

## From Seed to Sale



Research & Development  
Analytical Testing

### Challenges faced

You care about the science and your data that can advance research in cannabis-based medicines and health products.

### Your needs – our solutions

You need accurate and reproducible standards and best in class sample prep for your analytical testing methods like HPLC and GC.

### Your product highlights

Quintix® and Secura® balances, Picus® and Tacta® pipettes, arium® lab water, Minisart® syringe filters



Grower  
Mass & Moisture Testing

### Challenges faced

Mold is the most problematic pathogen that not only poses health risks to consumers but financial risks to growers and manufacturers. Early in-process moisture testing to recommended levels (6-12%) helps safeguard against contamination.

### Your needs – our solutions

Check dry weight and moisture content of your samples using ultra fast moisture analyzers:

- small sample sizes avoid waste and reduce testing costs
- accurate and precise temperature control prevents the loss of terpenes and other essential volatile compounds

### Your product highlights

MA160, MA35 and MA37 moisture analyzers



Extractor  
Quality Control

### Challenges faced

When producing medicinal cannabis and hemp-derived products you need to perform many tests to ensure product safety and reduce liability concerns with tests for moisture content, bioburden, terpene profiling, residual solvents, heavy metals, potency and pesticide.

### Your needs – our solutions

Measure, weigh, dilute, filter and prepare accurate samples and standards for HPLC, ICPMS, GC analysis:

- reliable and intuitive sample prep tools to help meet standards for safety and labelling.
- economical sample prep solutions for cost-effective quality control

### Your product highlights

Picus® and Tacta® pipettes, Midi Plus™ Pipetting Controller, Bottle-top dispensers, Minisart® syringe filters, arium® lab water, Centrisart® centrifuges



Testing Labs  
Quality Control

### Challenges faced

Consumers need to know you are serious about their safety. Your data is your integrity. You need reproducible and accurate results that meet the standards for safety and labelling.

### Your needs – our solutions

Get best in class standards preparation with defined concentrations for quantitative HPLC by using Cubis® MSA Individual balance with Q-App software:

- intuitive navigation and full automation reduces risks of user inaccuracy
- automatic documentation keeps you compliant
- reliable planning keeps processes simple and efficient

### Your product highlights

Cubis® Dosing System with Q-App software for HPLC standards



Sartorius helps you build efficient work-flows, so you can concentrate on achieving high-quality results in your medicinal cannabis analysis. For further information visit: [www.sartorius.com/cannabis-testing](http://www.sartorius.com/cannabis-testing)