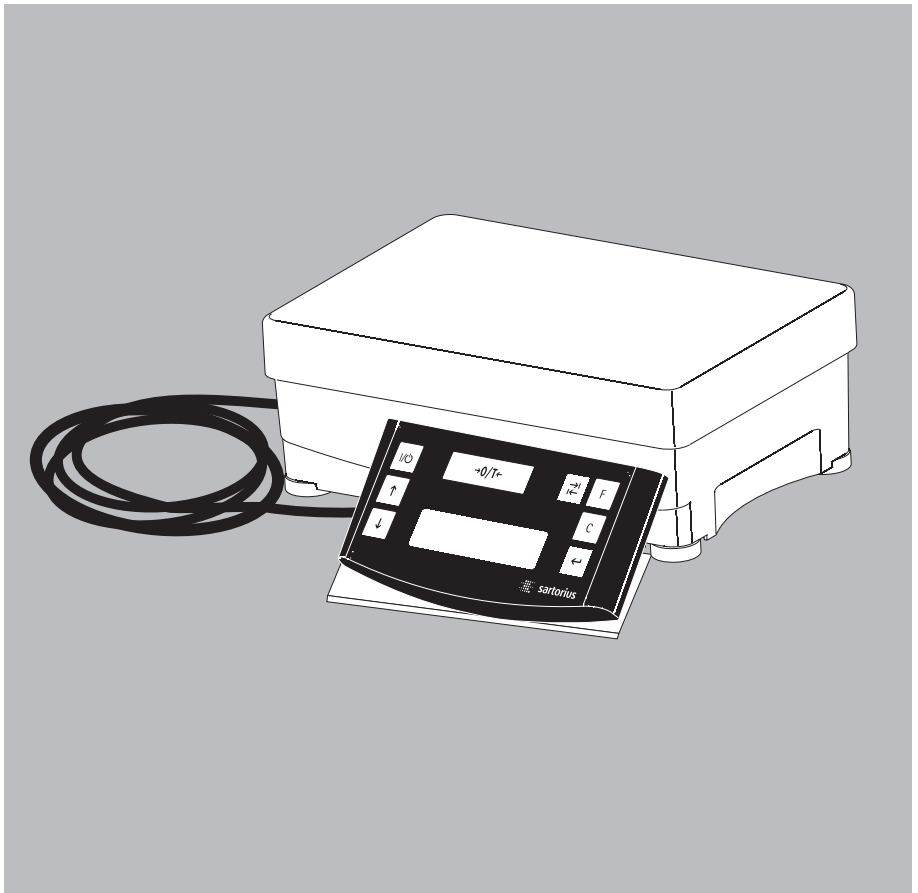
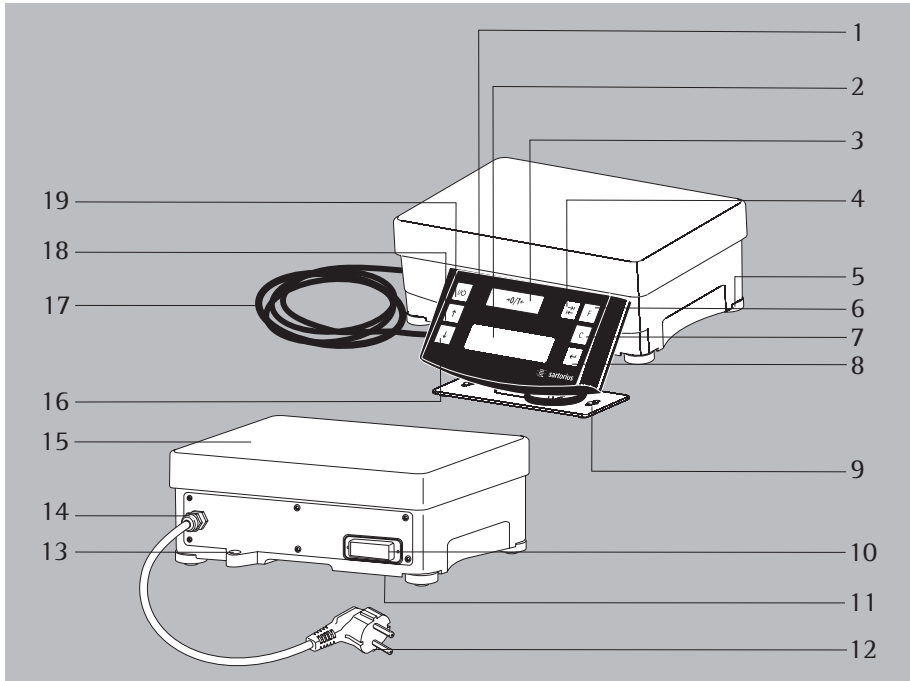


