

PRODUCT INFORMATION



Tricosanoic Acid methyl ester

Item No. 26868

CAS Registry No.: 2433-97-8

Synonyms: C23:0 methyl ester,
Methyl Tricosanoate, SFE 24:0

MF: $C_{24}H_{48}O_2$

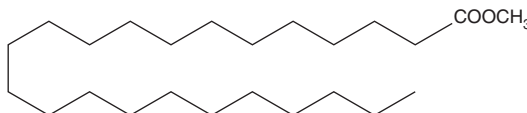
FW: 368.6

Purity: $\geq 98\%$

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

Tricosanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the tricosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Tricosanoic acid methyl ester is soluble in the organic solvent chloroform at a concentration of approximately 50 mg/ml.

Description

Tricosanoic acid methyl ester is an ester form of tricosanoic acid (Item No. 26867). It has been found in peanut butter and in latent fingerprint residue samples.^{1,2}

References

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. Michalski, S., Shaler, R., and Dorman, F.L. The evaluation of fatty acid ratios in latent fingerprints by gas chromatography/mass spectrometry (GC/MS) analysis. *J. Forensic Sci.* **58(Suppl 1)**, S215-S220 (2013).

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, 02/14/2024

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