

PRODUCT INFORMATION



Deuterated Saturated/Monounsaturated Fatty Acid MaxSpec® LC-MS Mixture

Item No. 30652

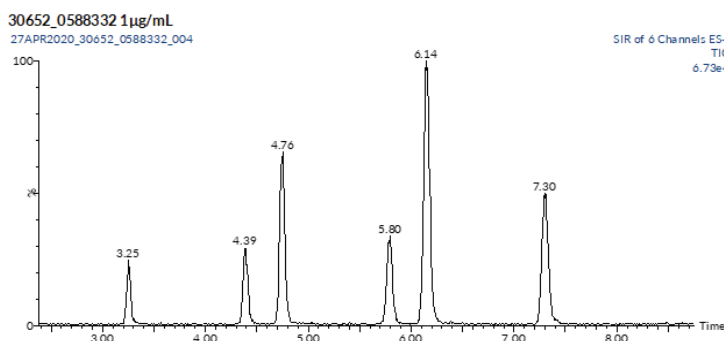
Supplied as: A solution in ethanol (10 µg/ml of each compound)
Fill volume: >1 ml
Storage: -20°C
Stability: ≥5 years

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

Description and Contents

The Deuterated Saturated/Monounsaturated Fatty Acid MaxSpec® LC-MS Mixture contains deuterated forms of select saturated or monounsaturated fatty acids. 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. This mixture is ideally suited for use as an internal standard in mass spectrometry applications. 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 the deuterium-labeled saturated fatty acids lauric acid-d₂₃ (Item No. 28080), myristic acid-d₂₇ (Item No. 9003317), palmitic acid-d₃₁ (Item No. 16497), stearic acid-d₃₅ (Item No. 9003318) and the deuterium-labeled monounsaturated fatty acids oleic acid-d₁₇ (Item No. 9000432) and palmitoleic acid-d₁₄ (Item No. 9000431).



Item Number: 30652		Item Name: Deuterated Saturated/Monounsaturated Fatty Acid MaxSpec® LC-MS Mixture		
Item Number	Item Name	Formula Weight:	MS/MS Transition:	RT (min)
28080	Lauric Acid-d ₂₃	223.5	222.3	3.25
9003317	Myristic Acid-d ₂₇	255.5	254.4	4.39
9000431	Palmitoleic Acid-d ₁₄	268.4	267.3	4.76
16497	Palmitic Acid-d ₃₁	287.6	286.4	5.80
9000432	Oleic Acid-d ₁₇	299.6	298.4	6.14
9003318	Stearic Acid-d ₃₅	319.7	318.5	7.30
LC-MS Conditions: Waters Acquity UPLC – Xevo TQD				
Mobile Phase A: 60/40/0.1 ACN/H ₂ O/Formic Acid + 10 mM Ammonium Formate				
Mobile Phase B: 10/90/0.1 ACN/IPA/Formic Acid + 10 mM Ammonium Formate				
Column: Waters BEH C8, 2.1x100 mm, 1.7 µm, 30°C			Flow Rate: 250 µl/min	
LC Gradient: Isocratic 25%B to 50%B over 7 minutes				
Negative Electrospray Ionization		SIR Scan		

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