

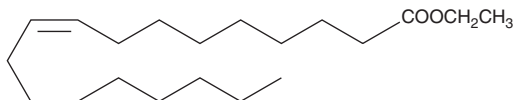
Product Information



Oleic Acid ethyl ester

Catalog No. 10008201

CAS Registry No.: 111-62-6
Formal Name: 9Z-octadecenoic acid, ethyl ester
Synonym: Ethyl Oleate
MF: C₂₀H₃₈O₂
FW: 310.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that oleic acid ethyl ester be stored as supplied at -20°C. It should be stable for at least one year.

Oleic acid ethyl ester 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 oleic acid ethyl ester in these solvents is approximately 100 mg/ml.

Oleic acid ethyl ester is sparingly soluble in aqueous buffers. Further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Oleic acid is a monounsaturated fatty acid and is one of the major components of membrane phospholipids. It contributes about 17% of the total fatty acids esterified to phosphatidylcholine in porcine platelets.¹ Oleic acid ethyl ester is a neutral, more lipid-soluble form of oleic acid. As the free acid, it inhibits collagen-stimulated platelet aggregation by approximately 90% at a concentration of 10 µg/ml.¹ It inhibits fMLF-induced neutrophil aggregation and degranulation by 55% and 68%, respectively, at 5 µM.² Oleic acid, whether applied extracellularly (EC₅₀ = ~60 µM) to human platelets or released from membrane phospholipids, causes an increase in intracellular calcium levels.³

References

1. Wahle, K.W.J. and Peacock, L.I.L. Effects of isomeric *cis* and *trans* eighteen carbon monounsaturated fatty acids on porcine platelet function. *Biochim. Biophys. Acta* **1301**, 141-149 (1996).
2. Naccache, P.H., Moiski, T.F.P., Volpi, M., *et al.* Modulation of rabbit neutrophil aggregation and degranulation by free fatty acids. *J. Leukoc. Biol.* **36**, 333-340 (1984).
3. Sifafaka-Kapadai, A., Hanahan, D.J., and Javors, M.A. Oleic acid-induced Ca²⁺ mobilization in human platelets: Is oleic acid an intracellular messenger? *J. Lipid Mediat. Cell Signal.* **15**, 215-232 (1997).

Related Products

Elaidic Acid - Cat. No. 90250 • Oleic Acid - Cat. No. 90260 • Oleic Acid-2,6-diisopropylanilide - Cat. No. 10006782 • Fatty Acid ethyl ester Standard-PAK - Cat. No. 10008188 • Stearic Acid ethyl ester - Cat. No. 10008196 • Myristic Acid ethyl ester - Cat. No. 10008197 • Linoleic Acid ethyl ester - Cat. No. 10008198 • Linolenic Acid ethyl ester - Cat. No. 10008199 • Arachidonic Acid ethyl ester - Cat. No. 10008200 • Palmitic Acid ethyl ester - Cat. No. 10008202 • Lauric Acid ethyl ester - Cat. No. 10008203 • Palmitoleic Acid ethyl ester - Cat. No. 10008204

Cayman Chemical

Mailing addressPhoneFaxE-MailWeb

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our **website and in our catalog**.

Copyright Cayman Chemical Company, 02/25/2010