



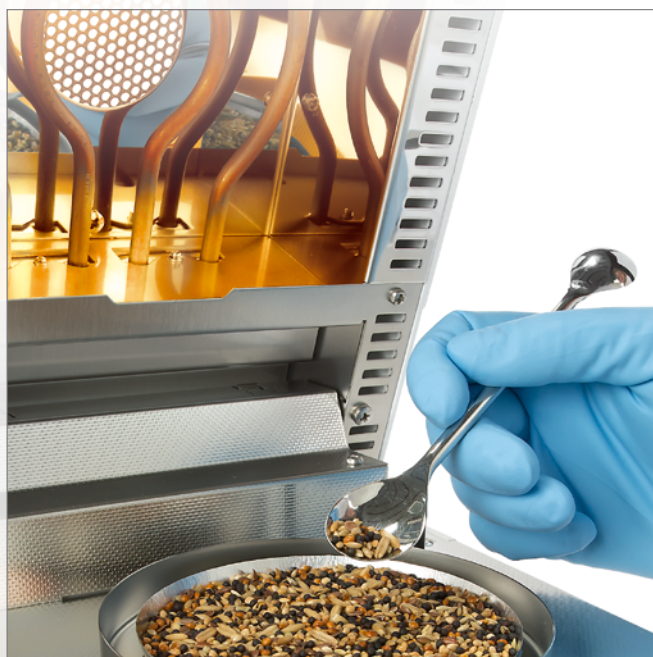
sartorius

MA35

## Rugged, Fully Automatic Infrared Moisture Analyzer for Use in Routine Operation

### Advantages:

- ▶ Time-saving alternative to a drying oven – quick measurement results in only 10-20 minutes instead of several hours
- ▶ Focused on the simple routine check
- ▶ Rugged device components, for use even in harsh environments
- ▶ Simple operation
- ▶ Best price-performance ratio with Sartorius quality – Made in Germany



sartorius  
130°C 10min  
50.36 g

## MA35 Infrared Moisture Analyzer

The MA35 is the ideal basic model for quick and reliable analysis of the material moisture of liquid, pasty and solid substances using the thermogravimetric method. Sartorius's decades of experience in the area of moisture analysis guarantees quick and reliable measurement results and high user-friendliness.



### Your Advantages at a Glance:

- ▶ **Ruggedness for all areas of application** The compact and rugged design, combined with great precision and measurement accuracy, make the MA 35 the ideal device for uniform measuring tasks. The selection of long-lasting and rugged components impervious to vibration and dirt guarantees reliable measurement results, even in a harsh environment. An ideal partner for use in production, incoming goods or in the laboratory.
- ▶ **Simple use** Sartorius's unique, fully automatic endpoint determination makes programming a switch-off criterion unnecessary. This saves time because, on the one hand, time-consuming optimizations are no longer required, and on the other, operation of the device is simplified so that it can be used competently by anyone. This is made possible by continuous monitoring of the drying cycle by the MA35. This monitoring then stops the measurement fully automatically using a switch-off criterion that is calculated based on the drying curve as soon as the sample weight is constant, and where despite heating, no additional loss of weight is detectable. The measurement accuracy required for this is provided by a Sartorius weighing system with readability to 1 mg, which has been specially optimized for use in high temperature ranges.
- ▶ **Quick measurement results** For sample heating, the MA35 features two powerful metal tubular-shaped heating elements generating 360 watts of power. These elements facilitate quick and uniform heating of samples and thus short measurement times, which means a significant amount of time is saved compared to the drying oven because the finished end result can already simply be read within a few minutes.



### Your Contact

[www.sartorius.com](http://www.sartorius.com)

Specifications subject to change without notice.  
Copyright Sartorius Lab Instruments GmbH & Co. KG.  
Printed in the EU on paper bleached without chlorine.  
Publication No.: W--1524-e170401  
Order No.: 98649-022-97  
Ver. 04 | 2017