

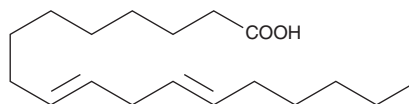
PRODUCT INFORMATION



Linoelaidic Acid

Item No. 90160

CAS Registry No.: 506-21-8
Formal Name: 9E,12E-octadecadienoic acid
Synonym: FA 18:2
MF: $C_{18}H_{32}O_2$
FW: 280.5
Purity: $\geq 98\%$
Supplied as: A solution in ethanol
Storage: -20°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

Linoelaidic acid 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of linoelaidic acid in these solvents is approximately 100 mg/ml.

Linoelaidic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of linoelaidic acid should be diluted with the aqueous buffer of choice. The solubility of linoelaidic acid in PBS (pH 7.2) is approximately 0.1 mg/ml and approximately 1 mg/ml in 0.15 M Tris-HCl (pH 8.5). We do not recommend storing the aqueous solution for more than one day.

Description

Linoelaidic acid is the all-trans fatty acid homolog of linoleic acid. Trans-fatty acids are generally not detected in raw food oils, but constitute up to 25% of the unsaturated fat in partially hydrogenated oils.¹

References

1. Ratnayake, W.M.N., Chen, Z.Y., Pelletier, G., *et al.* Occurrence of 5c,8c,11c,15t-eicosatetraenoic acid and other unusual polyunsaturated fatty acids in rats fed partially hydrogenated canola oil. *Lipids* **29**(10), 707-714 (1994).

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, 02/12/2024

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