



**Syntax Description**

**Sartorius**

**SICS Interface**



# Contents

1. Intended Use	2
2. Settings for Cubis MSA and MSU Balances	5
3. Syntax of SICS Commands	6
4. Syntax of SICS Responses	7
5. Description of SICS Commands	8

## 1. Intended Use

SICS (Standard Interface Common Set) consists of commands that are used to control Cubis MSA and MSU balances through interfaces.

The command scope covers functions for reading measurement data and triggering weighing commands (e.g., tare), and functions for remote access to the user interface (writing text in displays, activating keys, reading out key functions).

Use a program such as the SartoTerminal from Sartorius to communicate with the balance.

## The Following Commands Are Available:

### ► Level 0, Version 2.3x

- @ - Reset all SICS commands
- I0 - List all available commands
- I1 - Send information about the level and its versions
- I2 - Query balance model
- I3 - Query software version of balance (BAC)
- I4 - Query balance serial number
- I5 - Query software version of display (APC)
- S - Send weight value at stability
- SI - Send weight value without stability
- SIR - Send automatic weight values at and without stability
- Z - Zero the balance at stability
- ZI - Zero the balance without stability

### ► Level 1, Version 2.2x

- T - Tare the balance at stability
- TI - Tare the balance without stability
- D - Write text in display
- DW - Delete text from display
- K - Key control
- TA - Query and allocate tare memory
- TAC - Delete tare memory
- SR - Send weight value if there is a weight change

### ► Level 2

- SU - Send weight value at stability with current weight unit (with motorized draft shield)
- WS - Query door position of motorized draft shield; open or close door
- PWR - Turn balance on/off (standby)
- I10 - Query/set the balance ID
- I11 - Query balance type
- I14 - Query balance components
- M01 - Query/set application filter
- M02 - Query/set filter adjustment
- M03 - Query/set automatic zeroing
- M04 - Query/set I/O inputs
- M07 - Query/activate/deactivate automatic draft shield
- M12 - Acoustic signal (beep)
- M13 - Activate/deactivate touchscreen softkeys
- M24 - Query/activate/deactivate "Print" key; print stable or unstable weight values
- M39 - Query/activate bar graph display in working environment
- P100 - Send text line to printer
- C1 - Execute calibration/adjustment (as set in menu)

## ► Remote Control

- P112 – Write text in selected line on display
- P113 – Delete text from selected line in display
- P114 – Overwrite task or user names
- P120 – Turn off bar graph in checkweigher
- P121 – Turn on bar graph in checkweigher
- RM20 – Activate/deactivate user input
- RM30 – Assign new function to softkeys
- RM32 – Assign new order to softkeys
- RM34 – Create a dynamic parameter
- RM35 – Immediately change softkey designations
- RM36 – Assign/query function for multiple softkey lines
- RM37 – Prepare preset softkey designations for display
- RM38 – Activate RM36-assigned softkey lines
- RM39 – Activate/deactivate RM30-assigned softkey functions
- RM44 – Query/set input with barcode scanner
- RM48 – Change order of standard keys
- RM49 – Activate/deactivate info text
- RM51 – Activate/deactivate selection window
- RM52 – Define properties for a window with info text
- RM53 – Activate/deactivate window with info text
- RM54 – Activate/deactivate window with info

## ► Additional Sartorius Commands

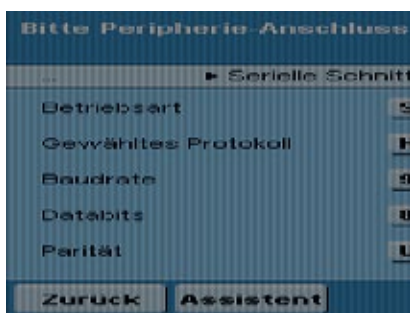
- SA – Send weight value at stability and store in Alibi memory (with optional ID)
- CMD – Execute application command
- PAR – Query parameter
- MN36 – Assign a function to several menus
- MN38 – Display or hide MN36-assigned menus
- TX36 – Assign text to several text pages
- TX37 – Overwrite a line on a text page
- TX38 – Activate/deactivate TX38-assigned text pages

## 2. Settings for Cubis MSA and MSU Balances

### SICS Commands Can Be Used via the Following Interfaces:

- ▶ Serial (RS-232)
- ▶ USB
- ▶ Bluetooth
- ▶ Ethernet

Go to Menu > Device parameters > Configure ports. The interface must be set to "SICS" mode:



### Additional Settings:

- ▶ Handshake:
  - No handshake
  - Software handshake: XON/OFF
  - Hardware handshake: RTS/CTS
- ▶ Baud rate:
  - 300 baud
  - 600 baud
  - 1200 baud
  - 2400 baud
  - 4800 baud
  - 9600 baud
  - 19200 baud
  - 38400 baud
  - 57600 baud
  - 115200 baud
- ▶ Data bits:
  - 7 data bits
  - 8 data bits
- ▶ Parity:
  - No
  - Odd
  - Even
- ▶ Stop bits:
  - 1 stop bit
  - 2 stop bits
- ▶ Log data:
  - Switch off
  - Switch on

### 3. Syntax of SICS Commands

An SICS command consists of an identifier (ID) and optional parameters.

