

Moisture Analysis Individual – Method Development

The Convenient and
Reliable Way to Obtain
Correct Results



Benefits

- Individual method development in the Sartorius Application Laboratory based on your requirements and samples
- Save time by bypassing method optimization and improving analysis times
- Reliability through guaranteed accurate measurements and optimal traceability

Product Information

The moisture content has a major impact on the quality (shelf life, suitability for storage) and on the usability of materials. Therefore, determining the moisture content is mandatory in many different areas. Oven drying is the most widespread method for measuring moisture content. This is a standard method in many in-house standard operating procedures (SOPs) and is incorporated into various statutory guidelines and regulations.

Description

The use of moisture analyzers is advisable to accelerate measurement times and to determine moisture content as a key criterion in supporting a process workflow. Because these analyzers work with infrared or microwave radiation to remove the moisture from a sample, they take only a few minutes to deliver a reading of the moisture result. To obtain reliably reproducible and traceable results, these analyzers should be aligned with your existing reference method (oven drying | Karl Fischer titration). Our specialists at the Sartorius Application Laboratories based in the USA and in Germany will be happy to optimize your parameters.

Easy

If you consult with us first, you can just send us your sample – we will take care of the rest. The Sartorius Application Laboratory will use your samples to develop individual methods that deliver optimal results for your measurements. Simply enter the pre-configured methods in your moisture analyzer and you can get started right away.

Fast

Avoid time-consuming reference measurements with a drying oven or Karl Fischer titration, and long-drawn-out method developments. You will also benefit over the long term from methods optimized in terms of measurement times, reproducibility and traceability, which will make your process more efficient.

Reliable

Whether your samples need to comply with individual test specifications, German Industrial Standards (DIN), ASTM standards or national pharmacopeia requirements – you can completely rely on our methods. We will correlate the results of your moisture analyzer with those of the standard method you require, and thus guarantee dependable, accurate results and traceability. This will also be documented in our test report.

Rely on our expertise. We have already successfully aligned over 15,000 different samples with the results of existing reference methods for our customers.

Description of the Individual Services Provided

- Detailed definition of requirements and consultation provided by our product specialists
- Analysis of your samples in the Sartorius Application Laboratory
 - Determination of reference values according to the relevant applicable standard (DIN, Section 64 of the German Food and Feed Code LFBG, ASTM, pharmacopeia, etc.) or according to your individual SOP using the following methods:
 - Oven drying method
 - Karl Fischer titration, coulometric, with or without a drying oven
- Method development using your samples to correlate a reference method with moisture analyzer results
 - Optimization of sample preparation and processing
 - Optimization of your sample application (application type and sample quantity)
 - Determination of the optimal parameters, depending on your need for accuracy or speed
- Evaluation of measurement results according to your specific requirements
- Preparation of a detailed test report



Services Available

Individual method development and correlation with your required reference method for guaranteed accurate measurement results and optimal traceability.

Order Number	MA35	MA37	MA100	MA160	Mark3	LMA200
YMD01B (incl. short report)	▪	▪				
YMD02B (incl. short report)				▪		
YMD01P (incl. premium report)			▪		▪	▪

Prerequisites and General Terms


- Following consultation, the customer sends the samples to our Sartorius Application Laboratory in New York, USA
- The customer provides full details of the particular sample and, if necessary, safety data sheets
- The customer accepts return of excess sample material, if applicable
- Sartorius reserves the right to refuse to accept delivery of individual hazardous materials and active ingredients
- Any test equipment provided may only be brought into contact with known sample material in order to prevent contamination by unknown materials
- Test equipment may be returned only in its original packaging
- This offer is available only in selected countries

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com