

CERTIFICATE of ANALYSIS



ACCREDITED
ISO 17034 #AR-1774

Kava Component Mixture 9 (CRM)

Certified Reference Material

Item No.:	41260
Batch No.:	0803531
Expiry Date:	10MAR2027 (valid from date of certification)
Supplied as:	A 250 µg/ml (nominal) multi-component solution in acetonitrile
Volume per Ampule:	Not less than 1 ml. Ampules are overfilled.
Storage:	Unopened at -20°C ± 10°C.
Safety:	Refer to Safety Data Sheet
Intended Use:	For analytical testing purposes only, not intended for human or animal use.
Instructions for Use:	This product is designated for one-time use and should be used immediately after opening. It is advised that laboratories warm the vial to room temperature prior to opening and use measured volumes.

Certified Concentration

Compound	Corrected Purity*	Concentration
Desmethoxyyangonin	98.34%	250.0 µg/ml ± 2.5 µg/ml
Dihydrokavain	99.70%	250.0 µg/ml ± 3.0 µg/ml
Dihydromethysticin	99.26%	250.1 µg/ml ± 3.3 µg/ml
(±)-Kavain	99.44%	250.0 µg/ml ± 3.3 µg/ml
Methysticin	97.55%	250.2 µg/ml ± 3.3 µg/ml
Yangonin	98.57%	250.1 µg/ml ± 3.0 µg/ml
Flavokawain A	98.52%	249.9 µg/ml ± 3.3 µg/ml
Flavokawain B	98.82%	250.0 µg/ml ± 3.3 µg/ml
Flavokawain C	99.00%	249.9 µg/ml ± 1.3 µg/ml

Concentration is calculated based on product mass, solution mass, corrected purity, and density at 20°C. It is traceable to SI units through an unbroken chain of measurements. Uncertainty of concentration is expressed as an expanded uncertainty in accordance with ISO standards for Testing Laboratories and Reference Material Producers at the approximate 95% confidence interval using a coverage factor of k=2 and incorporates uncertainties from the corrected purity, solution preparation, homogeneity, and long- and short-term stability. Concentration was verified by comparison to an independently prepared calibration standard.

*Corrected purity is determined as follows:

Corrected Purity = [(100 - % LOD - % ROI)*Chromatographic Purity/100] or [(100 - % KF - % RS - % ROI)*Chromatographic Purity/100].

Where applicable, optical rotation, chiral purity, and/or isotopic purity testing are performed to support the identification of the reference material, therefore the uncertainty is considered null.

Approval: 
Roxanne Franckowski

Title: Senior Manager of ISO Quality

Certification Date: 10MAR2025

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



CERTIFICATE of ANALYSIS



CRM Assay

Method Parameters													
Cayman Method	F0594												
Column	3.0 x 75 mm, 2.2 µm ShimPack ODS II												
Mobile Phase	A: Water B: Isopropanol												
Gradient	<table><tr><td>Time (min)</td><td>%B</td></tr><tr><td>0-2</td><td>20%</td></tr><tr><td>2-5.5</td><td>20-27.5%</td></tr><tr><td>5.5-8</td><td>27.5-72.5%</td></tr><tr><td>8-10</td><td>72.5%</td></tr><tr><td>10.1-12</td><td>20%</td></tr></table>	Time (min)	%B	0-2	20%	2-5.5	20-27.5%	5.5-8	27.5-72.5%	8-10	72.5%	10.1-12	20%
Time (min)	%B												
0-2	20%												
2-5.5	20-27.5%												
5.5-8	27.5-72.5%												
8-10	72.5%												
10.1-12	20%												
Flow Rate	0.7 ml/min												
Column Temp	50°C												
Wavelength	UV monitored at 245 nm												

Neat Material Quality Information

Qualifier	CAS No.	Item No.	Batch No.
Desmethoxyyangonin	15345-89-8	40313	0701292
Dihydrokavain	587-63-3	40406	0707850
Dihydromethysticin	19902-91-1	39944	0700696
(±)-Kavain	3155-48-4	39943	0701886
Methysticin	495-85-2	40311	0702568
Yangonin	500-62-9	40312	0700458
Flavokawain A	37951-13-6	40316	0702575
Flavokawain B	1775-97-9	40416	0702665
Flavokawain C	37308-75-1	40317	0702666

Homogeneity

A minimum sample size of 0.3 µg was used to determine homogeneity. Homogeneity was determined by HPLC using ampules selected from a random sampling plan from the early, middle, and late fill positions.

