

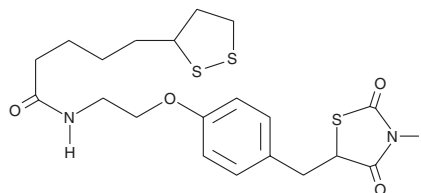
PRODUCT INFORMATION



CAY10506

Item No. 10009079

CAS Registry No.: 292615-75-9
Formal Name: N-[2-[4-[(2,4-dioxo-5-thiazolidinyl)methyl]phenoxy]ethyl]-1,2-dithiolane-3-pentanamide
MF: C₂₀H₂₆N₂O₄S₃
FW: 454.6
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

CAY10506 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of CAY10506 in ethanol is approximately 3 mg/ml and approximately 5 mg/ml in DMSO and DMF.

CAY10506 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of CAY10506 should be diluted with the aqueous buffer of choice. CAY10506 has a solubility of 0.15 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Anti-diabetic drugs of the thiazolidinedione (TZD) structural class as well as α -lipoic acid activate peroxisome proliferator-activated receptor γ (PPAR γ), a nuclear transcription factor that controls glucose metabolism and lipid homeostasis. CAY10506 is a hybrid lipoic acid-TZD derivative that transactivates human PPAR γ with an EC₅₀ value of 10 μ M.¹

Reference

1. Chittiboyina, A.G., Venkatraman, M.S., Mizuno, C.S., *et al.* Design and synthesis of the first generation of dithiolane thiazolidinedione- and phenylacetic acid-based PPAR γ agonists. *J. Med. Chem.* **49**(14), 4072-4084 (2006).

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, 11/29/2021

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