

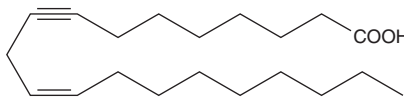
PRODUCT INFORMATION



8,11-Eicosadiynoic Acid

Item No. 90100

CAS Registry No.: 82073-91-4
Formal Name: 8,11-Eicosadiynoic acid
Synonyms: EDYA, FA 20:4
MF: $C_{20}H_{32}O_2$
FW: 304.5
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 1 year



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

Laboratory Procedures

8,11-eicosadiynoic acid (EDYA) is supplied as a crystalline solid. A stock solution may be made by dissolving the EDYA in the solvent of choice, which should be purged with an inert gas. EDYA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of EDYA in these solvents is approximately at least 100 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of EDYA can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of EDYA in 0.15 M Tris-HCl (pH 8.5) is at least 1 mg/ml, whereas the solubility in PBS (pH 7.2) is less than 100 $\mu\text{g}/\text{ml}$. For greater aqueous solubility, EDYA can be directly dissolved in 0.1 M Na_2CO_3 and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

Description

EDYA is an acetylenic fatty acid which inhibits arachidonic acid metabolism at several unrelated enzymatic stages. EDYA inhibits arachidonate uptake by platelets, arachidonoyl synthase, acyl-CoA synthase, $\Delta 5$ -desaturase, and COX.¹

Reference

1. Laposata, M., Kaiser, S.L., Reich, E.L., *et al.* Eicosadiynoic acid: A non-toxic inhibitor of multiple enzymatic steps in the production of eicosanoids from arachidonic acid. *Prostaglandins* **33**, 603-614 (1987).

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, 02/20/2024

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