

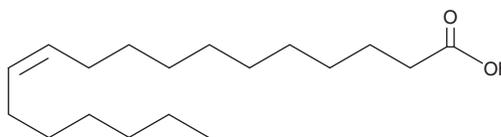
PRODUCT INFORMATION



cis-Vaccenic Acid

Item No. 20023

CAS Registry No.: 506-17-2
Formal Name: 11Z-octadecenoic acid
Synonyms: C18:1 *n*-7, C18:1(11Z),
FA 18:1, *cis*-11-Octadecenoic Acid
MF: C₁₈H₃₄O₂
FW: 282.5
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

cis-Vaccenic 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 *cis*-vaccenic acid in these solvents is approximately 30 mg/ml.

cis-Vaccenic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of *cis*-vaccenic acid should be diluted with the aqueous buffer of choice. *cis*-Vaccenic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method.

Description

cis-Vaccenic acid is an ω-7 fatty acid that has been found in mango pulp.¹ It induces differentiation of, and γ-globin synthesis in, K562 and JK-1 cells, as well as isolated sickle cell transgenic mouse bone marrow erythroid progenitor cells (TMbmEPSCs).²

References

1. Shibahara, A., Yamamoto, K., Nakayama, T., *et al.* *cis*-Vaccenic acid in mango pulp lipids. *Lipids* **21**(6), 388-394 (1986).
2. Aimola, I.A., Inuwa, H.M., Nok, A.J., *et al.* *Cis*-vaccenic acid induces differentiation and up-regulates gamma globin synthesis in K562, JK1 and transgenic mice erythroid progenitor stem cells. *Eur. J. Pharmacol.* **776**, 9-18 (2016).

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, 01/24/2024

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