

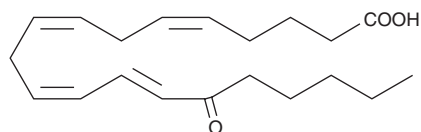
PRODUCT INFORMATION



15-OxoETE

Item No. 34730

CAS Registry No.: 81416-72-0
Formal Name: 15-oxo-5Z,8Z,11Z,13E-eicosatetraenoic acid
Synonym: 15-KETE
MF: $C_{20}H_{30}O_3$
FW: 318.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 279 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥ 2 years



Special Conditions: Oxygen and light sensitive

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

Laboratory Procedures

15-OxoETE is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. 15-OxoETE is miscible in these solvents.

15-OxoETE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 15-OxoETE should be diluted with the aqueous buffer of choice. The solubility of 15-OxoETE in PBS (pH 7.2) is approximately 0.80 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

15-OxoETE is produced by oxidation of the 15-hydroxyl of 15-HETE.¹ Whether this is a major pathway of 15-HETE metabolism is yet to be clearly established.

Reference

1. O'Flaherty, J.T., Kuroki, M., Nixon, A.B., *et al.* 5-Oxo-eicosanoids and hematopoietic cytokines cooperate in stimulating neutrophil function and the mitogen-activated protein kinase pathway. *The Journal of Biological Chemistry* **271**(30), 17821-17828 (1996).

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, 08/10/2022

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