

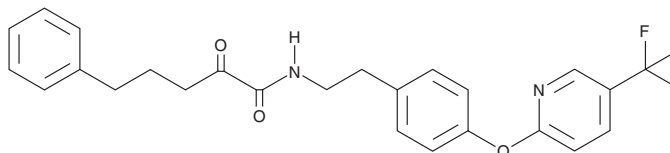
PRODUCT INFORMATION



LEI-110

Item No. 36789

CAS Registry No.: 2313525-90-3
Formal Name: α -oxo-N-[2-[4-[[5-(trifluoromethyl)-2-pyridinyl]oxy]phenyl]ethyl]-benzenepentanamide
MF: $C_{25}H_{23}F_3N_2O_3$
FW: 456.5
Purity: $\geq 98\%$
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

LEI-110 is supplied as a solid. A stock solution may be made by dissolving the LEI-110 in the solvent of choice, which should be purged with an inert gas. LEI-110 is soluble in ethanol.

Description

LEI-110 is an inhibitor of adipose phospholipase A_2 (AdPLA $_2$; $IC_{50} = 100$ nM).¹ It also inhibits HRas-like suppressor 2 (HRASL2), retinoid acid receptor responder protein 3 (RARRES3), and calcium-independent N-acyltransferase (iNAT; IC_{50} s = 158, 158, and 25 nM, respectively). LEI-110 (10 μM) decreases arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) levels in U2OS cells expressing human AdPLA $_2$ and prevents oleic acid-induced increases in lipid droplet formation in HepG2 cells.

Reference

1. Zhou, J., Mock, E.D., Martella, A., *et al.* Activity-based protein profiling identifies α -ketoamides as inhibitors for phospholipase A_2 group XVI. *ACS Chem. Biol.* **14**(2), 164-169 (2019).

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, 12/21/2022

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