

PRODUCT INFORMATION



Stearic Acid ethyl ester

Item No. 10008196

CAS Registry No.: 111-61-5
Formal Name: octadecanoic acid, ethyl ester
Synonyms: Ethyl Octadecanoate, Ethyl Stearate, NSC 8919, Octadecanoic Acid ethyl ester

MF: C₂₀H₄₀O₂

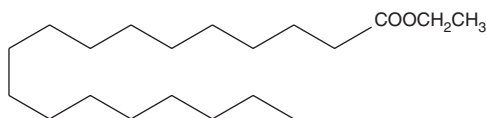
FW: 312.5

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥1 year



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

Laboratory Procedures

Stearic acid ethyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvent such as dimethyl formamide purged with an inert gas can be used. The solubility of stearic acid ethyl ester in this solvent is approximately 30 mg/ml.

Description

Stearic acid is a saturated fatty acid commonly found in animal and vegetable fats that is frequently used in cosmetics, candles, soaps, plastics, oil pastels, and for softening rubber. Stearic acid ethyl ester (ethyl stearate) is the neutral, more lipid soluble form of the free acid. It perturbs the cell cycle and induces apoptosis in HepG2 cells and is a marker of excessive alcohol consumption that can be isolated from an individual's hair.^{1,2}

References

1. Aydin, A., Celik, H.A., Deveci, R., *et al.* Induction of apoptosis by fatty acid ethyl esters in HepG2 cells. *Food Chem. Toxicol.* **43(1)**, 139-145 (2005).
2. Hartwig, S., Auwärter, V., and Pragst, F. Fatty acid ethyl esters in scalp, public, axillary, beard and body hair as markers for alcohol misuse. *Alcohol Alcohol.* **38(2)**, 163-167 (2003).

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, 02/20/2019

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