

PRODUCT INFORMATION

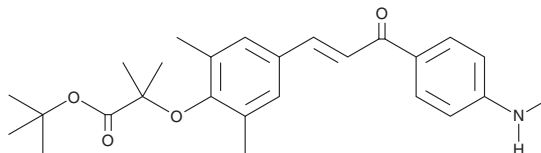


Anti-NASH Agent 1

Item No. 42614

CAS Registry No.: 2409685-41-0
Formal Name: 2-[2,6-dimethyl-4-[3-[4-(methylamino)phenyl]-3-oxo-1-propen-1-yl]phenoxy]-2-methyl-propanoic acid, 1,1-dimethylethyl ester
Synonym: Anti-Non-Alcoholic Steatohepatitis Agent 1

MF: C₂₆H₃₃NO₄
FW: 423.6
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

Anti-NASH agent 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the anti-NASH agent 1 in the solvent of choice, which should be purged with an inert gas. Anti-NASH agent 1 is soluble (≥10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Anti-NASH agent 1 is an anti-non-alcoholic steatohepatitis (NASH) agent and a derivative of the PPARα and PPARδ agonist GFT505 (elafibranor; Item No. 23508).¹ It inhibits free fatty acid-induced triglyceride accumulation and aspartate aminotransferase (AST) activity in HepG2 cells when used at a concentration of 40 μM. Anti-NASH agent 1 (3-30 mg/kg) decreases disease severity, reduces serum triglyceride, total cholesterol, LDL, Ast, alanine transaminase (Alt), TNF-α, IL-6, and IL-8 levels, and increases serum HDL and liver glutathione (GSH) levels in a mouse model of NASH induced by a methionine- and choline-deficient diet.

Reference

1. Xiang, C., Chen, X., Yao, J., et al. Design, synthesis and anti-NASH effect evaluation of novel GFT505 derivatives *in vitro* and *in vivo*. **257**, 115510 (2023).

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, 02/05/2025

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