

# Mark 3 LTE Moisture Analyzer

Application Examples  
Chemicals &  
Minerals Industry



## Benefits

- Thousands of materials already correlated to reference methods
- No pre-programming required on the part of the operator
- One button operation to perform a test
- Two year complete parts and labor warranty
- 60 programs can be stored and recalled by name
- Hard copy results with optional internal printer

## Product Information

The Mark 3 LTE is the perfect analyzer for your production floor. Fast, accurate moisture measurement allows adjustments to be made in the production cycle in minutes rather than hours. Reliable and affordable, the Mark 3 LTE can save you time and money.

# Correlation to Oven Data

Type Material	Oven Reference	Mark 3 LTE Results	Test Times	STD DEV
TiO2	79.72% S	79.65% S	6.0 Min.	.048
PVA (Polyvinyl Acetate)	7.05% S	7.11% S	12.0 Min.	.069
Carbon Black Solution	7.20% S	7.26% S	9.9 Min.	.102
Sodium Phosphate	0.945% M	0.95% M	4.1 Min.	.010
Cathode Powder	4.25% M	4.28% M	3.9 Min.	.009
Cathode Slurry	44.60% S	44.56% S	3.5 Min.	.170
Soda Ash	0.261% M	0.27% M	5.9 Min.	.021
EVOH	0.130% M	0.12% M	12.7 Min.	.005
Urea Prills	0.240% M	0.24% M	10.3 Min.	.006
Calcium Citrate	4.91% M	4.81% M	4.7 Min.	.037
Ammonium Sulfate	0.30% M	0.32% M	7.8 Min.	.045
Solvent-Based Caulk	65.52% M	65.58% M	5.6 Min.	.142
Lime Cake	30.53% M	30.44% M	14.0 Min.	.091
Coal	6.75% M	6.71% M	10.8 Min.	.201
Magnesium Stearate	3.79% M	3.76% M	7.8 Min.	.030


Table demonstrates correlation results between convection or vacuum oven and the Mark 3 LTE 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](http://www.sartorius.com)