Installation Instructions**Sartorius PMA.Power
Model PMA35001**

Electronic Paint-mixing Scales



General View of the PMA35001



- | | | | |
|---|--|----|--|
| 1 | Display and control unit | 9 | Mounting bracket for display and control unit |
| 2 | Display | 10 | Interface (25-contact D-Sub socket) |
| 3 | Zero/tare key | 11 | Grounding terminal |
| 4 | Toggle key | 12 | Power cord with plug |
| | You can configure the PMA35001 menu to enable toggling between grams (g) and parts per pound (p) | | |
| 5 | Leveling foot | 13 | Level indicator |
| 6 | Factor key | 14 | Cable gland |
| 7 | Clear key and [REC] key for paint-mixing applications; depends on the menu setting | 15 | Load plate |
| 8 | [ENTER] key and [MEM] key for paint-mixing applications | 16 | 'Down' key |
| | | 17 | Power supply and data cable for the display and control unit |
| | | 18 | 'Up' key for paint-mixing applications |
| | | 19 | On/off (standby) key |

Contents

General View of the PMA35001	2
Contents	3
Intended Use	3
Warnings and Safety Precautions	4
Getting Started	5
Operation	7
Applications	8
Calibration/Adjustment	11
Operating Menu Configuration	12
Troubleshooting Guide	17
Care and Maintenance	18
Recycling	19
Interface Description	20
Specifications	21
Accessories	21
EU Declaration of Conformity	22
FCC Supplier's Declaration of Conformity	23

Symbols

The following symbols are used in these instructions:

- indicates required steps
- indicates steps required only under certain conditions
- > describes what happens after you have performed a particular step
- indicates an item in a list
- △ indicates a hazard

Intended Use

The PMA35001 scale is specially designed for use in paint-mixing applications. This scale can be operated through the keypad in stand-alone operation or using application software (such as a paint-mixing program from a paint manufacturer) installed on a connected PC.

If you wish to create your own application software, Sartorius can supply the required drivers for DOS or Windows operating systems.

Note:

- Make sure to read these operating instructions carefully before installing and operating your PMA35001 scale.
- As described in the manufacturer's declaration, type PMA35001 electronic scales (only with the power cord supplied by Sartorius, type H03 or comparable) may be used in flammable environments if the room is not at risk of inflammation due to excessive dust or fiber deposits. Paint-mixing rooms are generally free of accumulated dust or fibers.

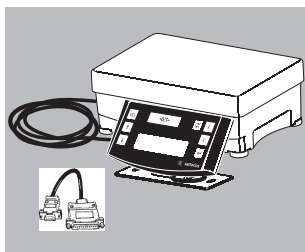
Warnings and Safety Precautions

The weighing instrument complies with the European Council Directives as well as international regulations and standards for electrical equipment, electromagnetic compatibility, and the stipulated safety requirements. Improper use or handling, however, can result in damage and/or injury. To prevent damage to the equipment, read these operating instructions carefully before using your scale.

- Keep these instructions in a safe place.
- Do not expose the weighing instrument to aggressive chemical vapors or to extreme temperatures, moisture, shocks, or vibration.
- Exposure to excessive electromagnetic interference can cause the readout value to change. Once the disturbance has ceased, the instrument can be used again in accordance with its intended purpose.
- Do not use electrical equipment out of doors.
- The scale may not be operated in hazardous areas or areas containing potentially explosive agents. If you use electrical equipment in installations and under ambient conditions subject to stricter safety standards than those described in the manual, you must comply with the provisions as specified in the applicable regulations for installation in your country.
- Make absolutely sure to disconnect the scale from power (unplug the power cord) before you connect or disconnect a peripheral device (printer or PC) to or from the interface port.
- If you use cables purchased from another manufacturer, check the pin assignments in the cable against those specified by Sartorius before connecting the cable to Sartorius equipment, and disconnect any wires that are assigned differently. The operator shall be solely responsible for any damage or injuries that occur when using cables not supplied by Sartorius. Connect only Sartorius accessories and options, as these are optimally designed for use with your PMA scale.
- The power connection must be made in accordance with the regulations applicable in your country. If you need assistance, contact your Sartorius dealer or the Sartorius Service Center. Any installation work that does not conform to the instructions in this manual will result in forfeiture of all claims under the manufacturer's warranty.
- If there is any indication that safe operation of the equipment is no longer warranted; for example, if the AC adapter is defective or shows sign of damage, disconnect the equipment from power and lock it in a safe place to ensure that it cannot be used for the time being.
- Disconnecting the grounding conductor is prohibited. Make sure the applicable accident prevention regulations are observed by all operating personnel.
- Always make sure the equipment is disconnected from AC power before performing any installation, cleaning, maintenance or repair work on the scale or AC adapter. If the equipment housing is opened by anyone other than persons authorized by Sartorius, all claims under the manufacturer's warranty are forfeited.
- The casing on all connecting cables, as well as the casing on wires inside the equipment housing, is made of PVC. Chemicals that corrode this material must be kept away from these cables.
- Do not leave the interface port uncovered. If the port is not used, or in preparation for shipping, replace the cap to protect the data interface from vapors, moisture and dust or dirt.
- Never use a hammer to close a paint can while it is still on the load plate, as this will damage the weighing system.

