

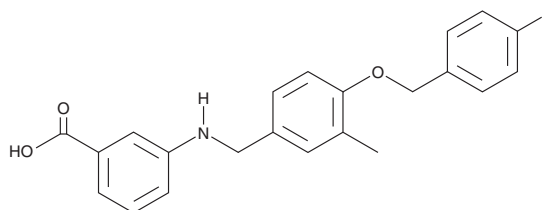
PRODUCT INFORMATION



CAY10767

Item No. 30605

CAS Registry No.: 2376590-40-6
Formal Name: 3-[[[4-[(4-fluorophenyl)methoxy]-3-methylphenyl]methyl]amino]-benzoic acid
MF: C₂₂H₂₀FNO₃
FW: 365.4
Purity: ≥98%
UV/Vis.: λ_{max}: 228 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

CAY10767 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10767 in the solvent of choice, which should be purged with an inert gas. CAY10767 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10767 in these solvents is approximately 12, 5, and 20 mg/ml, respectively.

CAY10767 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10767 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10767 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10767 is an agonist of peroxisome proliferator-activated receptor α (PPAR α ; EC₅₀ = 37 nM in a reporter assay).¹ It is greater than 2,700-fold selective for PPAR α over PPAR γ or PPAR δ .

Reference

1. Dou, X., Nath, D., Shin, H., *et al.* Evolution of a 4 benzyloxy-benzylamino chemotype to provide efficacious, potent, and isoform selective PPAR α agonists as leads for retinal disorders. *J. Med. Chem.* **63**(6), 2854-2876 (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, 07/02/2020

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