

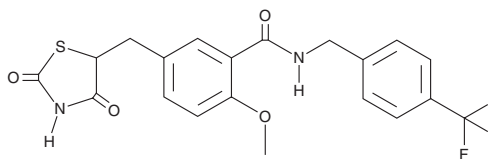
PRODUCT INFORMATION



KRP 297

Item No. 33699

CAS Registry No.: 213252-19-8
Formal Name: 5-[(2,4-dioxo-5-thiazolidinyl)methyl]-2-methoxy-N-[[4-(trifluoromethyl)phenyl]methyl]-benzamide
Synonym: MK-767
MF: C₂₀H₁₇F₃N₂O₄S
FW: 438.4
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

KRP 297 is supplied as a solid. A stock solution may be made by dissolving the KRP 297 in the solvent of choice, which should be purged with an inert gas. KRP 297 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

KRP 927 is a thiazolidinedione dual agonist of PPAR α and PPAR γ (EC₅₀s = 149 and 83 nM, respectively, in transactivation assays).¹ It decreases hyperglycemia and hyperinsulinemia in *ob/ob* mice in a dose-dependent manner. KRP 927 (10 or 30 mg/kg) reduces serum total cholesterol and triglyceride levels in hamsters.

Reference

1. Doebber, T.W., Kelly, L.J., Zhou, G., *et al.* MK-0767, a novel dual PPAR α / γ agonist, displays robust antihyperglycemic and hypolipidemic activities. *Biochem. Biophys. Res. Commun.* **318**(2), 323-328 (2004).

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, 11/04/2022

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