

PRODUCT INFORMATION



Hexacosanoic Acid

Item No. 13354

CAS Registry No.: 506-46-7

Synonyms: C26:0, Ceric Acid, Cerotic Acid,
FA 26:0, NSC 4205

MF: $C_{26}H_{52}O_2$

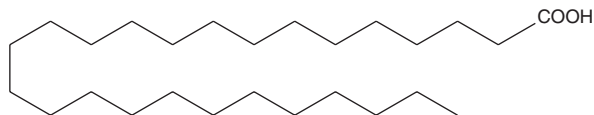
FW: 396.7

Purity: $\geq 98\%$

Supplied as: A crystalline 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

Hexacosanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the hexacosanoic acid in the solvent of choice, which should be purged with an inert gas. Hexacosanoic acid is soluble in the organic solvent chloroform at a concentration of approximately 2.5 mg/ml.

Description

Hexacosanoic acid is a 26-carbon saturated (26:0) very-long-chain fatty acid. Plasma levels of hexacosanoic acid are increased in a mouse model of adrenoleukodystrophy, a peroxisomal disorder.¹ Increased levels of hexacosanoic acid in red blood cells or in plasma positively correlated with atherosclerotic risk factors and metabolic syndrome, respectively.^{2,3}

References

1. Kobayashi, T., Shinnoh, N., Kondo, A., *et al.* Adrenoleukodystrophy protein-deficient mice represent abnormality of very long chain fatty acid metabolism. *Biochem. Biophys. Res. Commun.* **232(3)**, 631-636 (1997).
2. Antoku, Y., Tsukamoto, K., Miyoshi, Y., *et al.* Correlations of elevated levels of hexacosanoate in erythrocyte membranes with risk factors for atherosclerosis. *Atherosclerosis* **153(1)**, 169-173 (2000).
3. Kume, A., Miyazaki, T., Kitamura, Y., *et al.* High levels of saturated very long-chain fatty acid (hexacosanoic acid; C26:0) in whole blood are associated with metabolic syndrome in Japanese men. *Diab. Res. Clin. Pract.* **80(2)**, 259-264 (2008).

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, 02/07/2024

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