

# Mark 3 Moisture Analyzer

Application Examples  
Cosmetics Industry



## Benefits

- Over 10,000 different 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
- 300 materials can be stored and recalled by name
- Hard copy results with standard internal printer

## Product Information

Fast, accurate moisture measurement with the Mark 3 allows adjustments to be made in the processing cycle in minutes rather than hours.

# Correlation Results

Type Material	Oven Reference	Mark 3 Results	Test Times	STD DEV
Nail Polish	31.28% S	31.31% S	5.4 Min.	0.073
Nitro-base Solution	27.70% S	27.68% S	7.0 Min.	0.120
Polish Remover	0.273% S	0.268% S	15.3 Min.	0.034
Talc	0.070% M	0.069% M	2.4 Min.	0.004
Styling Spritz	9.43% S	9.44% S	5.5 Min.	0.070
Anti-Bacterial Soap	43.52% M	43.497% M	5.6 Min.	0.121
Body Lotion	17.51% M	17.45 % M	12.0 Min.	0.160
Collagen Slurry	93.54% M	93.555% M	11.6 Min.	0.030
Body Wash	21.57% S	21.62% S	11.5 Min.	0.151

Table demonstrates correlation results between convection or vacuum oven and the Mark 3 moisture analyzer (LMA100P).

## 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)