

PRODUCT INFORMATION



Kanna Alkaloids Mixture

Item No. 45074

Synonym: *S. tortuosum* Alkaloids Mixture
Supplied as: A solution in acetonitrile (100 µg/ml of each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥3 years

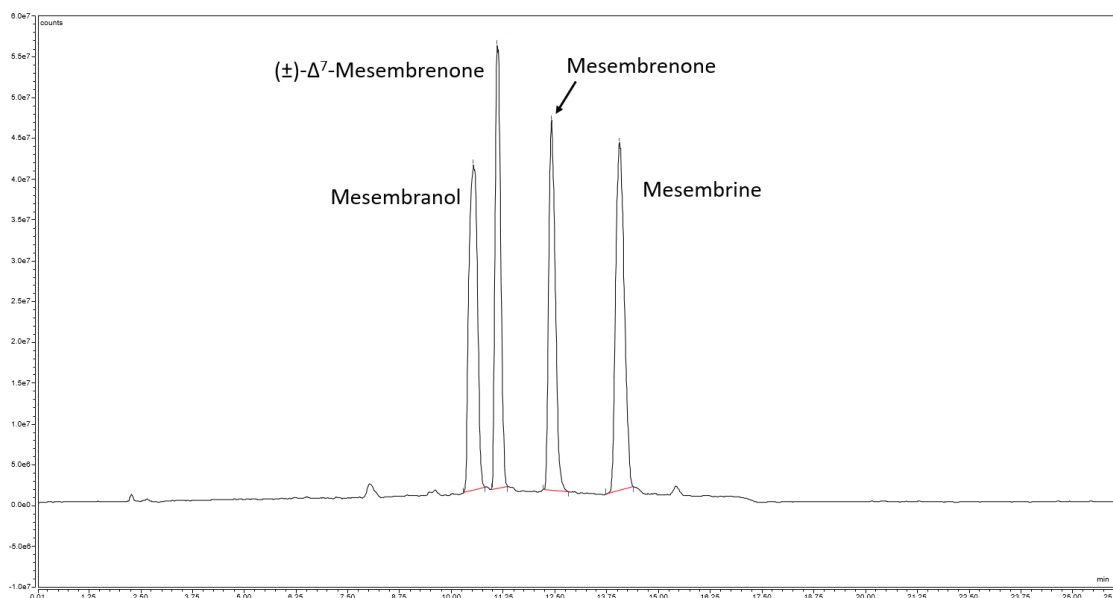
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Kanna Alkaloids Mixture (Item No. 45074) is an analytical reference standard mixture containing alkaloids found in the kanna (*S. tortuosum*) plant that can be used in the qualitative identification of key kanna alkaloids. The mixture contains mesembrine, mesembrenone, (±)- Δ^7 -mesembrenone, and mesembranol. 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 alkaloids found in the kanna (*S. tortuosum*) plant. Included are:

Item No. 45074	Kanna Alkaloids Mixture	
Item Name	CAS Number	Concentration (µg/ml)
Mesembrine	24880-43-1	100
Mesembrenone	468-54-2	100
(±)- Δ^7 -Mesembrenone	80287-15-6	100
Mesembranol	23544-42-5	100



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, 03/16/2026

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