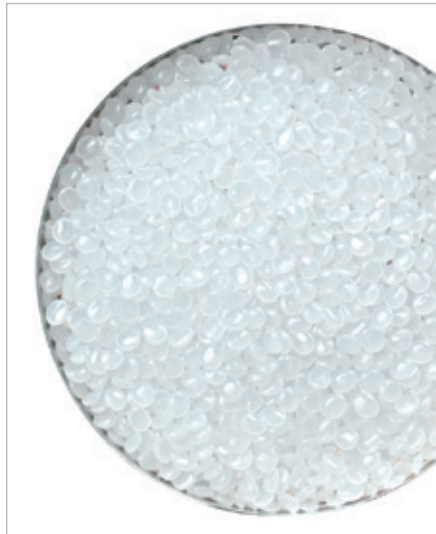




# Mark 3 Moisture Analyzer

Moisture Analysis



## 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	Cyclac 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](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  
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](http://www.sartorius.com)