

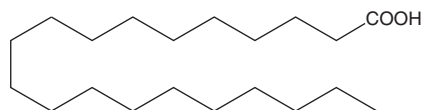
# PRODUCT INFORMATION



## Arachidic Acid

Item No. 9000339

**CAS Registry No.:** 506-30-9  
**Formal Name:** eicosanoic acid  
**Synonyms:** Arachic Acid, C20:0, *n*-Eicosanoic Acid, Icosanoic Acid, NSC 93983  
**MF:** C<sub>20</sub>H<sub>40</sub>O<sub>2</sub>  
**FW:** 312.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

Arachidic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the arachidic acid in an organic solvent purged with an inert gas. Arachidic acid is soluble in organic solvents such as ethanol and dimethyl formamide (DMF). The solubility of arachidic acid in these solvents is approximately 0.1 and 2 mg/ml, respectively.

Arachidic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, arachidic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Arachidic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Arachidic acid is a long-chain saturated fatty acid that has been found in peanut butter and anaerobic fungi.<sup>1,2</sup> It inhibits rabbit neutrophil aggregation induced by N-formyl-methionyl-leucyl-phenylalanine (fMLP; Item No. 21495) when used at a concentration of 5 μM.<sup>3</sup> Formulations containing arachidic acid have been used as surfactants in the manufacture of soaps and cosmetics.

### Reference

1. Negoita, M., Mihai, A.L., Adascalului, A., *et al.* Comparison of the fatty acid composition of peanut butter by applying different fat extraction procedures. *Rev. Chim. (Bucharest)* **69(11)**, 3023-3032 (2018).
2. Koppová, I., Novotná, Z., Štrosová, L., *et al.* Analysis of fatty acid composition of anaerobic rumen fungi. *Folia Microbiol. (Praha)* **53(3)**, 217-220 (2008).
3. Naccache, P.H., Moiski, T.F., Volpi, M., *et al.* Modulation of rabbit neutrophil aggregation and degranulation by free fatty acids. *J. Leukoc. Biol.* **36(3)**, 333-340 (1984).

#### 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, 11/04/2022

#### 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