

# PRODUCT INFORMATION



## New Mexico Pesticide Mixture 3 (CRM)

Item No. 39368

**Synonym:** NM Pesticide Mixture 3  
**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

New Mexico Pesticide Mixture 3 (CRM) (Item No. 39368) is a certified reference material mixture containing pesticides. The mixture contains bifentazate, etoxazole, imazalil, and imidacloprid, 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 in 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 seven pesticides used during routine analytical checks and action limit testing on *Cannabis* and hemp samples. Included are:

Item No. 39368	New Mexico Pesticide Mixture 3 (CRM)	
Item Name	CAS Number	Concentration (mg/ml)
Bifentazate	149877-41-8	1.0
Etoxazole	153233-91-1	1.0
Imazalil	35554-44-0	1.0
Imidacloprid	138261-41-3	1.0
Myclobutanil	88671-89-0	1.0
Paclobutrazol	76738-62-0	1.0
Piperonyl Butoxide	51-03-6	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