ID [Parameter<sub>1</sub>] [Parameter<sub>2</sub>] [Parameter<sub>3</sub>] ...[Parameter<sub>n</sub>] <CR><LF>

The identifiers (IDs) consist of ASCII characters and are written only in capital letters. The parameters must be separated with a space. If a text parameter contains at least one space, this parameter must be put in quotation marks. Each command must end with a carriage return and line feed (#0D#0A or <CR><LF>).

Example: Write text in the third line of the display

P112 3 "Place the second component on the balance"<CR><LF>

### 4. Syntax of SICS Responses

The balance sends a response to each SICS command.

The responses may contain one or more weight values and/or text.

ID Status [Parameter<sub>1</sub>] [Parameter<sub>2</sub>] [Parameter<sub>3</sub>] ... [Parameter<sub>n</sub>] <CR><LF>

"ID" corresponds to the command identifier that was sent to the balance (exceptions: S instead of SI, SIR, SR, and I4 instead of @)  
 "Status" provides feedback on how the command was executed:

- ▶ A – Command executed; no further response will be sent
- ▶ B – Command executed; a further response will be sent  
(example: I0 - List all commands)
- ▶ C – Key pressed, function was not executed, and response sent (e.g., K 3)
- ▶ D – Weight value without stability
- ▶ I – Command could not be executed (e.g., because balance is already tared)
- ▶ L – Command has a syntax error and could not be executed
- ▶ S – Weight value at stability
- ▶ + – Weight value too high
- ▶ - – Weight value too low

The return parameters are weight values with corresponding weight units, numerical parameters, or text.



## I1 – Send information about the level and its versions

Syntax:

Command: I1<CR><LF>

Response: I1 A "P<sub>1</sub>" "P<sub>2</sub>" "P<sub>3</sub>" "P<sub>4</sub>" "P<sub>5</sub>"<CR><LF>

P<sub>1</sub>: "01" SICS Level 0 and SICS Level 1 available

P<sub>2</sub>: Version from Level 0

P<sub>3</sub>: Version from Level 1

P<sub>4</sub> and P<sub>5</sub>: Empty, as Level 2 and Level 3 are not available

or

I1 I<CR><LF>

I: Command cannot currently be executed

Example:

Command: I1

Response: I1 A "01" "2.30" "2.20" "" ""

## I2 – Query balance model

Syntax:

Command: I2<CR><LF>

Response: I2 A "P<sub>1</sub>"<CR><LF>

P<sub>1</sub>: Balance model description

or

I2 I<CR><LF>

I: Command cannot currently be executed

Example:

Command: I2

Response: I2 A "MSA3203P"

## I3 – Query software version of balance (BAC)

Syntax:

Command: I3<CR><LF>

Response: I3 A "P<sub>1</sub>"<CR><LF>

P<sub>1</sub>: Software version of balance

or

I3 I<CR><LF>

I: Command cannot currently be executed

Example:

Command: I3

Response: I3 A "00-39-05"

## I4 – Query balance serial number

Syntax:

Command: I4<CR><LF>

Response: I4 A "P<sub>1</sub>"<CR><LF>

P<sub>1</sub>: Balance serial number

or

I4 I<CR><LF>

I: Command cannot currently be executed

Example:

Command: I4

Response: I4 A "23201202"



## I5 – Query software version of display (APC)

Syntax:

Command: I5<CR><LF>

Response: I5 A "P<sub>1</sub>"<CR><LF>

or

I5 I<CR><LF>

P<sub>1</sub>: Software version of display

I: Command cannot currently be executed

Example:

Command: I5

Response: I5 A "01-60-04"

## S – Send weight value at stability

If the balance has a motorized draft shield and is set to automatic draft shield, the draft shield is shut first and then the weight value is sent at stability. The draft shield may open after this command is executed, depending on the motorized draft shield setting (see command M07).