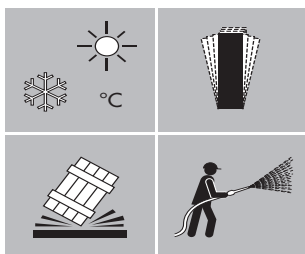
Getting Started

- Unpack the scale carefully.
- After unpacking the equipment, please check it immediately for any visible damage.



Equipment Supplied

- Scale
- Load plate
- Adapter cable



Setting up the Scale

Choose a suitable location where your scale will not be exposed to drafts, heat radiation, moisture or vibrations. Make sure to read these operating instructions carefully before connecting the scale to power.

- △ Make sure to observe the safety instructions.

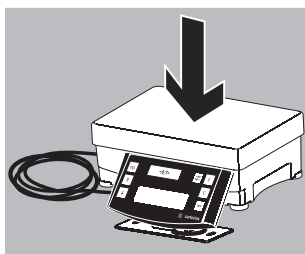
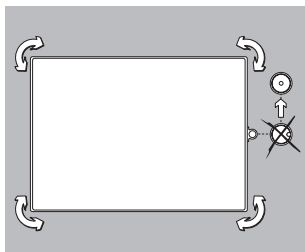
Leveling the Weighing Platform

Purpose:

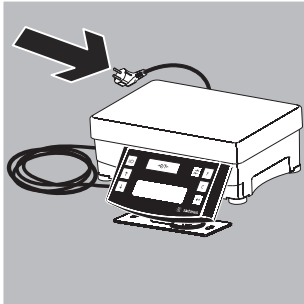
- To compensate for uneven areas at the place of installation
- To ensure that the equipment is placed in a perfectly horizontal position for consistently reproducible weighing results

Always level the weighing platform again any time after it has been moved to a different location.

- Level the weighing platform using the four leveling feet. Turn the feet until the air bubble is centered in the level indicator.
- Check to ensure that all leveling feet rest securely on the work surface.
- > Each of the leveling feet must support an equal load.
- > Adjusting the leveling feet:
 - To raise the weighing platform, extend the leveling feet (turn counterclockwise). To lower the weighing platform, retract the leveling feet (turn clockwise).

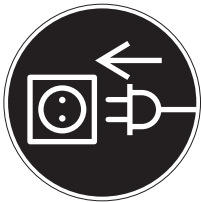


- Place the load plate on the scale.



Connecting the Scale to AC Power

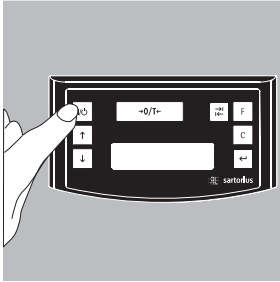
The display and control unit is powered through the scale's power cord and power PCB.




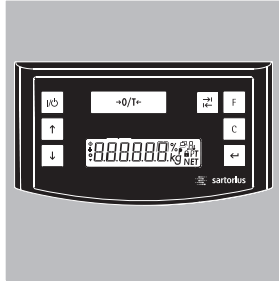
- Plug the power cord into a wall outlet (mains).

- ⚠ Make sure to observe the safety instructions. Make sure that the voltage rating printed on this unit is identical to your local line voltage. If the voltage specified on the label or the plug design of the power supply do not match the rating or standard you use, please contact your Sartorius office or dealer. Use only original Sartorius power cables. Use of cables from other manufacturers, even if these units have a registered approval rating from a national testing laboratory, requires the consent of a qualified electrician.
- If there is visible damage to the equipment or power cord, unplug the equipment and lock it in a secure place to ensure that it cannot be used for the time being.
 - Use only extension cords that meet the applicable standards and have a protective grounding conductor.
 - Disconnecting the grounding conductor is prohibited.

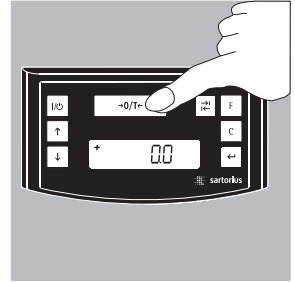
Operation

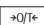


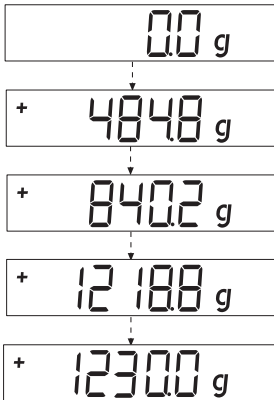
Press the  key (19) to switch on the scale.



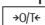
A self-test runs automatically, at the end of which the readout shows **0.0 g**.



If a different value is displayed, press  key to tare the scale.



Scale Functions

Center an empty paint can on the load plate. Press the  key (zero/tare; 3). The display shows **0.0 g**. Pour the first component of your formula into the can and read off the weight when the stability symbol (in this example, "g") is displayed. Add the other components up to the desired weight (formula). Remove the filled paint can from the load plate.



Never use a hammer to close a paint can while it is still on the load plate, as this will damage the weighing system.

