

PRODUCT INFORMATION



Colorado Pesticide Mixture Panel (CRM)

Item No. 39858

Synonym: CO Pesticide Mixture Panel
Supplied as: A 1 mg/ml solution (each pesticide) in acetonitrile
Fill volume: >1 ml
Storage: -20°C
Stability: ≥2 years
Special conditions: Vortex or sonicate before use

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

Description and Data

Colorado Pesticide Mixture Panel (CRM) (Item No. 39858) is a panel containing three certified reference material mixtures each composed of 9-10 pesticides for a total of 28 pesticides (>1 ml each ampule). The pesticides included in the panel correspond to the required list of pesticides by the state of Colorado for profiling of *Cannabis* and/or *Cannabis* products for residual pesticides. The panel is designed for use as a reference standard for these pesticides during routine analytical checks using GC- or LC-MS. The three mixtures are supplied in amber ampules in which the headspaces have been purged with argon. This product is intended for research and forensic applications.

Colorado Pesticide Mixture Panel (CRM) (Item No. 39858) includes acaricides, fungicides, and anthelmintics, as well as organophosphate, pyrethroid, and neonicotinoid insecticides, among others.

Item No.	Item Name	Pesticide	CAS No.
39493	Colorado Pesticide Mixture 1 (CRM)	Bifenthrin	82657-04-3
		Boscalid	188425-85-6
		Chlorpyrifos	2921-88-2
		lambda-Cyhalothrin	91465-08-6
		DDVP (Dichlorvos)	62-73-7
		Dimethoate	60-51-5
		Malathion	121-75-5
		Metalaxyl	57837-19-1
		Permethrins	52645-53-1
		Pyriproxyfen	95737-68-1
Item No.	Item Name	Pesticide	CAS No.
39494	Colorado Pesticide Mixture 2 (CRM)	Azoxystrobin	131860-33-8
		Bifenazate	149877-41-8
		Dinotefuran	165252-70-0
		Diuron (DCMU)	330-54-1
		Imazalil	35554-44-0
		Myclobutanil	88671-89-0
		Spinosad	168316-95-8
		Spiromesifen	283594-90-1
		Spirotetramat	203313-25-1

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, 10/27/2023

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

PRODUCT INFORMATION



Table continued:

Item No.	Item Name	Pesticide	CAS No.
39495	Colorado Pesticide Mixture 3 (CRM)	Abamectin	71751-41-2
		Carbaryl	63-25-2
		Clothianidin	210880-92-5
		Etoxazole	153233-91-1
		Imidacloprid	138261-41-3
		Propiconazole	60207-90-1
		Tebuconazole	107534-96-3
		Thiabendazole	148-79-8
Thiamethoxam	153719-23-4		

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