

PRODUCT INFORMATION



Cannabis Flavonoids Mixture

Item No. 37832

Supplied as: A solution in N,N-dimethylacetamide (100 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

Cannabis Flavonoids Mixture (Item No. 37832) is a mixture of analytical reference standards containing flavonoids. The mixture contains kaempferol, vitexin, cannflavin A, and isoquercetin, among others. 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 and LC-MS instrumentation. This product is intended for research and forensic applications.

Contents

Item Number	Item Name	CAS Registry No.	Concentration
25143	Orientin	28608-75-5	100 µg/ml
15116	Vitexin	3681-93-4	100 µg/ml
31212	Isovitexin	38953-85-4	100 µg/ml
24926	Isoquercetin	482-35-9	100 µg/ml
10012600	Myricetin	529-44-2	100 µg/ml
10005169	Quercetin (hydrate)	849061-97-8	100 µg/ml
10004161	Luteolin	491-70-3	100 µg/ml
11852	Kaempferol	520-18-3	100 µg/ml
10010275	Apigenin	520-36-5	100 µg/ml
33395	Chrysoeriol	491-71-4	100 µg/ml
21970	Cannflavin A	76735-57-4	100 µg/ml

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, 02/28/2025

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