

PRODUCT INFORMATION

ALA and GLA Oxylin MaxSpec® LC-MS Mixture

Item No. 22638

Synonym: α-Linolenic Acid and γ-Linolenic Acid Oxylin LC-MS Mixture
Supplied as: A solution in ethanol (1 µg/ml of each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥6 years

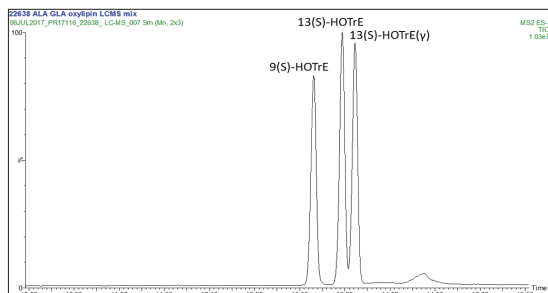
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

The ALA and GLA oxylin MaxSpec® LC-MS mixture contains oxylin metabolites derived from α-linolenic acid (ALA; Item No. 90210) and γ-linolenic acid (GLA; Item No. 90220). The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in LC-MS applications. The solution may be used as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

This mixture contains 9(S)-HOTrE (Item No. 39420), 13(S)-HOTrE (Item No. 39620), and 13(S)-HOTrE(γ) (Item No. 39610).

Contents



Item Number: 22638		ALA and GLA Oxylin MaxSpec® LC-MS Mixture			
Item Number	Item Name	Formula	Mass (Da)	[MH] ⁺ (m/z)	RT (min)
39420	9(S)-HOTrE	C ₁₈ H ₃₀ O ₃	294.4	293.3	13.16
39620	13(S)-HOTrE	C ₁₈ H ₃₀ O ₃	294.4	293.3	13.45
39610	13(S)-HOTrE(γ)	C ₁₈ H ₃₀ O ₃	294.4	293.3	13.62
LC-MS Conditions: Waters Acquity UPLC-Xevo TQD					
Mobile Phase A: Water + 0.1% Formic Acid					
Mobile Phase B: Acetonitrile + 0.1% Formic Acid					
Column: Waters BEH C8, 2.1 x 100 mm, 1.7 µm			Flow Rate: 400 µl/min		
Negative Electrospray Ionization			Selective Ion Monitoring		

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, 03/23/2026

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