



MA160

In 3 Steps to right and compliant results

Scope

Sartorius' almost 30 years proven fully automatic endpoint determination for our moisture analyzers enables the devices to stop the measurement automatically as soon as a constant sample weight is reached. Thus, it is no longer necessary to program a shutoff parameter. The new MA160 even goes one step further. The integrated Method-Development-Assistant suggests a suitable temperature depending on the characteristics of your sample.



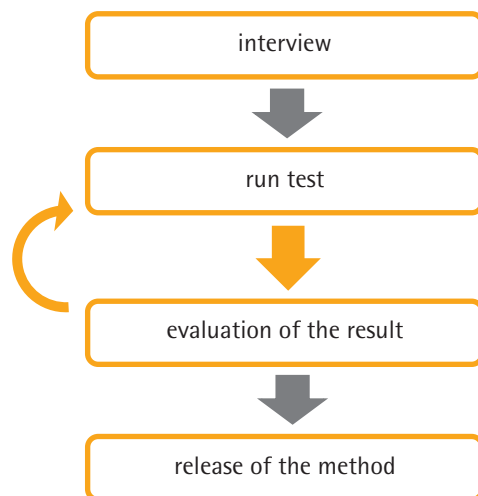
Technical Note

Executive Summary

The most common reference methods for the determination of the moisture content are the determination with the drying oven, vacuum oven or a titration procedure. No matter which method is used, it is very important to measure the right moisture content of each sample material. The biggest disadvantage of all these standard methods is that they are very time-consuming and sometimes difficult to proceed. Therefore, it is hard to measure the moisture content during production as you have a time delay of sometimes up to one day until you get your results. This you can overcome by using a MA160 rapid IR-Moisture Analyzer from Sartorius. To determine the moisture content in a standard complying way you have to adjust your MA160 method parameters to your standard reference method. This can be achieved through the modification of the measuring parameters. By doing this, you can be sure to get results comparable to your standard method in minimum time. The MA160 is the first Moisture Analyzer on the market with an included assistant function in order to make it as easy as possible to find the parameters giving you comparability to your standard method. On the basis of the main characteristics of your sample the MA160 supports you to find the right method parameters for your sample in just 3 easy steps.

Method development in 3 Steps

- ▶ Step 1: interview session
- ▶ Step 2: method test
- ▶ Step 3: release the new method



Learn more about the new MA160 Moisture Analyzer:
www.sartorius.com/en/moisture-analyzer/ma160/

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius.com

USA Toll-free +1.800.635.2906
UK +44.1372.737159
France +33.1.70.62.50.00
Italy +39.039.4659.1
Spain +34.913.586.095
Russian Federation +7.812.327.53.27
Japan +81.3.3740.5408

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--1140-e160801
Order No.: 98649-021-41
Ver. 08 | 2016