

LMA100 Moisture Analyzer

Operating Instructions



sartorius

1.

가

.

가

.

,

.

가

.

,

.

.

.

2.

Heating

,

1ea

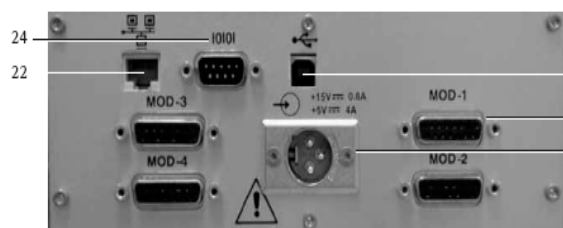
Tweezers

Operating instructions manual

LMA100 ()



Control Module connector panel



Heater Module connector panel



No. Designation

Heater Module

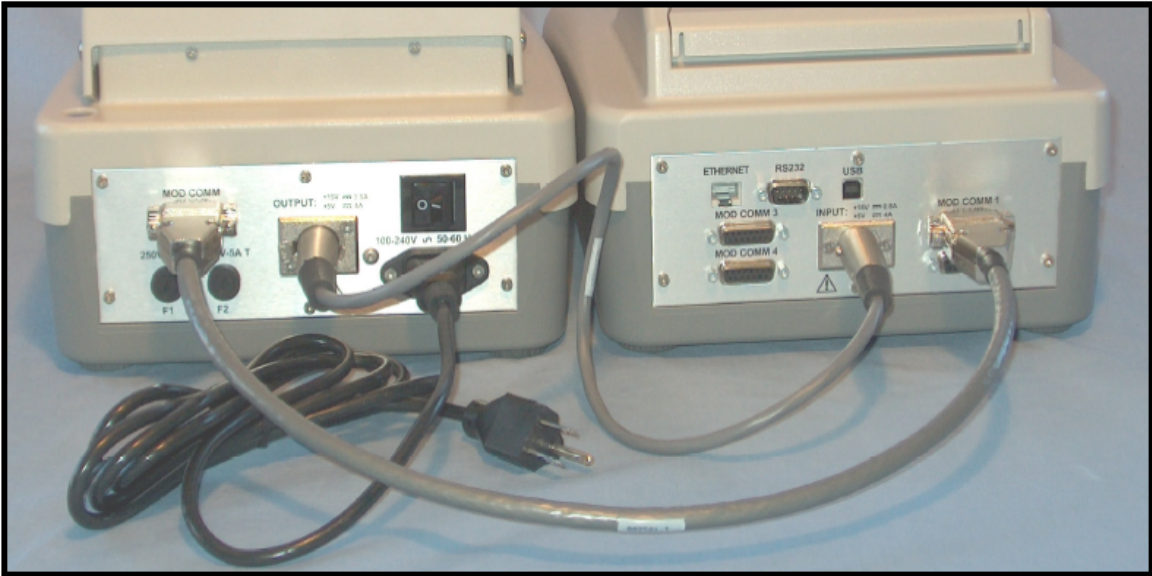
- 1 Pan shield
- 2 Pan support
- 3 Cylindrical quartz heating tubes
- 4 RTD temperature sensor
- 5 Heater adjustment tool connector
- 6 LED status lights
- 7 Service panel, heaters & RTD
- 8 Level vial
- 9 Adjustable leveling feet
- 10 Inter-module cable receptacle
- 11 Main power switch
- 12 Main cord receptacle
- 13 Fuse 1 (250 V~5A T)
- 14 Fuse 2 (250 V~5A T)
- 15 Module power cable receptacle

No. Designation

Control Module

- 16 Printer cavity
- 17 1/4 VGA display
- 18 Soft keys and display overlay
- 19 Keypad
- 20 Inter-module cable receptacle
- 21 Module power cable receptacle
- 22 Ethernet socket
- 23 USB socket

LMA100



Heating



MOD COMM

MOD COMM1

OUTPUT INPUT

Warm - up

15

Enter

LMA100P



Start/Stop

Enter

Print
Feed

Navigational keys

Soft keys

Print

Recall

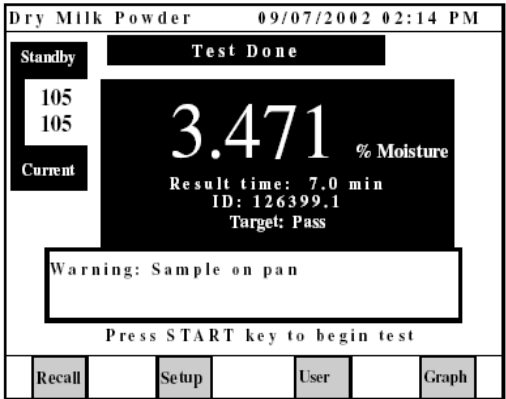
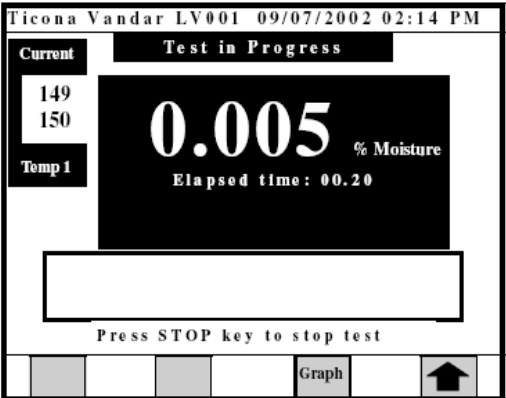
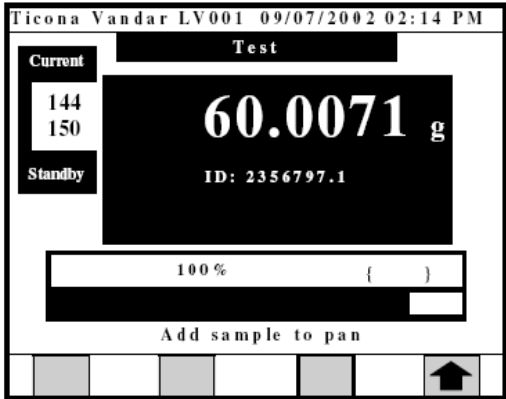
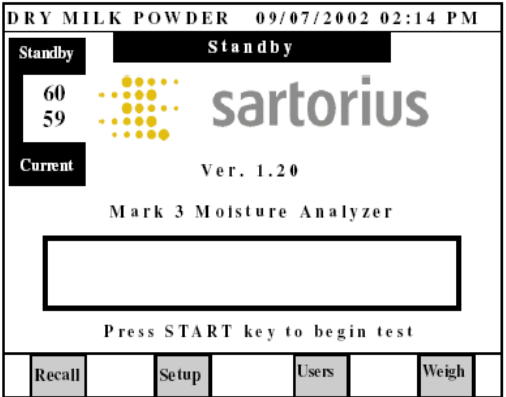
Setup

Users

Up arrow

Left arrow

LMA100P



Recall

가 Start/Stop

(, Zero .)

"Add sample to pan"

Beeper 가 ,

"Close heater hood"

"Sampling"

가

가

가

"Test Done"

*

tweezer

(

)