

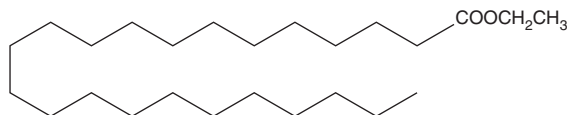
# Product Information



## Ethyl Tricosanoate

Catalog No. 10006874

**CAS Registry No.:** 18281-07-7  
**Formal Name:** tricosanoic acid, ethyl ester  
**MF:** C<sub>25</sub>H<sub>50</sub>O<sub>2</sub>  
**FW:** 382.7  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that ethyl tricosanoate be stored as supplied at -20°C. It will be stable for at least two years.

Ethyl tricosanoate is supplied as a crystalline solid. Ethyl tricosanoate is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of ethyl tricosanoate be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of ethyl tricosanoate in chloroform is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Ethyl tricosanoate, a long chain saturated fatty acid ethyl ester, is found in small quantities in liverwort, *C. conicum*.<sup>1</sup> The free acid decreases the hemolytic activity of *S. aureus* delta toxin on human erythrocytes.<sup>2</sup> Both the ethyl ester and the free acid can be used as standards for analysis of lipid mixtures.

### References

1. Matsuo, A., Nakayama, M., Hayashi, S., *et al.* Fatty acid ethyl esters in the liverwort *Conocephalum conicum*. *Phytochem.* **19**, 1848-1849 (1980).
2. Kapral, F.A. Effect of fatty acids on *Staphylococcus aureus* delta-toxin hemolytic activity. *Infect. Immun.* **13(1)**, 114-119 (1976).

### Related Products

Methyl Pentacosanoate - Cat. No. 10006451 • Palmitic Acid methyl ester - Cat. No. 10007358 • Fatty Acid ethyl ester Standard Mixture - Cat. No. 10008188 • Stearic Acid ethyl ester - Cat. No. 10008196 • Myristic Acid ethyl ester - Cat. No. 10008197 • Palmitic Acid ethyl ester - Cat. No. 10008202 • Lauric Acid ethyl ester - Cat. No. 10008203

**WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.**

#### 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 under separate cover to the MSDS supervisor at 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, 01/06/2010

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

www.caymanchem.com