Applications

Formulation (Calculation by a Factor)

The formulation program makes it easy to adjust your formulas for different final quantities of a paint-mixing recipe (for example, to make 250 ml of a 1-liter recipe). Press the **[F]** key (6) to select the desired factor. You can choose from the following: 0.5 1.0 1.5 2.0 2.5 3.0 3.5 4.0 4.5 5.0 10.0 15.0 20.0 25.0.

Press the **[↑]** key (18) (up) or **[↓]** key (16) (down) to modify the factor in 1.0 steps for factors of 5.0 or more, in 0.1 steps for factors of 1.0 to 5.0, or in 0.01 steps for factors of up to 1.0.

Note:

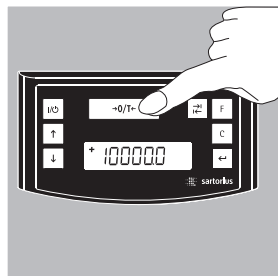
A flashing arrow **▼** on the display indicates that the weight value displayed is not valid in legal metrology (i.e., not legal for trade).

Example:

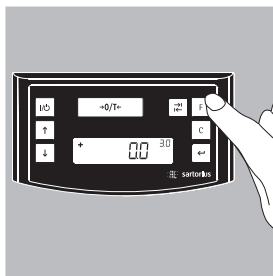
During formulation, the weight value is display in **g**. Let's suppose you want to weigh 3 kg of a basic formula that is for a total amount of 1 l. With the recalculation mode, you do not need to manually recalculate the individual components.

The basic formula for 1 liter is:

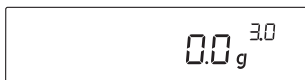
250 g first component
+ 250 g second component
+ 500 g third component
Total: 1000 g



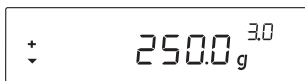
1. Place the empty paint can on the load plate and tare the scale.



2. Press the **[F]** key (6) repeatedly to set the factor to 3 for this example.



3. **3.0** next to the weight readout.



4. Slowly pour in the first component until the display shows **250 g**



5. Pour in the second component until **500 g** is displayed.



6. Pour in the last component **until 1000 g** is displayed



This concludes the recalculation example.

According to the display, exactly 1000 g was poured in, but the paint can actually contains 3 kg by weight in accordance with the factor you selected (3).

The procedure is the same for any other conversion factor.

Weighing Using the Recalculation Mode

Let's suppose that you poured in too much of one color component for a given formula (in this example, a 4-component recipe).

This example further assumes that you previously poured in all of the other amounts exactly according to each of the values you entered and saved by pressing the $\boxed{\leftarrow}$ key [MEM] (9). To correct for the overpour, press the $\boxed{\downarrow}$ key (16) to start the recalculation program; **C** flashes on the display. Use the $\boxed{\uparrow}$ key (18; up) or $\boxed{\downarrow}$ key (16; down) to correct the value so that it matches the amount called for in the formula. Then press $\boxed{\leftarrow}$ [MEM] (8) to have the scale calculate the amount to be added for each of the components that were already poured. The display shows the amounts required to correct the formula up to the point at which the overpour occurred. After the correction has been completed, you can continue filling the remaining components.

Note:

You can correct overpours as often as needed.

Keep in mind that the total quantity of paint at the conclusion of filling increases each time you correct a component. Press the $\boxed{\text{C}}$ key (7) to check how much the total quantity will be. The **C** stands for "Correction factor."

A flashing arrow \blacktriangledown on the display indicates that the weight value displayed is not valid in legal metrology (i.e., not legal for trade).

Example (Cumulative Weighing)

1. Center an empty paint can on the load plate.
+ 118.0 g

2. Press the \square key (zero/tare; 3).
0.0 g

3. Add first component
+ 50.0 g

4. Press the \square key [MEM] (8)
STO 01

5. Add the second component
+ 110.0 g

6. Press the \square key [MEM] (8)
STO 02

7. Add the third component
+ 203.0 g
The required amount of this component has been exceeded by 3 g (= overpour).

8. Press the \square key (16) to start the recalculation mode. A C ("Correction") flashes on the display.

9. Press the \square key (16) repeatedly to correct the value.
+ 200.0 g

10. Press the \square key [MEM] key (8)
COR 01

11. The readout shows the amount of the first component to be added, and "C1" is shown in the upper right-hand corner of the display.
- 1.7 g.

12. Add the first component until the readout shows 0.0 g

13. Press the \square key [MEM] key (8)
COR 02

14. The readout shows the amount of the second component to be added, and "C2" is displayed. - 2.0 g

15. Add until 0.0 is displayed.

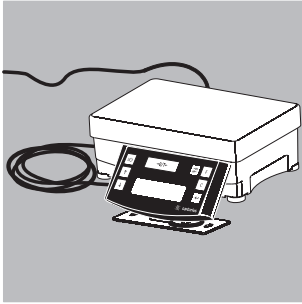
16. Press the \square key [MEM].
The scale returns to the formulation program and the "C" is no longer displayed.
+ 200.0 g.

