

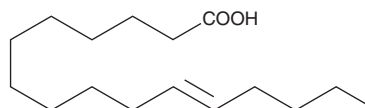
PRODUCT INFORMATION



11-Hexadecenoic Acid

Item No. 26730

CAS Registry No.: 2271-34-3
Formal Name: 11-hexadecenoic acid
Synonyms: C16:1 (*cis*-11/*trans*-11) Fatty Acid, FA 16:1
MF: C₁₆H₃₀O₂
FW: 254.4
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

11-Hexadecenoic acid is supplied as a liquid. A stock solution may be made by dissolving the 11-hexadecenoic acid in the solvent of choice which should be purged with an inert gas. 11-Hexadecenoic acid is soluble in organic solvents such as hexane, chloroform, and ethanol.

Description

11-Hexadecenoic acid is a monounsaturated fatty acid. 11-*cis*-Hexadecenoic acid and 11-*trans*-hexadecenoic acid have been found in ewe milk fat, and levels of each increase with dietary lipid supplementation of linseed, sunflower, olive, or fish oils.¹ 11-*trans*-Hexadecenoic acid has also been found in intramuscular fat samples isolated from male and female foals.² This product is a mixture of 11-*cis*- and 11-*trans*-hexadecenoic acids.

References

1. Luna, P., Rodríguez-Pino, V., and Angel de la Fuente, M. Occurrence of C16:1 isomers in milk fats from ewes fed with different dietary lipid supplements. *Food Chem.* **117(2)**, 248-253 (2009).
2. Sarriés, M.V., Murray, B.E., Troy, D., *et al.* Intramuscular and subcutaneous lipid fatty acid profile composition in male and female foals. *Meat Sci.* **72(3)**, 475-485 (2006).

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