



## Mark 3 LTE Feuchtebestimmer



### Anwendungs- beispiel

Papier und Zellstoff

Mark 3 ermöglicht eine schnelle und präzise Feuchtemessung und erlaubt eine Anpassungen der Produktionszyklen innerhalb von Minuten statt Stunden.

- Über 10.000 verschiedenen Proben sind bereits auf die Referenzmethode abgeglichen.
- Keine weitere Programmierung durch den Anwender notwendig
- Ein Knopfdruck zum Starten des Tests
- 2 Jahre Garantie
- 300 Programme können gespeichert werden
- Interner Drucker für direkten Ausdruck der Ergebnisse am Platz

Material	Ofen Referenz	Mark 3	Testzeit	Std. Abw.
Schwarzlauge	67,64% S	67,58 % S	6,3 Min.	0,154
Deckblatt, leicht	6,47% M	6,33 % S	3,2 Min.	0,097
Deckblatt, schwer	6,33% M	6,37 % M	3,3 Min.	0,041
Baumwollfasern, nass	61,66% M	61,59 % M	10,7 Min.	0,117
Klebestreifen	3,45% M	3,37 % M	6,6 Min.	0,050
Polypropylenfasern	0,45% M	0,44% M	7,6 Min.	0,071
Zeitungspapier	5,50% M	5,58% M	8,8 Min.	0,214
Polyesterfaser	1,47% M	1,54% M	1,6 Min.	0,030
Hartholz Zellstoff	10,18% M	10,01% M	12,2 Min.	0,152
Sandpapier, Körnung 80	2,01% M	2,10 % M	8,6 Min.	0,070
Sägemehl o. Holzmehl	5,76% M	5,64% M	4,0 Min.	0,103
Holzbrei	20,02% S	19,98% S	14,8 Min.	0,094
Zitronenholz	30,53% M	30,44% M	14,0 Min.	0,091
Holzspäne	5,61% M	5,54% M	5,0 Min.	0,056
Holzfaslerplatte	5,34% M	5,39% M	7,2 Min.	0,042

Tabelle zeigt die Korrelation der Ergebnisse zwischen Konvektions- oder Vakuum-Ofen und dem Mark 3 Feuchtebestimmer.

# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Weender Landstrasse 94-108  
37075 Goettingen  
Phone +49.551.308.0  
Fax +49.551.308.3289

### France & Suisse Romande

Sartorius France S.A.S.  
ZI Les Paluds  
Avenue de Jouques - CS 31090  
13781 Aubagne Cedex  
Phone +33.442.845600  
Fax +33.442.846545

### Austria

Sartorius Austria GmbH  
Franzosengraben 12  
1030 Vienna  
Phone +43.1.7965760.0  
Fax +43.1.7965760.24

### Belgium

Sartorius Belgium N.V.  
Leuvensesteenweg, 248/B  
1800 Vilvoorde  
Phone +32.2.756.06.71  
Fax +32.2.253.45.95

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki  
Phone +358.9.755.951  
Fax +358.9.755.95.292

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi  
Phone +3623.457.227  
Fax +3623.457.147

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11  
Phone +353.1.8089050  
Fax +353.1.8089388

### Italy

Sartorius Italy S.r.l.  
Via dell'Antella, 76/A  
50012 Antella-Bagno a Ripoli (FI)  
Phone +39.055.63.40.41  
Fax +39.055.63.40.526

### Poland

Sartorius Poland sp.z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn  
Phone +48.61.6473830  
Fax +48.61.6473839

### Russian Federation

LLC "Sartorius ICR"  
and LLC "Biohit"  
Uralskaya str. 4, Lit. B  
199155, Saint-Petersburg  
Phone +7.812.327.5.327  
Fax +7.812.327.5.323

### Spain & Portugal

Sartorius Spain S.A.  
Offices in Madrid:  
C/ Isabel Colbrand, 10-12, of. 70  
28050 Madrid  
Phone Spain +34.902.123.367  
Phone Portugal +351.800.855.800  
Fax Spain +34.91.358.84.85  
Fax Portugal +351.800.855.799

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)  
Phone +41.44.746.50.00  
Fax +41.44.746.50.50

### U.K.

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ  
Phone +44.1372.737159  
Fax +44.1372.729927

## America

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
Fax +1.631.254.4253

### Argentina

Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires  
Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Av. Dom Pedro I, 241  
Bairro Vila Pires  
Santo André  
São Paulo  
Cep 09110-001  
Phone +55.11.4451.6226  
Fax +55.11.4451.4369

### Canada

Sartorius Canada Inc.  
2179 Dunwin Drive #4  
Mississauga, ON L5L 1X2  
Phone +1.905.569.7977  
Toll-Free +1.800.668.4234  
Fax +1.905.569.7021

### Mexico

Sartorius de México S.A. de C.V.  
Circuito Circunvalación Poniente  
No. 149  
Ciudad Satélite  
53100, Estado de México  
México  
Phone +52.5555.62.1102  
Fax +52.5555.62.2942

## Asia | Pacific

### Australia

Sartorius Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175  
Phone +61.3.8762.1800  
Fax +61.3.8762.1828

### China

Sartorius Scientific  
Instruments (Beijing) Co., Ltd.  
33 Yu An Road, Airport Industrial Park  
Zone B,  
Shunyi District, Beijing 101300,  
P.R.China  
Phone +86.10.8042.6300  
Fax +86.10.8042.6486

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwung Tong  
Kowloon, Hong Kong  
Phone +852.2774.2678  
Fax +852.2766.3526

### India

Sartorius Weighing India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra,  
Nelamangala Tq  
562 123 Bangalore, India  
Phone +91.80.4350.5250  
Fax +91.80.4350.5253

### Japan

Sartorius Stedim Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan  
Phone +81.3.4331.4300  
Fax +81.3.4331.4301

### Malaysia

Sartorius Malaysia Sdn. Bhd  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia  
Phone +60.3.8996.0622  
Fax +60.3.8996.0755

### Singapore

Sartorius Singapore Pte. Ltd  
1 Science Park Road,  
The Capricorn, #05-08A,  
Singapore Science Park II  
Singapore 117528  
Phone +65.6872.3966  
Fax +65.6778.2494

### South Korea

Sartorius Korea Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, BunDang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400  
Phone +82.31.622.5700  
Fax +82.31.622.5799

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310  
Phone +66.2643.8361-6  
Fax +66.2643.8367



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