

CERTIFICATE of ANALYSIS



Deuterated Primary Prostaglandin Metabolite MaxSpec® LC-MS Mixture

Item No.: 19304
Batch No.: 0820750

Expiry Date: 13MAY2031 (valid from date of certification)
Supplied as: A solution in ethanol (1 µg/ml of each compound; nominal); in amber ampules

Contents: See table
Volume per Ampule: Not less than 1 ml. Ampules are overfilled. It is advised that labs use measured volumes.

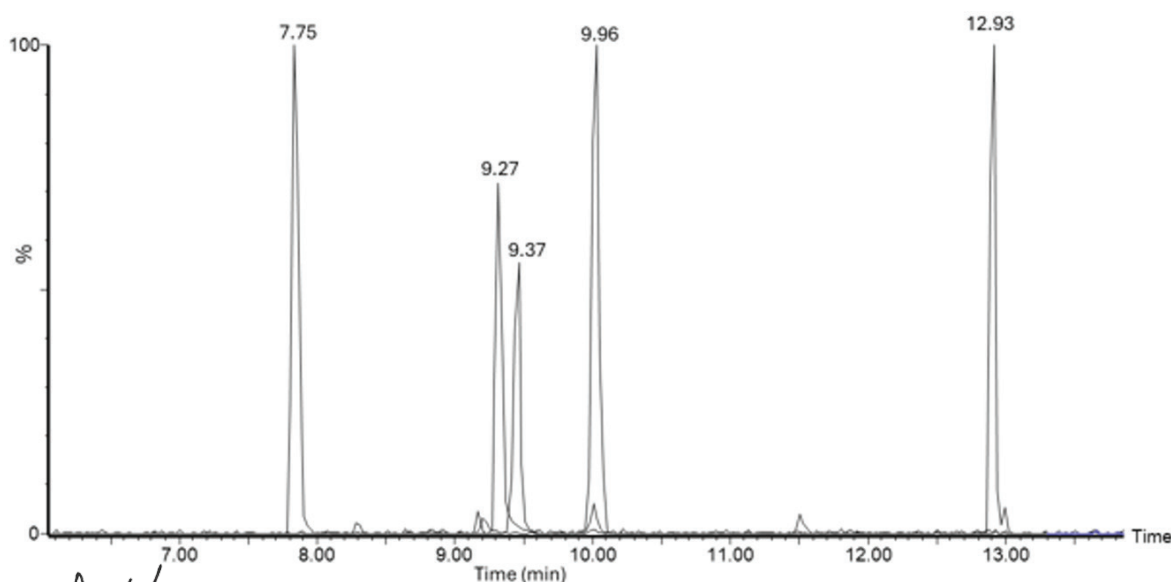
Storage & Handling: Store upright and unopened at -20°C. Warm to room temperature prior to opening.

Description

MaxSpec® LC-MS mixtures are purpose-built for qualitative mass spectrometry applications. This solution has been prepared using gravimetric and volumetric preparation procedures from a combination of neat and formulated pure standards, and is supplied in ampules sealed under argon. Ongoing stability tests are performed to ensure the presence of each analyte throughout the shelf life of the product. Follow recommended storage and handling conditions.

Quality Information

Test	Limit	Result	Meets Specification
Negative ESI-HRMS	Identities confirmed	Pass	✓



Approval:

Title: Manager, Analytical Chemistry

Certification Date: 13MAY2026

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiry date when stored unopened as recommended.

CERTIFICATE of ANALYSIS



Quality Control Table

Item No.	Item Name	Formula	RT (min)	Formula Weight	MS/MS Transition
10008989	11 β -Prostaglandin F _{2α} -d ₄	C ₂₀ H ₃₀ D ₄ O ₅	7.75	358.5	357>197
10007793	13,14-dihydro-15-keto Prostaglandin F _{2α} -d ₄	C ₂₀ H ₃₀ D ₄ O ₅	9.27	358.5	357>187
10010606	13,14-dihydro-15-keto Prostaglandin E ₂ -d ₄	C ₂₀ H ₂₈ D ₄ O ₅	9.37	356.5	355>113
10007978	13,14-dihydro-15-keto Prostaglandin D ₂ -d ₄	C ₂₀ H ₂₈ D ₄ O ₅	9.96	356.5	355>211
318570	15-deoxy- Δ 12,14-Prostaglandin J ₂ -d ₄	C ₂₀ H ₂₄ D ₄ O ₃	12.93	320.5	319>203
HRAM LC-MS/MS Conditions: Ultimate 3,000 UHPLC&Q-Exactive Triple Quadrupole-Orbitrap (Thermo Scientific)					
Mobile Phase A: Water with 0.1% Formic Acid				Time (min)	%A
Mobile Phase B: Acetonitrile with 0.1% Formic Acid				0.00	95
Column: Waters BEH C18 2.1 x 100 mm, 1.7 μ m				4.00	75
Flow Rate: 400 μ L/min				12.00	50
Negative Electrospray Ionization				12.01	2
SIR Scan				13.00	2
				13.01	95

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	13MAY2026	Initial version

Disclaimers

Material Safety Data

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some but not all of the information required for the safe and proper use of this material. Before use, review the complete Material Safety Data Sheet, which has been sent via email to your institution.

Warranty and Limitation of Remedy

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specification.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver of Buyer of all claims hereunder with respect to said material. For further details, please refer to our Warranty and Limitations of Remedy located on our website and in our catalog.

This Certificate shall not be reproduced except in full, without written approval from the Cayman Chemical Company.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com