

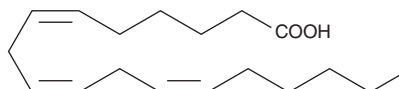
PRODUCT INFORMATION



γ -Linolenic Acid

Item No. 90220

CAS Registry No.: 506-26-3
Formal Name: 6Z,9Z,12Z-octadecatrienoic acid
Synonyms: FA 18:3, GLA
MF: C₁₈H₃₀O₂
FW: 278.4
Purity: \geq 98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

GLA 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 or dimethyl formamide purged with an inert gas can be used. The solubility of GLA in these solvents is approximately 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. If an organic solvent-free solution of GLA is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of GLA in 0.15 M Tris-HCl (pH 8.5) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GLA is an ω -6 fatty acid which can be elongated to arachidonic acid for endogenous eicosanoid synthesis. It is a weak LTB₄ receptor antagonist, inhibiting [³H]-LTB₄ binding to porcine neutrophil membranes with a K_i of 1 μ M. GLA produces 53% inhibition at a 1 mg/kg dose in an *in vivo* model of LTB₄-induced bronchoconstriction.¹

Reference

1. Yagaloff, K.A., Franco, L., Simko, B., *et al.* Essential fatty acids are antagonists of the leukotriene B₄ receptor. *Prostaglandins Leukot. Essent. Fatty Acids* **52(5)**, 293-297 (1995).

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, 02/12/2024

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