

CERTIFICATE of ANALYSIS



Cannabis Flavonoids Mixture

Item No.: 37832
Batch No.: 0671650

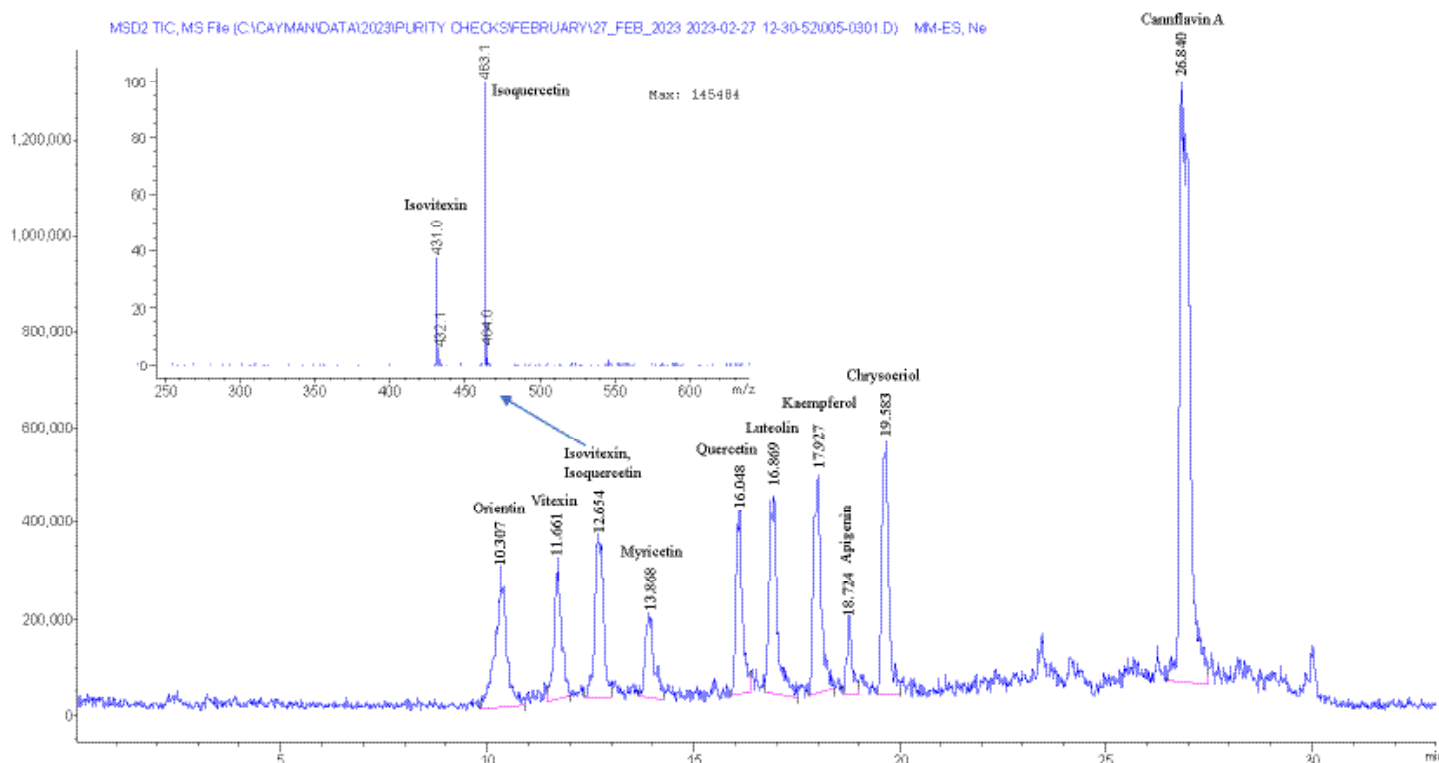
Expiry Date: 15MAR2027 (valid from date of certification)
Supplied as: Each analyte at 100 µg/ml solution in N,N-dimethylacetamide; in a deactivated glass ampule

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

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

Quality Information

Test	Specification	Result
Analyte Presence, LC-MS	All Analytes Identified	Pass



Approval:

Title: Manager, Analytical Chemistry

Certification Date: 08MAR2023

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	Exact Mass	RT (min)
25143	Orientin	448.1	10.307
15116	Vitexin	432.1	11.661
31212	Isovitexin	432.1	12.654
24926	Isoquercetin	464.1	12.654
10012600	Myricetin	318.0	13.868
10005169	Quercetin (hydrate)	302.0	16.048
10004161	Luteolin	286.1	16.869
11852	Kaempferol	286.1	17.927
10010275	Apigenin	270.1	18.724
33395	Chrysoeriol	300.1	19.583
21970	Cannflavin A	436.2	26.840
LC-MS Conditions: Agilent HPLC-Quadrupole MS			
Mobile Phase A: 90:10:0.1 Water:Methanol:Acetic Acid			
Mobile Phase B: 10:90:0.1 Water:Methanol:Acetic Acid			
Column: Kinetex Biphenyl, 250 x 4.6 mm, 5 µm, Ambient		Flow Rate: 1 ml/min	
Gradient: 25% B to 100% B over 20 minutes			
Negative APCI/ESI Ionization		TIC Scan	

Mixture was prepared using gravimetric solution preparation procedures from neat analytical standards. The concentration of this mixture was prepared with an estimated error of $\pm 10\%$.

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	08MAR2023	Initial version
02	03MAR2025	Expiry date extension

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