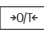
17. Press the \square key (7) [REC] to view the factor by which the total weight will exceed the original target defined in the recipe. (C = "Correction;" in this example, 1.03).
Total weight = original target \times correction factor)

18. Add remaining components
+ 1000.0 g

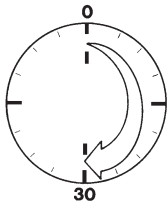
This concludes the recalculation example.

Calibration/Adjustment

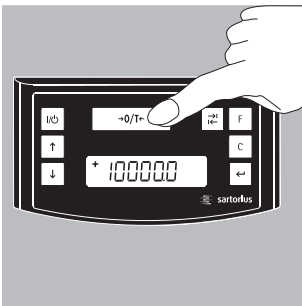


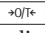
You can calibrate/adjust the scale by pressing the  key (3).

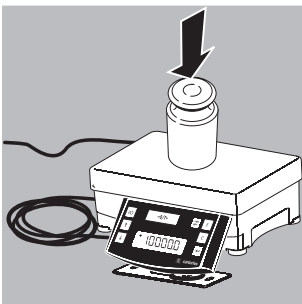
Calibration weight: 10,000 g; resolution: ± 0.016 g.



Always allow approximately 30 minutes for the scale to warm up after connecting it to AC power and before performing calibration/adjustment.



Press the  key (zero/tare) for 2 seconds; "10000" is displayed. Release the key.



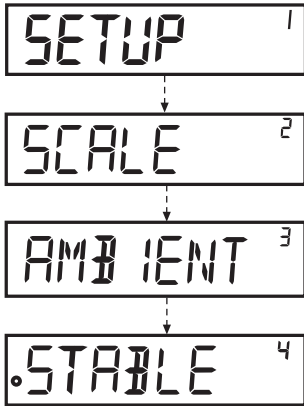
Center the calibration weight on the load plate (15). Adjustment is performed automatically. Remove the calibration weight.

Operating Menu Configuration

Accessing the Operating Menu

Example:

Adapting the Scale to Ambient Conditions



- Hold down the key [ENTER] for approximately 2 seconds; "SETUP" is displayed.
- Use the keys to select the desired menu item on this level.
- Press the key [Enter] to select the next menu level.
- The second menu level is displayed. Use the keys to select the desired menu item on this level.
- Press the key [ENTER] to select the next menu level.
- The third menu level is displayed. Use the keys to select the desired menu item on this level.
- Press the key [ENTER] to select the next menu level.
- The fourth menu level is displayed. Use the keys to select the desired menu item on this level.


This concludes the example.

- Press the key [ENTER]; is displayed, indicating that this item is set.
- Press (Clear) repeatedly to exit the menu.

Note:

Contact your local Sartorius office for a detailed list of the menu codes.

Configuring the Main Menu Settings

- Hold down the  key [ENTER] for approximately 2 seconds; "SETUP" is displayed.

Level 1






SETUP

Language Setting

Level 1 Level 2

LANGUAGE

- o GERMAN
- ENGLISH
- FRENCH
- ITALIAN
- etc.

- Press the  key to select LANGUAGE
- Press the  key [ENTER]
- Press the  keys to select a language
- Press the  key [ENTER];
o indicates the active setting
- Press the  key repeatedly to exit the menu

Default Unit: Grams or Parts per Pound

The default settings is active when the scale is switched on; defined under "SETUP > SCALE > UNIT":



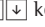


Level 1 Level 2 Level 3 Level 4

SETUP


SCALE

UNIT

- o GRAMS
- PT./P.D.

- Press the  key [ENTER]
- Press the  key [ENTER]
- Use the  keys to select the desired setting; e.g., "UNIT"
- Press the  key [ENTER]
- Use the  keys to select the desired unit; e.g., "GRAMS"

Activating the Toggle Key

When the toggle key  (4) is active, you can configure it to toggle the weight unit between grams and parts per pound. The unit is toggled when the key is pressed.

Level 1 Level 2 Level 3 Level 4




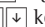




SETUP

APPLICATION

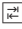
TOGGLE

OFF











- o ON

- Press the  key [ENTER]
- Use the  keys to select "APPLICATION"
- Press the  key [ENTER]
- Use the  keys to select "TOGGLE"
- Press the  key [ENTER]
- Use the  keys to select "ON"
- Press the  key [ENTER]; o indicates the active setting
- Press the  key (Clear) repeatedly to exit the menu.

Configuring the Toggle-key Function

Pressing the toggle key  (4) toggles the scale between the default unit (defined under SETUP > SCALE > UNIT; see page 13) and the unit defined as follows under SETUP > APPLICATION > UNIT.












