

## Moisture Analyzers in 2020 | 2021

### Overview of Standard Analyzers



Model	MA 35	MA 37	MA 160
Features	<ul style="list-style-type: none"> <li>▪ Rugged entry-level model</li> <li>▪ Easy operation</li> <li>▪ For repetitive measurements</li> </ul>	<ul style="list-style-type: none"> <li>▪ Faster analysis (due to more powerful heating unit than that in the MA35)</li> <li>▪ Touch screen</li> <li>▪ Easy to clean</li> <li>▪ Direct data transfer</li> <li>▪ Function testing</li> </ul>	<ul style="list-style-type: none"> <li>▪ AURI metal tube heater for the highest measurement speeds</li> <li>▪ Touch screen</li> <li>▪ Password protection</li> <li>▪ Memory for methods</li> <li>▪ Method Development Assistant</li> <li>▪ Memory for results</li> <li>▪ Easy to clean</li> <li>▪ Direct data transfer</li> <li>▪ Function testing</li> </ul>
Moisture measurement range in %	>1	>1	>1
Typical uses	<ul style="list-style-type: none"> <li>▪ Moisture analysis of recurrent samples</li> </ul>	<ul style="list-style-type: none"> <li>▪ Moisture analysis of recurrent samples</li> </ul>	<ul style="list-style-type: none"> <li>▪ Moisture analysis of recurrent samples</li> <li>▪ Moisture analysis of the widest variety of samples</li> </ul>
Typical industrial sectors	<ul style="list-style-type: none"> <li>▪ Sewage treatment plants</li> <li>▪ Biogas facilities</li> <li>▪ Manufacturing</li> <li>▪ Paints, lacquers, varnishes, dyes</li> </ul>	<ul style="list-style-type: none"> <li>▪ Food and beverage (incoming quality control)</li> <li>▪ Building materials and wood industries</li> <li>▪ Animal feed</li> </ul>	<ul style="list-style-type: none"> <li>▪ Chemical analytical labs</li> <li>▪ QA   QM</li> <li>▪ Food &amp; beverage</li> <li>▪ Manufacturing</li> <li>▪ Research and development</li> </ul>

# Overview of Premium Analyzers



Model	MA 100	LMA 100	LMA 200
Features	<ul style="list-style-type: none"> <li>▪ Analytical weighing system (0.1 mg)</li> <li>▪ Different, interchangeable heat sources</li> <li>▪ Memory for methods</li> <li>▪ Easy operation</li> <li>▪ Password protection</li> <li>▪ Various heating programs</li> <li>▪ Various shutoff criteria</li> <li>▪ The most flexible range of applications</li> <li>▪ Statistics function</li> </ul>	<ul style="list-style-type: none"> <li>▪ Analytical weighing system (0.1 mg)</li> <li>▪ Modular design consisting of operating and heating units (up to 4)</li> <li>▪ User management</li> <li>▪ Memory for methods</li> <li>▪ Large sample chamber</li> <li>▪ Special weighing pan for plastic samples</li> <li>▪ Syringe and concentration modes</li> <li>▪ Statistics function and memory for results</li> </ul>	<ul style="list-style-type: none"> <li>▪ Analytical weighing system (0.1 mg)</li> <li>▪ User management</li> <li>▪ Microwave drying</li> <li>▪ The fastest measurements (many take less than 1–2 min.)</li> <li>▪ Carousel-type sample holder</li> <li>▪ Memory for methods</li> <li>▪ Statistics function and memory for results</li> </ul>
Moisture measurement range in %	> 0.01	> 0.01	> 8
Typical uses	<ul style="list-style-type: none"> <li>▪ Analysis of samples with a low moisture content (below 1%), such as plastic</li> <li>▪ Moisture analysis of the widest variety of samples</li> </ul>	<ul style="list-style-type: none"> <li>▪ Analysis of samples with a low moisture content (below 1%), such as plastic</li> <li>▪ Special applications (e.g., highly volatile substances)</li> </ul>	<ul style="list-style-type: none"> <li>▪ Analysis of samples with a high moisture content (liquids, pastes, etc.)</li> </ul>
Typical industrial sectors	<ul style="list-style-type: none"> <li>▪ Processing of plastics</li> <li>▪ Pharma (granules and powder)</li> <li>▪ Food and beverage</li> <li>▪ R&amp;D</li> <li>▪ Paper and textile industries</li> </ul>	<ul style="list-style-type: none"> <li>▪ Processing of plastics</li> <li>▪ Pharma (granules and powder)</li> <li>▪ Food and beverage</li> <li>▪ Chemical industry</li> <li>▪ Automotive industry</li> </ul>	<ul style="list-style-type: none"> <li>▪ Food and beverage (dairies)</li> <li>▪ Pharma (extracts, homeopathic products)</li> <li>▪ Water-based paints and other coatings</li> <li>▪ Cosmetics industry</li> </ul>

## Germany

Sartorius Lab Instruments GmbH & Co. KG  
 Otto-Brenner-Straße 20  
 37079 Göttingen  
 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](http://www.sartorius.com)