

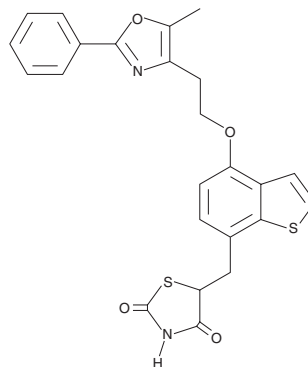
PRODUCT INFORMATION



Edaglitazone

Item No. 19117

CAS Registry No.: 213411-83-7
Formal Name: 5-[[4-[2-(5-methyl-2-phenyl-4-oxazolyl)ethoxy]benzo[b]thien-7-yl]methyl]-2,4-thiazolidinedione
Synonym: BM 13.1258
MF: C₂₄H₂₀N₂O₄S₂
FW: 464.6
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 271 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

Edaglitazone is supplied as a crystalline solid. A stock solution may be made by dissolving the edaglitazone in the solvent of choice, which should be purged with an inert gas. Edaglitazone is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

Edaglitazone is an agonist of peroxisome proliferator-activated receptor γ (PPAR γ).¹ It activates PPAR γ in a reporter assay using CV-1 cells expressing murine PPAR γ 2 when used at concentrations ranging from 0.1 to 1 μ M. Edaglitazone (1, 5, and 25 μ M) increases [³H]2-deoxyglucose transport in the presence and absence of insulin and increases glycogen synthesis in the presence of insulin in soleus muscle strips isolated from obese rats.

Reference

1. Fürnsinn, C., Brunmair, B., Meyer, M., *et al.* Chronic and acute effects of thiazolidinediones BM13.1258 and BM15.2054 on rat skeletal muscle glucose metabolism. *Br. J. Pharmacol.* **128(6)**, 1141-1148 (1999).

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, 11/09/2022

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