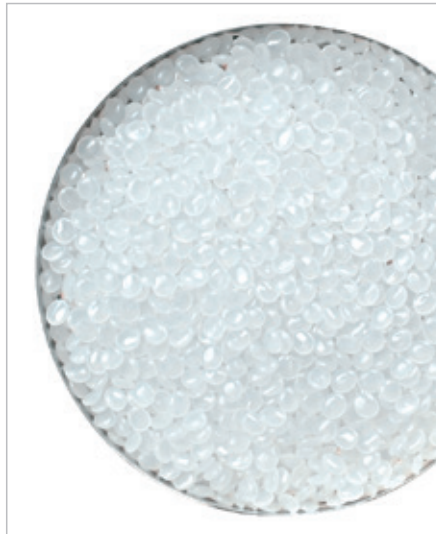




Mark 3 Moisture Analyzer



Application Examples

Plastics Industry

The Mark 3 conforms to ASTM standard D6980, "Standard Test Method for Determination of Moisture in Plastics by Loss in Weight". Fast, accurate moisture measurement with the Mark 3 allows adjustments to be made in the processing cycle in minutes rather than hours.

- Over 4000 different resin grades already correlated to Karl Rscher Titration
- No pre-programming required on the part of the operator
- One button operation to perform a test
- Two year complete parts and labor warranty
- Verifies performance of dryer systems
- 300 resin programs can be stored and recalled by name
- Hard copy results with standard internal printer

Generic	Specific Grade	KF Titration	Mark 3 Results	Test Times	STD DEV
ABS	Cycolac FR15	.034%	.032%	5.0 Min.	.002
Acetal	Delrin 500P	.020%	.020%	4.0 Min.	.001
ABS	Lustran LK279	.061%	.060%	10.0 Min.	.001
Acrylic	Acrylite S101	.045%	.042%	7.0 Min.	.001
Nylon 6	Akulon Ultraflow K	.031%	.031%	4.3 Min.	.002
Nylon 6	Capron 8234	.088%	.081%	9.0 Min.	.003
Nylon 6	Durethan BG30X	.071%	.072%	5.0 Min.	.004
Nylon 6/6	Zytel ST 801	.115%	.114%	12.5 Min.	.006
Nylon 6/6	Chemlon 133 GH	.064%	.063%	8.0 Min.	.004
Nylon 11	Rilsan Besno	.020%	.018%	3.3 Min.	.003
PC	Lexan 500	.015%	.016%	4.0 Min.	.002
PC	Makrolon RX-2530	.013%	.014%	4.0 Min.	.001
LCP	Zenite 55201	.009%	.010%	5.0 Min.	.001
LCP	Vectra E473i	.016%	.016%	4.7 Min.	.002
PET	Kosa 1101	48 ppm	49 ppm	5.0 Min.	.001
PET	Traytuff	.015%	.014%	5.0 Min.	.001
PET	RyniteRE 5231	.009%	.010%	4.9 Min.	.001
PBT	Celanex 6500	.022%	.021%	3.7 Min.	.002
PBT	Valox 325	.010%	.010%	4.0 Min.	.001
HDPE	Petrothene LR7	.005%	.006%	4.0 Min.	.001
PP	Thermofil P6-30	.020%	.019%	7.4 Min.	.001
PS	Styron 6079	.009%	.011%	4.0 Min.	.002
PSU	Udel GF-120	.031%	.031%	6.8 Min.	.002
PU	Isoplast 2303C	.015%	.017%	5.0 Min.	.001

Table demonstrates correlation results between Karl Fischer coulometric titration and the Mark 3 moisture analyzer.

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
Franzengraben 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