Level 1	Level 2	Level 3	Level 4
SETUP			
	APPLICATION		
		UNIT	
			o PT./PD.
			GRAMS

- Press the  key [ENTER]
- Use the   keys to select “APPLICATION”
- Press the  key [ENTER], press  to select “UNIT” and press  to confirm.
- Use the   keys to select the desired unit; e.g., “PT./PD.”
- Press the  key [ENTER]; o indicates the active setting.
- Press the  key (Clear)

Activating the “LOCK” Function

The LOCK function protects the scale from unauthorized use. When this function is active, the scale readout shows weight values only when there is active communication between the scale and a PC. If communication is interrupted, the readout goes blank and the display shows a padlock symbol. Activation of the LOCK function is configured under “EXTRAS.”

Level 1	Level 2	Level 3	Level 4
SETUP			
	EXTRAS		
		LOCK	
			OFF
			o ON

- Press the  key [ENTER]
- Use the   keys to select “EXTRAS”
- Press the  key [ENTER]
- Use the   keys to select “LOCK”
- Press the  key [ENTER]
- Use the   keys to select ON and press  to confirm
- Press the  key (Clear) repeatedly to exit the menu.

Configuring Password Protection

In addition to the LOCK function, you can configure password protection for additional security. With this feature, the LOCK function can be deactivated only by entering the password you configure. The password is numeric and can have up to 6 digits. Use the $\uparrow\downarrow$ keys to select the digits (0 through 9) for your password. The password is hidden on the readout; only dashes (“-----”) are shown. The first dash flashes to prompt input. Press the $\uparrow\downarrow$ keys as needed to select the desired digit (0 to 9) and then press the \leftarrow key [ENTER] key. The digit is stored and the second dash flashes on the display. Repeat the input procedure as described for the first digit. To store a space as a character in the password, press the \leftarrow key [ENTER] while the corresponding dash is flashing. Once all 6 characters have been stored, press the \leftarrow key [ENTER] to store the password.

Note:

Keep a copy of your password in a safe place.

The LOCK function can be deactivated only with this password.

Level 1 Level 2 Level 3

INPUT

PASSWORD

PW.NEW

- Use the $\uparrow\downarrow$ keys to select “INPUT”
- Press the \leftarrow key [ENTER]
- Press the \leftarrow key [ENTER]
- Use the $\uparrow\downarrow$ keys to select “PW.NEW”
- Enter the desired password and press the \leftarrow key [ENTER]
- Press the \leftarrow key [CLEAR] repeatedly to exit the menu

Changing the Password

To define a new password, the existing password must be entered first in the SETUP menu under “PASSWORD.” “PW.OLD” prompts this input. Once the old password is entered, the “PW.NEW” prompt is shown automatically. Enter the new password or press \leftarrow at each position to enter spaces. The display shows spaces.

Note:

Entering 6 spaces deletes the password, which deactivates the password function.

Level 1 Level 2 Level 3

INPUT

PASSWORD

PW.OLD

PW.NEW

- Use the $\uparrow\downarrow$ keys to select “INPUT”
- Press the \leftarrow key [ENTER]
- Press the \leftarrow key [ENTER]
- Enter the existing (old) password
- > After the old password have been entered correctly, “PW.NEW” is displayed.
- Enter the desired password and press the \leftarrow key [ENTER]
- Press the \leftarrow key [CLEAR] to return to the menu

Configuring Text Length ("LONG" or "SHORT")

You can define the length of the operator guidance texts shown on the display.

Level 1	Level 2	Level 3	Level 4
SETUP			
	EXTRAS		
		TEXTS	
			LONG
			o SHORT

- Press the key [ENTER]
- Use the keys to select "EXTRAS"
- Press the key [ENTER]
- Use the keys to select "TEXTS"
- Press the key [ENTER]
- Use the keys to select "SHORT" and press to confirm
- Press the key (Clear) repeatedly to exit the menu

Resetting the Scale: "RESET"

You can restore the factory settings in the scale.

Note:

If you have activated the password function, this feature is password-protected.

Level 1	Level 2	Level 3	Level 4
SETUP			
	RESET		
		MENU	
			YES
			o NO

- Press the key [ENTER]
- Use the keys to select "REST"
- Press the key [ENTER]
- Use the keys to select "MENU"
- Press the key [ENTER]
- Use the keys to select "YES"
- Press the key. Factory settings are restored. "MENU" is displayed
- Press the key (Clear) repeatedly to exit the menu

Code Settings

Select the "CODES" menu item to have menu items identified by numeric codes rather than texts.


Level 1	Level 2
LANGUAGE	
	GERMAN
	etc.
	o CODES

- Press the key to select LANGUAGE
- Press the key [ENTER]
- Use the keys to select "CODES"
- Press the key [ENTER]; **o** indicates the active setting
- Press the key (Clear) repeatedly to exit the menu

Note:

Contact your local Sartorius office for a detailed list of the menu codes.

