



Mark 3 LTE Feuchtebestimmer



Anwendungs- beispiel

Lebensmittelindustrie
– Snacks

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.
Brezel	2,62% M	2,61% M	3,4 Min.	0,037
Schweineschwarte	3,77% M	3,77% M	5,9 Min.	0,220
Kartoffelchips	1,41% M	1,40% M	6,9 Min.	0,035
Käseflips	3,07% M	3,05% M	5,6 Min.	0,041
Maiskörner	1,26% M	1,25% M	4,0 Min.	0,032
Minisalami	24,92% M	25,02% M	8,9 Min.	0,181
Maischips	2,24% M	2,22% M	8,5 Min.	0,074
Erdnussriegel	2,21% M	2,25% M	9,2 Min.	0,064
Pistazien	2,73% M	2,70% M	8,4 Min.	0,039
Erdbeergebäck	13,50% M	13,74% M	4,0 Min.	0,179
Käsecracker	3,88% M	3,85% M	4,5 Min.	0,068
Bananenchips	1,43% M	1,49% M	10,5 Min.	0,031
Schokoladenkuchen	6,69% M	6,73% M	7,0 Min.	0,119
Salzcracker	5,21% M	5,22% M	7,9 Min.	0,021

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

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