

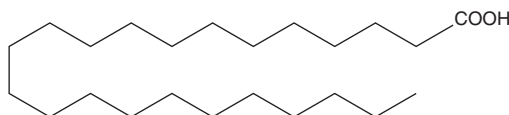
PRODUCT INFORMATION



Tricosanoic Acid

Item No. 26867

CAS Registry No.: 2433-96-7
Synonyms: C23:0 Fatty Acid, FA 23:0,
Tricosylic Acid
MF: $C_{23}H_{46}O_2$
FW: 354.6
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Tricosanoic acid is a naturally occurring saturated fatty acid that has been found in various edible mushroom species.¹ It inhibits lysis of human erythrocytes induced by *S. aureus* delta toxin in a concentration-dependent manner.² Tricosanoic acid levels are decreased in the sebum of patients with papulopustular rosacea and increased in malignant, but not adjacent nonmalignant, prostate tissue in patients with prostate cancer.^{3,4} Tricosanoic acid has been used as an internal standard for the quantification of fatty acids in chicken egg yolk.⁵

References

1. Ribeiro, B., Guedes de Pinho, P., Andrade, P.B., *et al.* Fatty acid composition of wild edible mushrooms species: A comparative study. *Microchem. J.* **93(1)**, 29-35 (2009).
2. Kapral, F.A. Effect of fatty acids on *Staphylococcus aureus* delta-toxin hemolytic activity. *Infect. Immun.* **13(1)**, 114-119 (1976).
3. Ní Raghallaigh, S. Bender, K., Lacey, N., *et al.* The fatty acid profile of the skin surface lipid layer in papulopustular rosacea. *Br. J. Dermatol.* **166(2)**, 279-287 (2012).
4. Jung, K., Reszka, R., Kamlage, B., *et al.* Tissue metabolite profiling identifies differentiating and prognostic biomarkers for prostate carcinoma. *Int. J. Cancer* **133(12)**, 2914-2924 (2013).
5. Wang, Y., Sunwoo, H., Cherian, G., *et al.* Fatty acid determination in chicken egg yolk: A comparison of different methods. *Poult. Sci.* **79(8)**, 1168-1171 (2000).

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, 03/07/2024

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