Syntax:

Command: S<CR><LF>

Response: S S w<sub>1</sub> u<sub>1</sub><CR><LF>

S +<CR><LF>

S -<CR><LF>

or

S I<CR><LF>

S: Weight value at stability

w<sub>1</sub>: Weight value

u<sub>1</sub>: Unit of weight

+: Balance overload

-: Balance underload

I: Command cannot currently be executed

Example:

Command: S

Response: S S 99.528 g

Current weight value at stability is 99.528 g

## SI – Send weight value without stability

Syntax:

Command: SI<CR><LF>

Response: S S w<sub>1</sub> u<sub>1</sub><CR><LF>

S D w<sub>1</sub> u<sub>1</sub><CR><LF>

S +<CR><LF>

S -<CR><LF>

or

S I<CR><LF>

S: Weight value at stability

D: Weight value without stability

w<sub>1</sub>: Weight value

u<sub>1</sub>: Unit of weight

+: Balance overload

-: Balance underload

I: Command cannot currently be executed

Example:

Command: SI

Response: S S 99.528 g

S D 362.359 g

Current weight value at stability is 99.528 g

Current weight value without stability is 362.359 g

## SIR – Send automatic weight values at and without stability

The SIR command is used by the balance to query weight values cyclically.

The frequency for weight value queries is set as part of the task ("Print Settings" menu, with parameter "Interval For Automatic Printout"): once, twice, five times, or ten times per second.

If no task is activated, set this parameter in Menu > Configure device > Configure data output

SIR is terminated with the commands @, S, SI, and SR or by turning the balance off.

Syntax:

Command: SIR<CR><LF>

Response: S S w<sub>1</sub> u<sub>1</sub><CR><LF>

S D w<sub>1</sub> u<sub>1</sub><CR><LF>

S +<CR><LF>

S -<CR><LF>

or

S I<CR><LF>

S: Weight value at stability

D: Weight value without stability

w<sub>1</sub>: Weight value

u<sub>1</sub>: Unit of weight

+: Balance overload

-: Balance underload

I: Command cannot currently be executed

Example:

Command: SIR

Response: S D 594.821 g

S D 228.896 g

S D 228.885 g

S S 228.890 g

## Z – Zero the balance at stability

If the Z command is sent to the balance, the balance waits for stability to be achieved and is then zeroed.

Syntax:

Command: Z<CR><LF>

Response: Z A<CR><LF>

or

Z I<CR><LF>

A: Balance zeroed

I: Command cannot currently be executed

Example:

Command: Z

Response: Z A

Balance zeroed

## ZI – Zero the balance without stability

If the ZI command is sent to the balance, the balance is zeroed immediately (even without stability).

Syntax:

Command: ZI<CR><LF>

Response: ZI D<CR><LF>

or

ZI I<CR><LF>

D: Balance zeroed without stability

I: Command cannot currently be executed

Example:

Command: ZI

Response: ZI D

Balance zeroed

## Level 1 Version 2.2x

### T – Tare the balance at stability

If the T command is sent to the balance, the balance waits for stability to be achieved and is then tared. The tare memory is overwritten with the new tare value. If the current weight value is less than zero, the balance cannot be tared (but can be zeroed). If the "Second tare" application is active, tare memory T1 is overwritten with this command. If the "Second tare" application is not active, the balance tare memory is overwritten with this command. In this case, it is not possible to preset the balance tare memory with an entered (non-weighed) weight value.

Syntax:

Command: T<CR><LF>

Response: T S  $w_1$   $u_1$ <CR><LF>

S: Balance tared at stability

$w_1$ : Weight value

$u_1$ : Unit of weight

or

T I<CR><LF>

I: Command cannot currently be executed

Example:

Command: T

Response: T S 29.817 g

Balance tared, and weight value 29.817 g recorded in the tare memory

### TI – Tare the balance even without stability

The balance is tared immediately. The tare memory is overwritten with the new tare value. If the current weight value is less than zero, the balance cannot be tared (but can be zeroed).

If the "Second tare" application is active, tare memory T1 is overwritten with this command.

If the "Second tare" application is not active, the balance tare memory is overwritten with this command.

Syntax:

Command: TI<CR><LF>

Response: TI D  $w_1$   $u_1$ <CR><LF>

D: Balance tared without stability

$w_1$ : Weight value

$u_1$ : Unit of weight

or

TI I<CR><LF>

I: Command cannot currently be executed

Example:

Command: TI

Response: TI D 29.817 g

Balance tared, and weight value 29.817 g recorded in the tare memory

## D – Write text in display

Text is written in the working environment; the weight value is not overwritten. If the text is too long and cannot be shown entirely in one line, it is cut off at the end of the line.

