



Moisture Analyzer Portfolio

Overview of Standard Analyzers



Model	MA35	MA37	MA160
Features	<ul style="list-style-type: none"> Rugged entry-level model Easy operation For repetitive measurements 	<ul style="list-style-type: none"> Faster analysis (due to more powerful heating unit than that in the MA35) Touch screen Easy to clean Direct data transfer Function testing 	<ul style="list-style-type: none"> AURI metal tube heater for the highest measurement speeds Touch screen Password protection Memory for methods Method Development Assistant Memory for results Easy to clean Direct data transfer Function testing
Moisture measurement range in %	> 1	> 1	> 1
Typical uses	Moisture analysis of recurrent samples	Moisture analysis of recurrent samples	<ul style="list-style-type: none"> Moisture analysis of recurrent samples Moisture analysis of the widest variety of samples
Typical industrial sectors	<ul style="list-style-type: none"> Sewage treatment plants Biogas facilities Manufacturing Paints, lacquers, varnishes, dyes 	<ul style="list-style-type: none"> Food and beverage (incoming quality control) Building materials and wood industries Animal feed 	<ul style="list-style-type: none"> Chemical analytical labs QA QM Food & beverage Manufacturing Research and development

Overview of Premium Analyzers



Model	MA100	LMA100	LMA200
Features	<ul style="list-style-type: none"> Analytical weighing system (0.1 mg) Different, interchangeable heat sources Memory for methods Easy operation Password protection Various heating programs Various shutoff criteria The most flexible range of applications Statistics function 	<ul style="list-style-type: none"> Analytical weighing system (0.1 mg) Modular design consisting of operating and heating units (up to 4) User management Memory for methods Large sample chamber Special weighing pan for plastic samples Syringe and concentration modes Statistics function and memory for results 	<ul style="list-style-type: none"> Analytical weighing system (0.1 mg) User management Microwave drying The fastest measurements (many take less than 1 – 2 min.) Carousel-type sample holder Memory for methods Statistics function and memory for results
Moisture measurement range in %	>0.01	>0.01	>8
Typical uses	<ul style="list-style-type: none"> Analysis of samples with a low moisture content (below 1%), such as plastic Moisture analysis of the widest variety of samples 	<ul style="list-style-type: none"> Analysis of samples with a low moisture content (below 1%), such as plastic Special applications (e.g., highly volatile substances) 	<ul style="list-style-type: none"> Analysis of samples with a high moisture content (liquids, pastes, etc.)
Typical industrial sectors	<ul style="list-style-type: none"> Processing of plastics Pharma (granules and powder) Food and beverage R&D Paper and textile industries 	<ul style="list-style-type: none"> Processing of plastics Pharma (granules and powder) Food and beverage Chemical industry Automotive industry 	<ul style="list-style-type: none"> Food and beverage (dairies) Pharma (extracts, homeopathic products) Water-based paints and other coatings Cosmetics industry

Germany
 Sartorius Lab Instruments
 GmbH & Co. KG
 Otto-Brenner-Strasse 20
 37079 Goettingen
 Phone +49.551.308.0
 Fax +49.551.308.3289
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--1521-e170602
 Order No.: 98649-020-30
 Ver. 06 | 2017