Troubleshooting Guide

Problem	Cause	Solution
No segments appear on the display	- No AC power is available	- Check the AC power supply
The weight readout shows "LOW"	- No load plate on the scale	- Position the load plate
The weight readout shows "HIGH"	- Weighing capacity exceeded	- Unload the scale
The weight readout changes constantly	- Unstable environment - Too much vibration, or the scale is exposed to a draft	- Change the place of installation - Access the menu to select the appropriate code for the ambient conditions (see "Operating Menu Configuration")
The weight readout is obviously wrong	- The sample is not stable - Scale not tared before weighing	- Tare the scale before weighing
No weight value is shown and the padlock symbol is displayed 	- Communication between scale and PC has been interrupted, activating the LOCK function	- Access the menu settings to deactivate the LOCK function - Check the connection

Care and Maintenance

Cleaning

- △ Never use concentrated acids, alkali solutions or pure alcohol to clean the equipment.
- △ Do not allow liquids to penetrate the equipment housing.
- Use a brush or a soft, dry, lint-free cloth to clean the scale.

Cleaning Stainless Steel Surfaces

Clean all stainless steel parts regularly. Remove the stainless steel load plate and thoroughly clean it separately. Use a damp cloth or sponge to clean stainless steel parts on the scale. You can use any commercially available household cleaning agent that is suitable for use on stainless steel. Clean stainless steel surfaces by wiping them down. Then clean the load plate thoroughly, making sure to remove all residues. Afterwards, wipe down stainless steel parts again using a clean, damp cloth or sponge and allow the equipment to dry. If desired, you can apply oil to the cleaned surfaces as additional protection.

- △ Solvents are permitted only for cleaning stainless steel parts.

Corrosive Environment

- Remove all traces of corrosive substances from the scale on a regular basis.

Storage and Shipping Conditions

- The packaging used for shipping your Sartorius equipment is optimally designed to prevent damage during transport. It is a good idea to save the box and all parts of the packaging for future storage or shipment of the equipment. Only the original packaging provides the best protection for shipment.
- Allowable storage temperature: -20 °C to $+75\text{ °C}$ (-4 °F to $+167\text{ °F}$)
- Allowable humidity during storage: up to 90%
- Please refer to the information under “Safety Inspection” below.

Safety Inspection

Safe operation of the scale is no longer ensured in the following cases:

- If there is visible damage to the connecting cable
- If the device no longer functions properly
- If the equipment has been stored for a relatively long period under unfavorable conditions
- If the equipment has been subjected to rough handling during shipment
- Make sure all warnings and safety precautions are observed, and notify your nearest Sartorius Service Center or the International Technical Support Unit based in Goettingen, Germany, if any of the above occurs. Maintenance and repair work may be performed only by authorized Sartorius service technicians who have access to the required maintenance manuals and have received the necessary training.

- △ Only authorized service technicians may open the equipment and perform maintenance work; this ensures that operation of the equipment is safe and trouble-free and the warranty remains in effect.

Recycling



The packaging is made from environmentally-friendly materials that can be used as secondary raw materials. If you no longer need this packaging, bring it to your local recycling and waste

disposal facility according to the regulations applicable in your country. In Germany, you can dispose of this material using the VfW dual system (contract number D-59101-2009-1129). The equipment, including accessories and batteries, must not be disposed of in general household waste, and must be recycled similar to electrical and electronic devices. For further information about disposal and recycling options, please contact your local service staff. The partners listed on the following website can be used for disposals within the EU:

- 1) Go to <http://www.sartorius.com>.
- 2) Select the summary under “Service.”
- 3) Then select “Information on Disposal.”
- 4) Addresses for local Sartorius disposal contacts can be found in the PDF files given on this webpage.



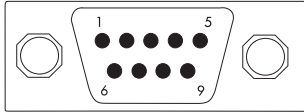
Sartorius will not take back equipment contaminated with hazardous materials (ABC contamination) either for repair or disposal.

Insert heading: “Service Address for Disposal”

Please refer to our website (www.sartorius.com) or contact the Sartorius Service Department for more detailed information regarding repair service addresses or the disposal of your device.

Interface Description

Connector on adapter cable



Pin Assignments

9-contact female connector:

Pin 2: (RxD) Receive data

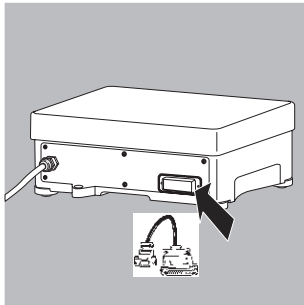
Pin 3: (TxD) Transmit data

Pin 4: (DTR) Data terminal ready

Pin 5: (GND) Ground

Pin 6: BPI bridge

Pin 8: Clear to send (CTS)



Installing the Adapter Cable

The adapter cable is required for operation of the scale with a PC. If you wish to create your own application software, Sartorius can supply the required drivers for DOS or Windows operating systems.

