

PRODUCT INFORMATION

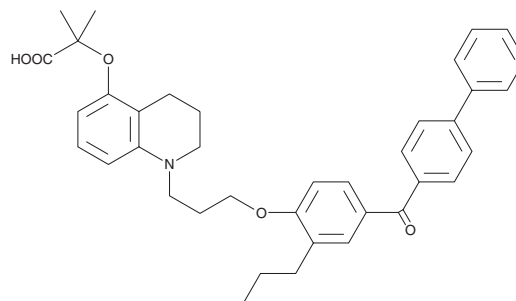


CAY10599

Item No. 13282

CAS Registry No.: 1143573-33-4
Formal Name: 2-[[1-[3-[4-([1,1'-biphenyl]-4-ylcarbonyl)-2-propylphenoxy]propyl]-1,2,3,4-tetrahydro-5-quinolinyl]oxy]-2-methyl-propanoic acid

MF: C₃₈H₄₁NO₅
FW: 591.7
Purity: ≥98%
UV/Vis.: λ_{max}: 299 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CAY10599 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10599 in the solvent of choice, which should be purged with an inert gas. CAY10599 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10599 in these solvents is approximately 25 mg/ml.

Description

CAY10599 is an agonist of peroxisome proliferator-activated receptor γ (PPAR γ ; EC₅₀ = 0.05 μ M in a transactivation assay).¹ It is selective for PPAR γ over PPAR α and PPAR δ (EC₅₀s = 3.99 and >10 μ M, respectively, in transactivation assays).

Reference

1. Lin, C.-H., Peng, Y.-H., Coumar, M.S., *et al.* Design and structural analysis of novel pharmacophores for potent and selective peroxisome proliferator-activated receptor γ agonists. *J. Med. Chem.* **52**(8), 2618-2622 (2009).

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, 09/13/2021

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