

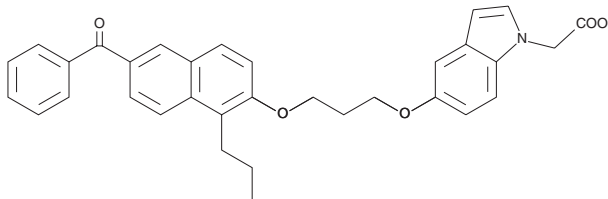
# PRODUCT INFORMATION



**CAY10573**

Item No. 10008846

**CAS Registry No.:** 853652-40-1  
**Formal Name:** 5-[3-[(6-benzoyl-1-propyl-2-naphthalenyl)oxy]propoxy]-1H-indole-1-acetic acid  
**MF:** C<sub>33</sub>H<sub>31</sub>NO<sub>5</sub>  
**FW:** 521.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 274 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

CAY10573 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10573 in an organic solvent purged with an inert gas. CAY10573 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10573 in these solvents is approximately 0.3 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

## Description

The peroxisome proliferator-activated receptors (PPARs) are lipid-activated transcription factors often used as drug targets for the treatment of metabolic disorders. CAY10573 is a PPAR agonist that displays potent binding at PPARα, γ, and δ with IC<sub>50</sub> values of 113, 50, and 223 nM, respectively. It potently transactivates PPARα, γ, and δ with EC<sub>50</sub> values of 8, 70, and 500 nM, respectively.<sup>1</sup> CAY10573 demonstrates stronger binding and functional activity for PPARγ than the antidiabetic compound rosiglitazone (IC<sub>50</sub> = 92 nM; EC<sub>50</sub> = 220 nM).<sup>1</sup>

## Reference

1. Mahindroo, N., Wang, C.-C., Liao, C.-C., *et al.* Indol-1-yl acetic acids as peroxisome proliferator-activated receptor agonists: Design, synthesis, structural biology, and molecular docking studies. *J. Med. Chem.* **49**(3), e1212-1216 (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, 01/24/2024

## 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