

Mark 3 Moisture Analyzer

Application Examples
Plastics Industry



Benefits

- Over 4000 different resin grades already correlated to Karl Fischer 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

Product Information

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.

Correlation to Karl Fischer Titration

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 66	Zytel ST 801	.115%	.114%	12.5 min.	.006
Nylon 66	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.

Germany

Sartorius Lab Instruments GmbH & Co. KG
 Otto-Brenner-Straße 20
 37079 Göttingen
 Phone +49 551 308 0

USA

Sartorius Corporation
 565 Johnson Avenue
 Bohemia, NY 11716
 Phone +1 631 254 4249
 Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com