



Mark 3 Feuchtebestimmer



Anwendungs- beispiel

Kunststoffe

Der Mark 3 Feuchtebestimmer entspricht dem ASTM Standard D6980 „Standard Testmethode für die Bestimmung von Feuchte durch Gewichtsabnahme“. Eine schnelle und präzise Feuchte- oder Feststoffmessung von lösemittelhaltigen oder wasserbasierten Materialien mit dem Mark 3 erlaubt eine Anpassungen der Produktionszyklen innerhalb von Minuten statt Stunden.

- Über 4.000 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

Generisch	Spez. Grad	KF Titration	Mark 3	Testzeit	Std. Abw.
ABS	Cyclac FR15	0,034%	0,032%	5,0 Min.	0,002
Acetal	Delrin 500P	0,020%	0,020%	4,0 Min.	0,001
ABS	Lustran LK279	0,061%	0,060%	10,0 Min.	0,001
Acryl	Acrylite S101	0,045%	0,042%	7,0 Min.	0,001
Nylon 6	Akulon Ultraflow K	0,031%	0,031%	4,3 Min.	0,002
Nylon 6	Capron 8234	0,088%	0,081%	9,0 Min.	0,003
Nylon 6	Durethan BG30X	0,071%	0,072%	5,0 Min.	0,004
Nylon 6/6	Zytel ST 801	0,115%	0,114%	12,5 Min.	0,006
Nylon 6/6	Chemlon 133 GH	0,064%	0,063%	8,0 Min.	0,004
Nylon 11	Rilsan Besno	0,020%	0,018%	3,3 Min.	0,003
PC	Lexan 500	0,015%	0,016%	4,0 Min.	0,002
PC	Makrolon RX-2530	0,013%	0,014%	4,0 Min.	0,001
LCP	L Zenite 55201	0,009%	0,010%	5,0 Min.	0,001
LCP	Vectra E473i	0,016%	0,016%	4,7 Min.	0,002
PET	Kosa 1101	48 ppm	49 ppm	5,0 Min.	0,001
PET	Traytuff	0,015%	0,014%	5,0 Min.	0,001
PET	RyniteRE 5231	0,009%	0,010%	4,9 Min.	0,001
PBT	Celanex 6500	0,022%	0,021%	3,7 Min.	0,002
PBT	Valox 325	0,010%	0,010%	4,0 Min.	0,001
HDPE	Petrothene LR7	0,005%	0,006%	4,0 Min.	0,001
PP	Thermofil P6-30	0,020%	0,019%	7,4 Min.	0,001
PS	Styron 6079	0,009%	0,011%	4,0 Min.	0,002
PSU	Udel GF-120	0,031%	0,031%	6,8 Min.	0,002
PU	Isoplast 2303C	0,015%	0,017%	5,0 Min.	0,001

Tabelle zeigt die Korrelation der Ergebnisse zwischen der Karl-Fischer-Titration 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