

PRODUCT INFORMATION

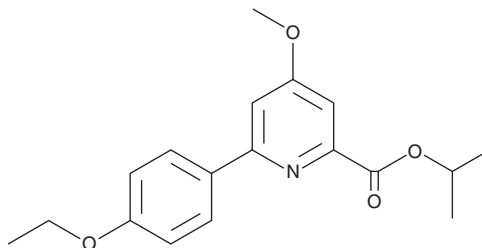


NG-497

Item No. 36886

CAS Registry No.: 2598242-66-9
Formal Name: 6-(4-ethoxyphenyl)-4-methoxy-2-pyridinecarboxylic acid, 1-methylethyl ester

MF: C₁₈H₂₁NO₄
FW: 315.4
Purity: ≥98%
UV/Vis.: λ_{max}: 256, 282 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

NG-497 is supplied as a crystalline solid. A stock solution may be made by dissolving the NG-497 in the solvent of choice, which should be purged with an inert gas. NG-497 is soluble in the organic solvent DMSO at a concentration of approximately 3 mg/ml.

Description

NG-497 is an inhibitor of adipose triglyceride lipase (ATGL; IC₅₀ = 1 μM), also known as patatin-like phospholipase domain-containing protein 2 (PNPLA2).¹ It is selective for ATGL over PNPLA1, PNPLA3, PNPLA4, and PNPLA5 at 100 μM. NG-497 selectively inhibits the triacylglycerol hydrolase activity of human and rhesus monkey ATGL over rat, dog, and marmoset ATGL, which it inhibits by less than 20% at 50 μM, and over mouse, goat, and pig ATGL at which it is inactive at 50 μM. NG-497 inhibits fatty acid and glycerol release induced by isoproterenol (Item No. 15592) in human Simpson-Golabi-Behmel syndrome (SGBS) adipocytes (IC₅₀ = 1.5 μM for both). It also inhibits forskolin-induced fatty acid release in primary human adipocytes.

Reference

1. Grabner, G.F., Gутtenberger, N., Mayer, N., *et al.* Small-molecule inhibitors targeting lipolysis in human adipocytes. *J. Am. Chem. Soc.* **144**(14), 6237-6250 (2022).

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, 10/09/2025

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