- △ Make sure to observe the safety instructions.
- Remove the protective cover from the interface port. Keep the protective cover in a safe place.
- Replace the interface cover when storing or shipping the scale.
- Plug in the adapter cable and tighten the retaining screws

Specifications

Model	PMA35001	
Weighing capacity	g	35000
Readability	g	0.1
Tare range (subtractive)	g	-35000
Linearity	g	<±0.2
Stability range (configured in operating menu)	digits	0.25 to 4
Humidity class	F	Anti-condensation
Ambient operating temperature range	°C	0 to +40 (+32 °F to +104 °F)
Load plate dimensions	mm	350 × 240
Scale housing (W × D × H)	mm	350 × 243 × 132.5
Net weight, approximate	kg	11.4
Calibration weight	kg	10; class E2 or better
Power consumption	VA	Average: 8; maximum: 16
Interface port	RS-232	
- Format	7-bit ASCII, 1 start bit, 1 or 2 stop bits	
- Parity	Even, odd, none	
- Transmission rates:	1,200 to 38,400 bit/s	
- Handshake mode	Software or hardware	

Accessories

Product	Order No.
Dust cover	YDC01PMA
RS-232 data cable, SBI (2 m)	YCC01-0027M2
RS-232 data cable, BPI (2 m)	YCC01-0028M2
RS-232 data cable, BPI (20 m)	YCC01-0028M20
USB/RS-232 data cable (SBI) (1.8 m)	YCO12
USB/RS-232 data cable (BPI) (1.8 m)	YCO13

EU Declaration of Conformity



CE EU-Konformitätserklärung EU Declaration of Conformity

Hersteller
Manufacturer Sartorius Lab Instruments GmbH & Co. KG
37070 Goettingen, Germany

erklärt in alleiniger Verantwortung, dass das Betriebsmittel
declares under sole responsibility that the equipment

Geräteart
Device type Hochlastige Farbmischwaage
High-capacity paint mixing scale

Modell
Model PMA35001

in der von uns in Verkehr gebrachten Ausführung allen einschlägigen Bestimmungen der folgenden Europäischen Richtlinien - einschließlich deren zum Zeitpunkt der Erklärung geltenden Änderungen - entsprecht und die anwendbaren Anforderungen folgender harmonisierter Europäischer Normen erfüllt:
in the form as delivered fulfils all the relevant provisions of the following European Directives - including any amendments valid at the time this declaration was signed - and meets the applicable requirements of the harmonized European Standards listed below:

2014/30/EU Elektromagnetische Verträglichkeit
Electromagnetic compatibility
EN 61326-1:2013

2014/35/EU Elektrische Betriebsmittel zur Verwendung innerhalb bestimmter Spannungsgrenzen
Electrical equipment designed for use within certain voltage limits
EN 61010-1:2010

2011/65/EU Beschränkung der Verwendung bestimmter gefährlicher Stoffe in Elektro- und Elektronikgeräten (RoHS)
Restriction of the use of certain hazardous substances in electrical and electronic equipment (RoHS)
EN 50581:2012

Jahreszahl der CE-Kennzeichenvergabe | *Year of the CE mark assignment:* 16

Sartorius Lab Instruments GmbH & Co. KG
Goettingen, 2016-04-20

Dr. Reinhard Baumfalk
Vice President R&D

Dr. Dieter Klausgrete
Head of International Certification Management

Diese Erklärung bescheinigt die Übereinstimmung mit den genannten EU-Richtlinien, ist jedoch keine Zusicherung von Eigenschaften. Bei einer mit uns nicht abgestimmten Änderung des Produktes verliert diese Erklärung ihre Gültigkeit. Die Sicherheitshinweise der zugehörigen Produktdokumentation sind zu beachten.

This declaration certifies conformity with the above mentioned EU Directives, but does not guarantee product attributes. Unauthorised product modifications make this declaration invalid. The safety information in the associated product documentation must be observed.

FCC Supplier's Declaration of Conformity

FCC Supplier's Declaration of Conformity



Device type Electronically laboratory balance

Model PMA35001

Party issuing Supplier's Declaration of Conformity / Responsible Party – U.S. Contact Information

Sartorius Corporation
5 Orville Dr Suite 200
11716 Bohemia, NY
USA
Telephone: +1.631.254.4249

FCC Compliance Statement

This device complies with Part 15 of the FCC Rules. Operation is subject to the following two conditions: (1) This device may not cause harmful interference, and (2) this device must accept any interference received, including interference that may cause undesired operation.

Information to the user

Note: This equipment has been tested and found to comply with the limits for a **class B** digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference in a residential installation. This equipment generates, uses and can radiate radio frequency energy and if not installed and used in accordance with the instructions, may cause harmful interference to radio communications. However, there is no guarantee that interference will not occur in a particular installation. If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try to correct the interference by one or more of the following measures:

- Reorient or relocate the receiving antenna.
- Increase the separation between the equipment and receiver.
- Connect the equipment into an outlet on a circuit different from that to which the receiver is connected.
- Consult the dealer or an experienced radio/TV technician for help.

Connections between the device and peripherals must be made using shielded cables in order to maintain compliance with FCC radio frequency emission limits.

Any modifications made to this device that are not approved by Sartorius may void the authority granted to the user by the FCC to operate this equipment.

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen, Germany

Phone: +49.551.308.0
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice. Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote the other gender as well.

Copyright notice:

This instruction manual, including all of its components, is protected by copyright. Any use beyond the limits of the copyright law is not permitted without our approval. This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated:
10 | 2018