR γ over PPAR α , PPAR β/δ , and retinoid X receptor α (RXR α) at 10 μ M. S26948 (30 mg/kg per day) decreases serum levels of glucose, triglycerides, non-esterified fatty acids, and insulin and increases hepatic palmitate oxidation in an *ob/ob* mouse model of type 2 diabetes. It decreases plasma levels of triglycerides and total and non-HDL cholesterol and the surface area of atherosclerotic lesions in the aortic sinus in an E2-KI transgenic mouse model of dyslipidemia and atherosclerosis fed a Western diet when administered at a dose of 30 mg/kg.

Reference

1. Carmona, M.C., Louche, K., Lefebvre, B., *et al.* S 26948: A new specific peroxisome proliferator-activated receptor γ modulator with potent antidiabetes and antiatherogenic effects. *Diabetes* **56(11)**, 2797-2808 (2007).

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, 11/18/2022

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