

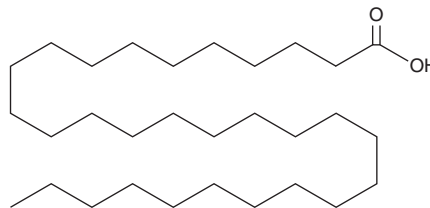
# PRODUCT INFORMATION



## Triacontanoic Acid

Item No. 34032

**CAS Registry No.:** 506-50-3  
**Synonyms:** C30:0, FA 30:0, NSC 53832, Myricic Acid, n-Triacontanoic Acid  
**MF:** C<sub>30</sub>H<sub>60</sub>O<sub>2</sub>  
**FW:** 452.8  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Triacontanoic acid is supplied as a solid. A stock solution may be made by dissolving the triacontanoic acid in the solvent of choice, which should be purged with an inert gas. Triacontanoic acid is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Triacontanoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triacontanoic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Triacontanoic acid has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Triacontanoic acid is a very long-chain saturated fatty acid. It is a component of D-003, a mixture of very-long chain aliphatic acids purified from sugar cane wax that has antiplatelet and cholesterol-lowering activities in animal models.<sup>1</sup>

### Reference

1. Gámez, R., Rodeiro, I., Fernández, I., *et al.* Preliminary evaluation of the cytotoxic and genotoxic potential of D-003: Mixture of very long chain fatty acids. *Teratog. Carcinoq. Mutagen.* **22(3)**, 175-181 (2002).

#### 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, 02/20/2024

#### 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