

PRODUCT INFORMATION



CAY10771

Item No. 31758

CAS Registry No.: 2522599-79-5
Formal Name: 4-[[[2-chloro-4-[[3-(2,6-dichlorophenyl)-5-(1-methylethyl)-4-isoxazolyl]methoxy]phenyl]methyl]amino]-benzeneacetic acid

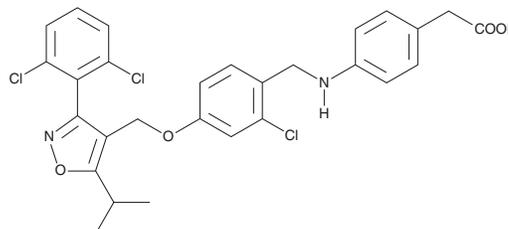
MF: C₂₈H₂₅Cl₃N₂O₄
FW: 559.9

Purity: ≥98%

Supplied as: A solution in acetonitrile

Storage: -20°C

Stability: ≥2 years



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

Description

CAY10771 is a dual agonist of farnesoid X receptor (FXR) and peroxisome proliferator-activated receptor δ (PPAR δ).¹ It activates FXR and PPAR δ in reporter assays using HEK293T cells (EC₅₀s = 0.94 and 1.5 μ M, respectively) and is selective for these receptors over retinoic acid receptor α (RAR α), retinoid X receptor α (RXR α), PPAR α , PPAR γ , and liver X receptor α (LXR α) at 10 μ M.

Reference

1. Schierle, S., Neumann, S., Heitel, P., *et al.* Design and structural optimization of dual FXR/PPAR δ activators. *J. Med. Chem.* **63**(15), 8369-8379 (2020).

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, 05/10/2024

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