

PRODUCT INFORMATION



Saroglitazar

Item No. 27851

CAS Registry No.: 495399-09-2

Formal Name: αS-ethoxy-4-[2-[2-methyl-5-[4-(methylthio)phenyl]-1H-pyrrol-1-yl]ethoxy]-benzenepropanoic acid

MF: C₂₅H₂₉NO₄S

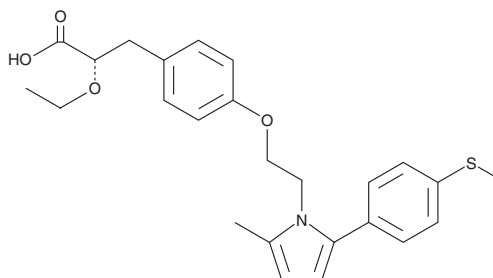
FW: 439.6

Purity: ≥90%

Supplied as: A neat oil

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Saroglitazar is supplied as a neat oil. A stock solution may be made by dissolving the saroglitazar in the solvent of choice, which should be purged with an inert gas. Saroglitazar is slightly soluble in methanol and chloroform.

Description

Saroglitazar is a free form of the peroxisome proliferator-activated receptor α (PPARα) and PPARγ dual agonist saroglitazar magnesium (Item No. 41694).

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, 09/26/2024

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