



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