

PRODUCT INFORMATION



10-oxo-12(Z)-Octadecenoic Acid ethyl ester

Item No. 37222

Synonyms: 10-keto-12Z-Octadecenoic Acid ethyl ester,
10-oxo-12-cis-Octadecenoic Acid ethyl ester

MF: C₂₀H₃₆O₃

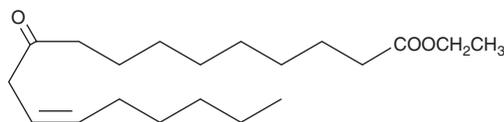
FW: 324.5

Purity: ≥95%

Supplied as: A 1 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

10-oxo-12(Z)-Octadecenoic acid ethyl ester is an ester form of 10-oxo-12(Z)-Octadecenoic acid (Item No. 36977).¹ It increases the expression of the gene encoding mitochondrial uncoupling protein 1 (*Ucp1*) in brown and white adipose tissue in a KKAY mouse model of obesity and diabetes.

Reference

1. Kim, M., Furuzono, T., Yamakuni, K., *et al.* 10-oxo-12(Z)-octadecenoic acid, a linoleic acid metabolite produced by gut lactic acid bacteria, enhances energy metabolism by activation of TRPV1. *FASEB J.* **3(11)**, 5036-5048 (2017).

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, 11/10/2025

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