

PRODUCT INFORMATION

Phytocannabinoid Qualitative Mixture 10

Item No. 18791

Purity: ≥98% for each compound
Supplied as: A solution in acetonitrile (100 µg/ml of each compound)
Storage: -20°C
Stability: ≥3 years

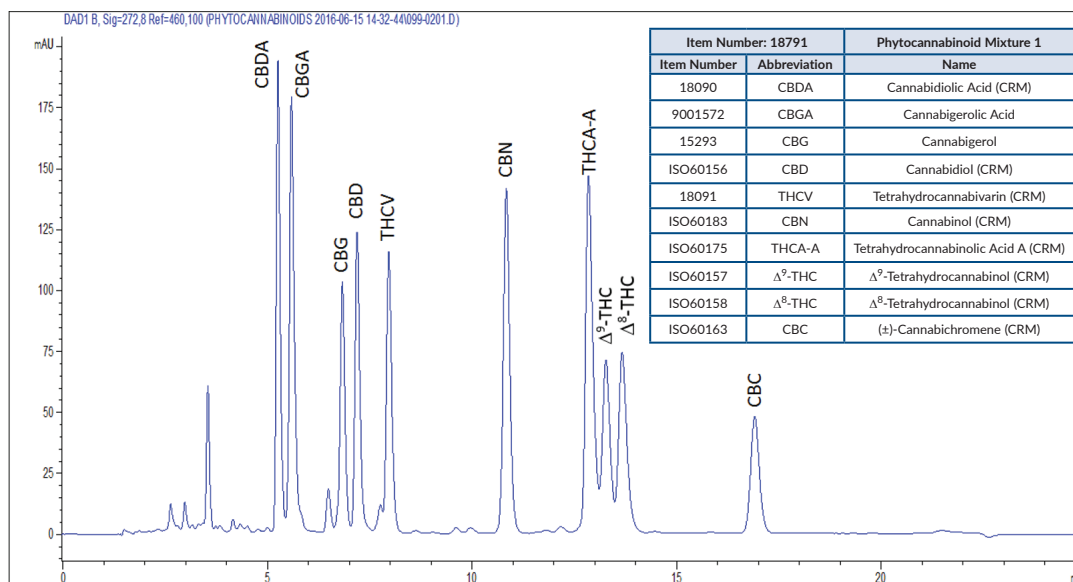
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Phytocannabinoid Qualitative Mixture 10 (Item No. 18791) is a mixture of compounds that can be isolated from plants in the genus *Cannabis*. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed to provide authentic standards for analyses using HPLC-UV detection; LC-MS analysis is also possible. This product is intended for forensic and research applications.

Contents

Each ampule contains at least 1 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, 08/16/2018

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