

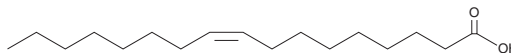
PRODUCT INFORMATION



cis-9-Heptadecenoic Acid

Item No. 36254

CAS Registry No.: 1981-50-6
Formal Name: 9Z-heptadecenoic acid
Synonyms: C17:1(9Z), FA 17:1, Margaroleic Acid
MF: $C_{17}H_{32}O_2$
FW: 268.4
Purity: $\geq 98\%$
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥ 4 years



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

Description

cis-9-Heptadecenoic acid is a monounsaturated fatty acid that is a major constituent of ruminant fat and milk.¹ The levels of *cis*-9-heptadecenoic acid in milk are negatively correlated with methane production in cows fed a high-lipid diet.²

References

1. Alves, S.P., Marcelino, C., Portugal, P.V., *et al.* Short communication: The nature of heptadecenoic acid in ruminant fats. *J. Dairy Sci.* **89**(1), 170-173 (2006).
2. Dijkstra, J., van Zijderveld, S.M., Apajalahti, J.A., *et al.* Relationships between methane production and milk fatty acid profiles in dairy cattle. *Anim. Feed Sci. Technol.* **166-167**, 23 (2011).

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, 06/17/2026

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