

PRODUCT INFORMATION



Nevada Pesticide Mixture 1 (CRM)

Item No. 39362

Synonym: NV Pesticide Mixture 1
Supplied as: A solution in acetonitrile (1 mg/ml 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

Nevada Pesticide Mixture 1 (CRM) (Item No. 39362) is a certified reference material mixture containing pesticides. The mixture contains abamectin, daminozide, myclobutanil, and etoxazole, 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 and action limit testing on *Cannabis* and hemp samples using LC- or GC-MS instrumentation. This product is intended for research and forensic applications.

This mixture contains nine pesticides used during routine analytical checks and action limit testing on *Cannabis* and hemp samples. Included are:

Item No. 39362	Nevada Pesticide Mixture 1 (CRM)	
Item Name	CAS Number	Concentration (mg/ml)
Abamectin	71751-41-2	1.0
Daminozide	1596-84-5	1.0
Etoxazole	153233-91-1	1.0
Imidacloprid	138261-41-3	1.0
Myclobutanil	88671-89-0	1.0
Paclobutrazol	76738-62-0	1.0
Spinosad	168316-95-8	1.0
Spirotetramat	203313-25-1	1.0
Trifloxystrobin	141517-21-7	1.0

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