

PRODUCT INFORMATION



Kava Lactones Mixture (CRM)

Item No. 39410

Supplied as: A solution in acetonitrile (1 mg/ml of each compound)

Fill volume: >1 ml

Storage: -20°C

Stability: ≥1 year

Special conditions: Vortex or sonicate before use

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

Description and Data

Kava Lactones Mixture (CRM) (Item No. 39410) is a certified reference material mixture containing kavalactones. The mixture contains desmethoxyyangonin, dihydrokavain, dihydromethysticin, (+)-kavain, methysticin, and yangonin. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and as a reference standard for these substances during routine analytical checks using HPLC-UV and LC-MS instrumentation. This product is intended for research and forensic applications.

This mixture contains 6 kavalactones. Included are:

Item No. 39410	Kava Lactones Mixture (CRM)	
Compound	CAS Number	Concentration (mg/ml)
Desmethoxyyangonin	15345-89-8	1.0
Dihydrokavain	587-63-3	1.0
Dihydromethysticin	19902-91-1	1.0
(±)-Kavain	3155-48-4	1.0
Methysticin	495-85-2	1.0
Yangonin	500-62-9	1.0

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, 06/25/2025

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