

PRODUCT INFORMATION



M. speciosa (Kratom) Alkaloid Mixture (CRM)

Item No. 39012

Synonym: *Mitragyna speciosa* (Kratom) Alkaloid Mixture

Supplied as: A solution of 0.1 N ammonia in methanol

Fill volume: >1 ml

Storage: -20°C

Stability: ≥1 year

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

Description and Data

M. speciosa (Kratom) Alkaloid Mixture (CRM) (Item No. 39012) is a certified reference material mixture containing alkaloids found in *M. speciosa* (Kratom in Thai). The mixture contains the opioids mitragynine and 7-hydroxy mitragynine, as well as paynantheine, speciogynine, and speciociliatine, which are structurally similar to known opioids. 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 five alkaloids found in *M. speciosa* (Kratom in Thai). Included are:

Item No. 39012	<i>M. speciosa</i> (Kratom) Alkaloid Mixture (CRM)	
Item Name	CAS Number	Concentration (µg/ml)
Mitragynine	4098-40-2	500
7-hydroxy Mitragynine	174418-82-7	100
Paynantheine	4697-66-9	500
Speciogynine	4697-67-0	500
Speciociliatine	14382-79-7	500

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, 01/16/2024

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