

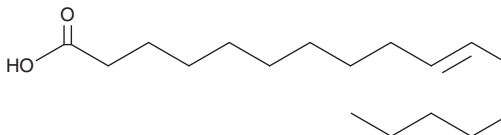
PRODUCT INFORMATION



trans-10-Heptadecenoic Acid

Item No. 22468

CAS Registry No.: 126761-43-1
Formal Name: 10E-heptadecenoic acid
Synonym: FA 17:1
MF: $C_{17}H_{32}O_2$
FW: 268.4
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 202 nm
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.

Laboratory Procedures

trans-10-Heptadecenoic acid 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. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of *trans*-10-heptadecenoic acid in these solvents is approximately 30 mg/ml.

trans-10-Heptadecenoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of *trans*-10-heptadecenoic acid should be diluted with the aqueous buffer of choice. *trans*-10-Heptadecenoic acid has a solubility of approximately 0.25 mg/ml in a 1:7 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

trans-10-Heptadecenoic acid is a C17:1 monounsaturated fatty acid that is a minor constituent of ruminant fats.¹ It is also found in the monounsaturated fatty acid pool of *C. atrox* (rattlesnakes) and remains at constant levels throughout periods of starvation.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. McCue, M.D. Western diamondback rattlesnakes demonstrate physiological and biochemical strategies for tolerating prolonged starvation. *Physiol. Biochem. Zool.* **80**(1), 25-34 (2007).

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, 02/03/2026

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