

Mycotoxin Analytical Standards

We understand that identifying mold and mycotoxins, such as aflatoxins and ochratoxin A, that can naturally contaminate hemp and *Cannabis* during its cultivation, transportation, storage, and even processing, is crucial to the assurance of your quality and safety testing. Linked to kidney and liver damage, reproductive disorders, and immune suppression as well as some being identified as potent carcinogens, mycotoxins are a known risk within the food safety industry. Cayman's Mycotoxin Mixture CRM and single-component standards have been designed with consumer safety in mind to help you carry out mycotoxin analysis as part of your essential testing program.



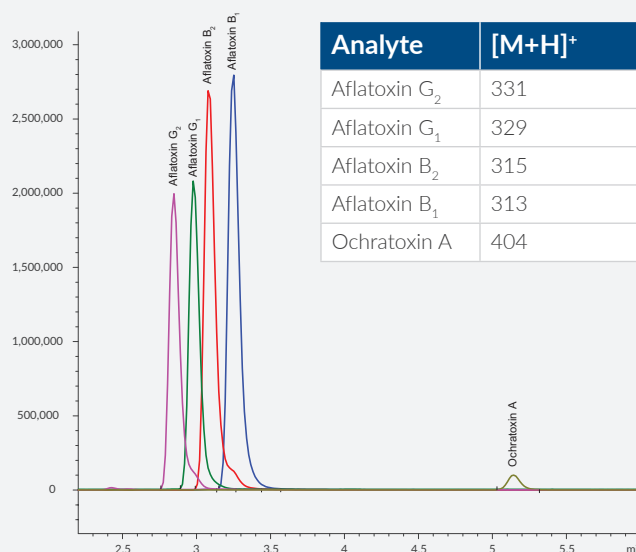
ACCREDITED
ISO/IEC 17025 #AT-1773
ISO 17034 #AR-1774

Simplify Mycotoxin Testing with Cayman's CRM Mixture

Mycotoxin Mixture 5 (CRM)

Item No. 28151

- Contains aflatoxin B₁, B₂, G₁, G₂, and ochratoxin A at 10 µg/ml of each compound
- Includes a certified concentration and uncertainty value for each analyte
- Intended for use as a calibration or secondary quality control standard
- Manufactured and tested to meet ISO 17034 and ISO/IEC 17025 international standards



Learn how pre-made mixtures improve quantitation accuracy in our app note at www.caymanchem.com/CRMmixtures

Single-Component Qualitative Standards

Item No.	Product Name
11293	Aflatoxin B ₁
11294	Aflatoxin B ₂
11295	Aflatoxin G ₁
11296	Aflatoxin G ₂
11297	Aflatoxin M ₁
11298	Aflatoxin M ₂

Item No.	Product Name
20437	Aflatoxin R ₀
11439	Ochratoxin A
16167	Ochratoxin B
20183	Ochratoxin C
11441	Sterigmatocystin
22070	O-methyl Sterigmatocystin

View a complete list of our analytical reference standards for hemp and *Cannabis* testing at www.caymanchem.com