

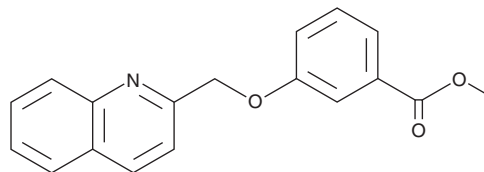
PRODUCT INFORMATION



CAY10788

Item No. 35324

CAS Registry No.: 119515-00-3
Formal Name: 3-(2-quinolinylmethoxy)-benzoic acid, methyl ester
MF: C₁₈H₁₅NO₃
FW: 293.3
Purity: ≥95%
UV/Vis.: λ_{max}: 229 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

CAY10788 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10788 in the solvent of choice, which should be purged with an inert gas. CAY10788 is soluble in organic solvents such as ethanol and DMSO. The solubility of CAY10788 in these solvents is approximately 1 mg/ml.

Description

CAY10788 is an antagonist of the cysteinyl leukotriene 1 (CysLT₁) receptor (IC₅₀ = 3.9 μM).¹

References

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