

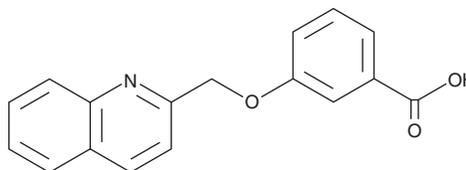
PRODUCT INFORMATION



CAY10790

Item No. 35326

CAS Registry No.: 107813-86-5
Formal Name: 3-(2-quinolinylmethoxy)-benzoic acid
MF: C₁₇H₁₃NO₃
FW: 279.3
Purity: ≥98%
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

CAY10790 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10790 in the solvent of choice, which should be purged with an inert gas. CAY10790 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10790 in these solvents is approximately 1 and 3 mg/ml, respectively.

Description

CAY10790 is a cysteinyl leukotriene 1 (CysLT₁) receptor antagonist (IC₅₀ = 2.8 μM) and G protein-coupled bile acid receptor 1 (GPBAR1) agonist.¹ It induces receptor transactivation in HEK293T cells expressing human GPBAR1 (EC₅₀ = 3 μM). CAY10790 (10 μM) reduces TNF-α and IL-1β levels in LPS-stimulated RAW 264.7 cells. It also reduces the proliferation of RAW 264.7 cells induced by leukotriene D₄ (LTD₄; Item No. 20310).

Reference

1. Fiorillo, B., Sepe, V., Conflitti, P., *et al.* Structural basis for developing multitarget compounds acting on cysteinyl leukotriene receptor 1 and G-protein-coupled bile acid receptor 1. *J. Med. Chem.* **64**(22), 16512-16529 (2021).

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, 03/26/2024

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