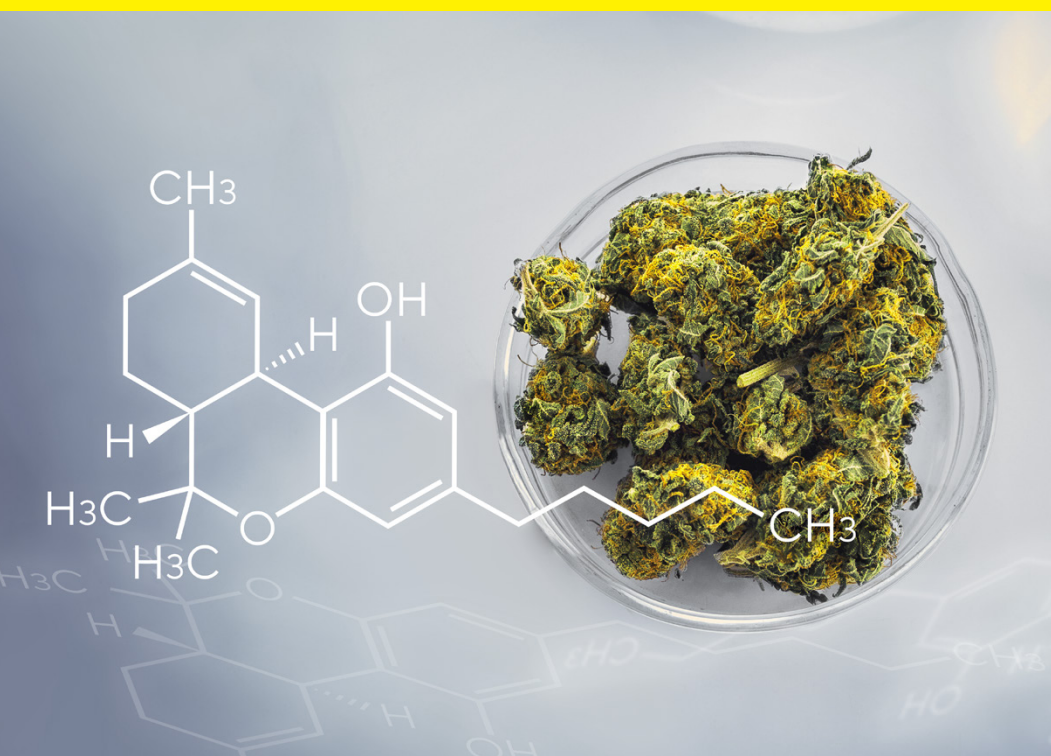


# SARTORIUS

## Make Quality Your Benchmark and Consumer Safety Your Priority

Sartorius Is Your Partner for Reliable Sample Preparation and Data Integrity for Cannabis Analysis.



Discover How  
Our Solutions  
Can Accelerate  
Your Analysis

States which have passed laws permitting usage of medicinal and recreational Cannabis products are also calling for rigorous testing of cannabis derivatives to ensure consumer safety.

Sartorius products not only support you in performing super-fast determination of your samples' moisture and weight, but also in the process of safely preparing your samples for analytical or microbiological analysis.

Get in touch with our specialists for expert advice on a specific need, or discover our solutions here:  
[www.sartorius.com/cannabis-testing](http://www.sartorius.com/cannabis-testing)

# Super-fast Determination of the Moisture Content of Cannabis as a Crucial Indicator of Potential Microbial Contamination

- Flexible, fully-automated, easy-to-use Infrared Moisture Analyzer MA160 for reliable and traceable results in the shortest possible time
- Get accurate measurements to determine the water content in cured cannabis, raw cannabis or edible cannabis products for quality control
- Quick and reliable measurement results in less than 30 minutes, rather than several hours of drying in the oven

## Analytical Cannabis Testing: Accurate and Consistent Lab Work Deserves Reliable and Precise Equipment and Best-in-Class Sample Prep

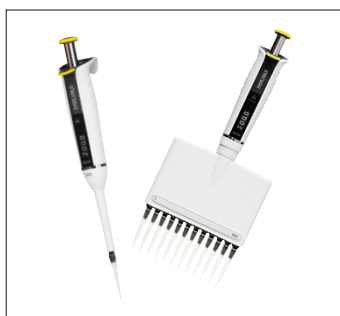
- Prepare your cannabis samples for analysis of pesticides, mycotoxins, residual solvents, heavy metals, terpenes and potency
- Get the highest sensitivity and performance from your analytical equipment with our state-of-the-art solutions

### Portfolio overview



#### Moisture Analyzers

Specify dry weight and moisture content of your samples using super-fast moisture analyzers.



#### Tacta® Pipettes

Enhance your workflow with robust and accurate pipettes, tips and bottle top dispensers.



#### Cubis® MSA Dosing System

Simplify your workflow with the semi-automatic preparation and documentation of highly reproducible calibration analytical standards for Cannabis testing.



#### Arium® Lab Water Systems

Use ultrapure water free of trace metals and organic components to avoid baseline noises and increase sensitivity of analytical experiments.



#### Minisart® Syringe Filters

Simply and safely remove solids from your sample with syringe filters.



#### Quintix® Balances

Choose from a range of intuitive operation lab balances to get the best weighing results and optimized workflows.