

PRODUCT INFORMATION



Benzodiazepine Mixture 6 (CRM)

Item No. 39011

Supplied as: A solution in acetonitrile (100 µg/ml of each compound)

Storage: -20°C

Stability: ≥1 year

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

Benzodiazepine Mixture 6 (CRM) (Item No. 39011) is a certified reference material mixture containing benzodiazepines. The mixture contains alprazolam, clonazepam, diazepam, lorazepam, temazepam, and oxazepam. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in calibration or control solutions, method development, and as a reference standard for these substances during routine analytical checks using HPLC-UV and LC-MS instrumentation. Benzodiazepine Mixture 6 (CRM) contains compounds that are regulated as Schedule IV compounds in the United States. Benzodiazepine Mixture 6 (CRM) (Item No. 39011) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

The mixture contains alprazolam, clonazepam, diazepam, lorazepam, temazepam, and oxazepam. It is designed for use in calibration or control solutions, method development, and as a reference standard for these substances during routine analytical checks using HPLC-UV and LC-MS instrumentation.

Included are:

Item No. 39011	Benzodiazepine Mixture 6 (CRM)	
Item Name	CAS Number	Concentration (µg/ml)
Alprazolam	28981-97-7	100
Clonazepam	1622-61-3	100
Diazepam	439-14-5	100
Lorazepam	846-49-1	100
Temazepam	846-50-4	100
Oxazepam	604-75-1	100

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, 02/19/2025

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