

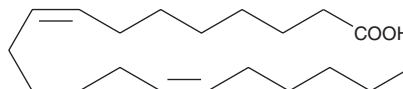
PRODUCT INFORMATION



8(Z),14(Z)-Eicosadienoic Acid

Item No. 16854

CAS Registry No.: 135498-07-6
Formal Name: (8Z,14Z)-eicosadienoic acid
Synonym: FA 20:2
MF: $C_{20}H_{36}O_2$
FW: 308.5
Purity: $\geq 95\%$
Supplied as: A solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

8(Z),14(Z)-Eicosadienoic 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 8(Z),14(Z)-eicosadienoic acid in these solvents is approximately 100 mg/ml.

8(Z),14(Z)-Eicosadienoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 8(Z),14(Z)-eicosadienoic acid should be diluted with the aqueous buffer of choice. The solubility of 8(Z),14(Z)-eicosadienoic acid in PBS (pH 7.2) and 0.15 M Tris-HCl (pH 8.5) is approximately 0.1 and 1 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

8(Z),14(Z)-Eicosadienoic Acid is an ω -8 $C_{20}:2$ fatty acid. The presence of 8(Z),14(Z)-eicosadienoic acid has been detected in human milk at a level of 0.19% (weight % total fatty acids).¹ Eicosadienoic acids are converted by desaturases, *in vivo*, to eicosatrienoic acids, which are potent vasodilators. The physiological effects of 8(Z),14(Z)-eicosadienoic acid are unstudied.

Reference

1. Ratnayake, W.M.N. and Chen, Z.-Y. *Trans*, n-3, and n-6 fatty acids in canadian human milk. *Lipids* **31(1)**, S279-S282 (1996).

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, 02/08/2024

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