

PRODUCT INFORMATION



Montana Pesticide Mixture Panel (CRM)

Item No. 39635

Synonym: MT Pesticide Mixture Panel
Supplied as: A solution in acetonitrile (1 mg/ml of each compound)
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

Montana Pesticide Mixture Panel (CRM) (Item No. 39635) is a panel containing three certified reference material mixtures composed of 2-9 pesticides and chlormequat (CRM) for a total of 17 pesticides (>1 ml each ampule). The pesticides included in the panel correspond to the required list of pesticides by the state of Montana 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 and chlormequat are supplied in amber ampules in which the headspaces have been purged with argon. This product is intended for research and forensic applications.

Montana Pesticide Mixture Panel (CRM) (Item No. 39635) includes fungicides, acaricides, and plant growth regulators, as well as neonicotinoid, pyrethroid, and carbamate insecticides.

Item No.	Item Name	Pesticide	CAS Number
39362	Nevada Pesticide Mixture 1 (CRM)	Abamectin	71751-41-2
		Daminozide	1596-84-5
		Etoxazole	153233-91-1
		Imidacloprid	138261-41-3
		Myclobutanil	88671-89-0
		Paclobutrazol	76738-62-0
		Spinosad	168316-95-8
		Spirotetramat	203313-25-1
		Trifloxystrobin	141517-21-7
Item No.	Item Name	Pesticide	CAS Number
39363	Nevada Pesticide Mixture 2 (CRM)	Acequinocyl	57960-19-7
		Bifenazate	149877-41-8
		Bifenthrin	82657-04-3
		Cyfluthrin	68359-37-5
		Pyrethrins	8003-34-7
Item No.	Item Name	Pesticide	CAS Number
39365	Montana Pesticide Mixture (CRM)	Fenoxycarb	72490-01-8
		Imazalil	35554-44-0
Item No.	Item Name	Pesticide	CAS Number
39360	Chlormequat (chloride) (CRM)	Chlormequat (chloride)	999-81-5

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, 10/13/2023

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