

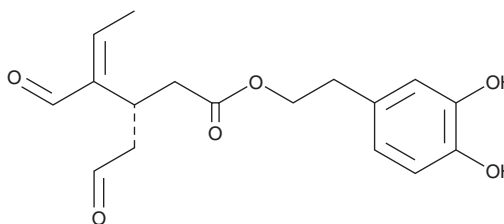
PRODUCT INFORMATION



Oleacein

Item No. 40719

CAS Registry No.: 149183-75-5
Formal Name: (3S,4E)-4-formyl-3-(2-oxoethyl)-4-hexenoic acid, 2-(3,4-dihydroxyphenyl) ethyl ester
MF: $C_{17}H_{20}O_6$
FW: 320.3
Purity: $\geq 95\%$
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Olea europaea*



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

Laboratory Procedures

Oleacein is supplied as a neat oil. A stock solution may be made by dissolving the oleacein in the solvent of choice, which should be purged with an inert gas. Oleacein is soluble in DMSO and acetonitrile.

Description

Oleacein is a secoiridoid polyphenol that has been found in *O. europaea* and has diverse biological activities.¹⁻⁴ It inhibits 5-lipoxygenase (5-LO; $IC_{50} = 2 \mu\text{M}$ for the human enzyme).¹ Oleacein (20 and 50 μM) reduces lipid accumulation and inhibits the transition into foam cells in human monocyte-derived macrophages.² It decreases lipid accumulation and levels of peroxisome proliferator-activated receptor γ (PPAR γ) and fatty acid synthase (FASN), as well as increases adiponectin levels, in 3T3-L1 adipocytes.³ *In vivo*, oleacein (20 mg/kg) reduces macrophage and lymphocyte infiltration in abdominal fat tissue and adipocyte size in a mouse model of high-fat diet-induced obesity. Oleacein also delays disease onset and progression in a mouse model of experimental autoimmune encephalomyelitis (EAE).⁴

References

1. Vougianniopoulou, K., Lemus, C., Halabalaki, M., *et al.* One-step semisynthesis of oleacein and the determination as a 5-lipoxygenase inhibitor. *J. Nat. Prod.* **77**(3), 441-445 (2014).
2. Filipek, A., Mikołajczyk, T.P., Guzik, T.J., *et al.* Oleacein and foam cell formation in human monocyte-derived macrophages: A potential strategy against early and advanced atherosclerotic lesions. *Pharmaceuticals (Basel)* **13**(4), 64 (2020).
3. Lepore, S.M., Maggisano, V., Bulotta, S., *et al.* Oleacein prevents high fat diet-induced adiposity and ameliorates some biochemical parameters of insulin sensitivity in mice. *Nutrients* **11**(8), 1829 (2019).
4. Gutiérrez-Miranda, B., Gallardo, I., Melliou, E., *et al.* Oleacein attenuates the pathogenesis of experimental autoimmune encephalomyelitis through both antioxidant and anti-inflammatory effects. *Antioxidants (Basel)* **9**(11), 1161 (2020).

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, 07/10/2024

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