Qualifier	%RSD Result	Acceptance Criteria
Desmethoxyyangonin	1.02%	≤3%
Dihydrokavain	0.85%	≤3%
Dihydromethysticin	0.83%	≤3%
(±)-Kavain	0.83%	≤3%
Methysticin	0.84%	≤3%
Yangonin	0.85%	≤3%
Flavokawain A	0.83%	≤3%
Flavokawain B	0.89%	≤3%
Flavokawain C	1.23%	≤3%

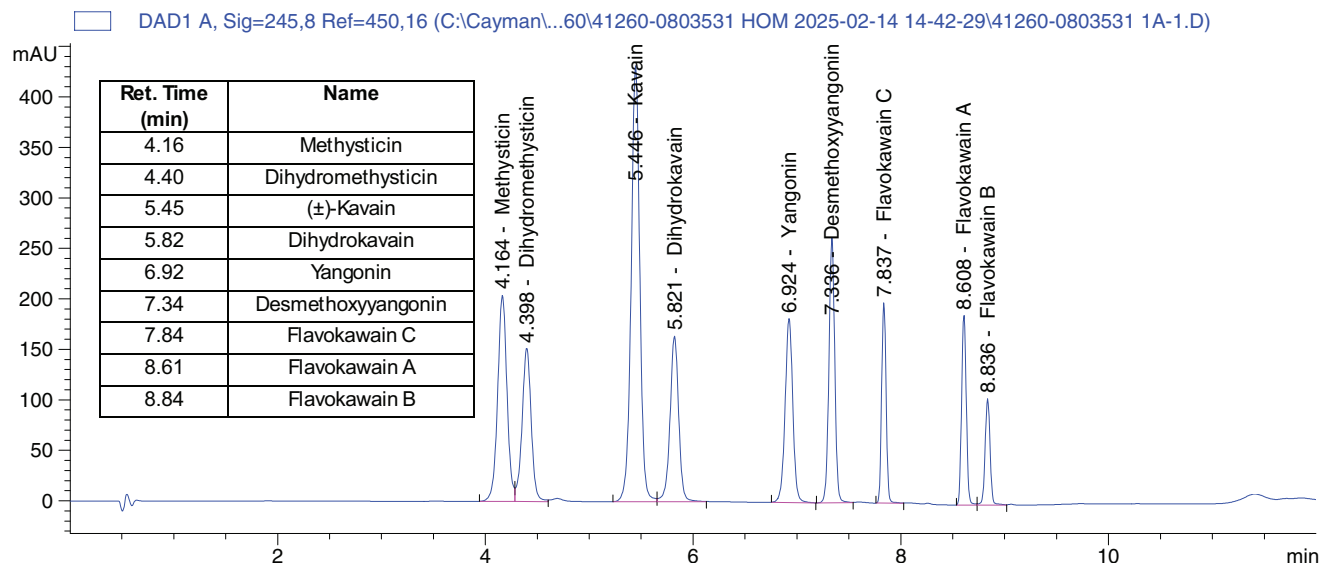
The recommended minimum quantity for use is 0.3 µg. Quantities below this have not been evaluated.

CERTIFICATE of ANALYSIS



Supplemental Data

HPLC-UV



Stability

The effect of the components of stability on the combined standard uncertainty of the CRM property value are considered negligible unless indicated in stability studies.

Short-Term Stability

Degradation was observed at 4°C and room temperature after 2 weeks. This data supports shipping of this product on dry ice.

Long-Term Stability

Long-term stability data predicts two years stability at the -20°C storage temperature. Long-term stability studies are ongoing and the Certificate of Analysis will be updated upon study completion.

CERTIFICATE of ANALYSIS



Quality Standard Documentation

The manufacturer of this Certified Reference Material is accredited under ISO 17034:2016 accreditation issued by ANAB. Refer to ANAB certificate and scope of accreditation AR-1774.

The manufacturer of this Certified Reference Material is accredited under ISO/IEC 17025:2017 accreditation issued by ANAB. Refer to the ANAB certificate and scope of accreditation AT-1773.

Revision History

Revision No.	Date	Reason for Revision
01	10MAR2025	Initial version
02	12FEB2026	Updated expiry date and certified concentration

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 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 ISO Quality Manager.

ISO CRT SD03 V. 3.1

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

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

crmquality@caymanchem.com
www.caymanchem.com