

PRODUCT INFORMATION



Kratom Alkaloids Mixture

Item No. 37649

Supplied as: A solution in 0.1 N ammonia in methanol (250 µg/ml of each compound)

Fill Volume: >1 ml

Storage: -20°C

Stability: ≥4 years

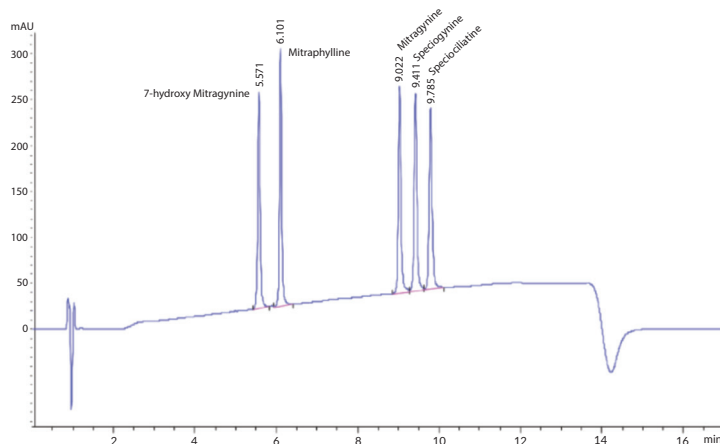
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Kratom alkaloids mixture (Item No. 37649) is a mixture of analytical reference standards containing alkaloids found in *M. speciosa* (Kratom in Thai).¹ The mixture contains the opioids mitragynine and 7-hydroxy mitragynine, as well as mitraphylline, 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 methods development and as a reference standard for these substances during routine analytical checks using HPLC or LC-MS instrumentation. This product is intended for research and forensic applications.

This mixture contains alkaloids found in *M. speciosa* (Kratom). Included are:

Item Number	Item Name	Concentration
11151	Mitragynine	250 µg/ml
13114	7-hydroxy Mitragynine	250 µg/ml
25794	Speciogynine	250 µg/ml
27246	Speciociliatine	250 µg/ml
25528	Mitraphylline	250 µg/ml



Item No.	Item Name	CAS Number	Concentration (µg/ml)
11151	Mitragynine	4098-40-2	250
13114	7-hydroxy Mitragynine	174418-82-7	250
25794	Speciogynine	4697-67-0	250
27246	Speciociliatine	14382-79-7	250
25528	Mitraphylline	509-80-8	250
HPLC Conditions			
Mobile Phase A: 90:100.1 Water:Acetonitrile:TFA			
Mobile Phase B: 10:900.1 Water:Acetonitrile:TFA			
Column: Kinetix Evo C18 2.7 µm, 4.6 x 100 mm		Flow Rate: 1.2 ml/min	
Wavelength: 210 nm			
Gradient: 0-10 min 0-40% B; 10-12 min hold 40% B; 12.1 min 40-0% B			

Reference

1. Basiliere, S. and Kerrigan, S. Identification of metabolites and potential biomarkers of kratom in urine. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **1140**, 121971 (2020).

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