

PRODUCT INFORMATION



CAY10794

Item No. 37585

CAS Registry No.: 1231243-91-6
Formal Name: N-[4-[(2-methoxyethyl)methylamino]phenyl]-2-phenyl-5-(trifluoromethyl)-4-oxazolecarboxamide

MF: C₂₁H₂₀F₃N₃O₃

FW: 419.4

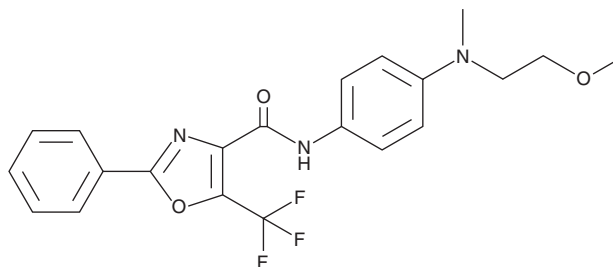
Purity: ≥98%

UV/Vis.: λ_{max}: 257 nm

Supplied as: A 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

CAY10794 is supplied as a solid. A stock solution may be made by dissolving the CAY10794 in the solvent of choice, which should be purged with an inert gas. CAY10794 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10794 in these solvents is approximately 2 and 3 mg/ml, respectively. CAY10794 is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CAY10794 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of CAY10794 in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10794 is an inhibitor of diacylglycerol acyltransferase 1 (DGAT-1; IC₅₀ = 0.325 μM).¹

Reference

1. Qian, Y., Wertheimer, S.J., Ahmad, M., *et al.* Discovery of orally active carboxylic acid derivatives of 2-phenyl-5-trifluoromethyloxazole-4-carboxamide as potent diacylglycerol acyltransferase-1 inhibitors for the potential treatment of obesity and diabetes. *J. Med. Chem.* **54**(7), 2433-2446 (2011).

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, 11/07/2022

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