Syntax:

Command: D "Text"<CR><LF>

Response: D A<CR><LF>

or

D I<CR><LF>

D L<CR><LF>

A: Text appears in the working environment

I: Command cannot currently be executed

L: Syntax error (check if the text is in quotation marks and that there is a space between D and "Text")

Example:

Command: D "Place the third component on the balance"

Response: D A

The text "Place the third component on the balance" appears in the working environment

## DW – Delete text from display

Text written in the working environment with the D command is deleted again. The original text – the corresponding task – appears again in the working environment.

Syntax:

Command: DW<CR><LF>

Response: DW A<CR><LF>

or

D I<CR><LF>

A: Text in the working environment deleted

I: Command cannot currently be executed

Example:

Command: DW

Response: DW A

Text deleted, and original text displayed again

## K – Key control

The K command can be used to lock the keys and/or query which key was pressed on the balance. Key control can be reset with commands "K 1," "@," or by turning the balance off and on again.

The return parameters for the keys (keycodes) for Cubis MSU and MSA balances are:

- ▶ Softkey 5 = 1
- ▶ Softkey 4 = 2
- ▶ Softkey 3 = 3
- ▶ Softkey 2 = 4
- ▶ Softkey 1 = 5
- ▶ TASK = 6
- ▶ USER = 7
- ▶ TARE = 8
- ▶ PRINT = 9

Syntax:

Command: K 1<CR><LF>

1: Balance keys and softkeys released; keycodes not sent in the response.  
– Press the keys and softkeys to execute the corresponding function (normal mode).

K 2<CR><LF>	2: Balance keys and softkeys locked; keycodes not sent in the response. – Pressing the keys and softkeys does not execute the corresponding function and does not send a response about which key or softkey was pressed. Softkeys are not displayed.
K 3<CR><LF>	3: Balance keys and softkeys locked; keycodes sent in the response. – Pressing the keys and softkeys does not execute the corresponding function, but does send a response about which key or softkey was pressed. The softkeys are overwritten in the balance display field with "EXTERNAL 1" to "EXTERNAL 5."
K 4<CR><LF>	4: Balance keys and softkeys released; keycodes sent in response. – Pressing the keys and softkeys executes the corresponding function and sends a response about which key or softkey was pressed. Command executed
Response: K A<CR><LF> or K I<CR><LF> K L<CR><LF>	I: Command cannot currently be executed L: Syntax error – command cannot be executed

## Example 1:

Command: K 3

Response: K A  
K C 8  
K C 6Command executed  
User pressed key 8 (TARE)  
User pressed key 6 (TASK)

## Example 2:

Command: K 4

Response: K A  
K A 9  
K A 2Command executed  
User pressed key 9 (PRINT); printout executed  
User pressed key 2 (softkey 2); function executed**TA – Query and allocate tare memory**

The TA command is used to read out the content of the tare memory. If a weight value is appended to the TA command as  $w_1$  and the "Second tare" application is active, tare memory T1 is overwritten with this value. Tare memory T on the balance can only be read out (if the "Second tare" application is not active), not overwritten.

The optional parameter  $u_2$  indicates the unit of weight. If this parameter is not added, the weight unit set on the balance is used.

## Syntax:

Command: TA  $w_1$   $u_1$ <CR><LF> $w_1$ : Weight value (optional) $u_1$ : Weight unit (optional)Response: TA A  $w_1$   $u_1$ <CR><LF> $w_1$ : Weight value $u_1$ : Unit of weight

or

TA I&lt;CR&gt;&lt;LF&gt;

I: Command cannot currently be executed

## Example 1:

Command: TA

Response: TA A 129.336 g

The tare memory records a weight value of 129.336 g

## Example 2:

Command: TA 130.56 g

Response: TA A 130.560 g

"Second tare" application is active; tare memory T1 will record value 130.56 g  
Weight value 130.56 g is recorded in the tare memory

## TAC – Delete tare memory

If the "Second tare" application is active, tare memory T1 is deleted, otherwise tare memory T is deleted.

Syntax:

Command: TAC<CR><LF>

Response: TAC A<CR><LF>

or

TAC I<CR><LF>

A: Tare memory deleted

I: Command cannot currently be executed

Example:

Command: TAC

Response: TAC A

If the "Second tare" application is active, tare memory T1 is deleted, otherwise tare memory T is deleted

## SR – Send weight value if there is a weight change

The SR command is used to monitor weight values and check that they are in a certain range. If the current weight value (at or without stability) changes by the preset deviation ( $w_1$ ) at least, a weight value is sent in the response (at or without stability). The new weight value is taken as the target value for the next measurements. If no weight value is specified (SR without  $w_1$ ), 12.5% of the current weight value is taken as the monitoring limit instead. The weight unit is not currently used; instead the specified weight value is taken to be in the current weight unit used by the balance.

Syntax:

Command: SR  $w_1$ <CR><LF>

Response: S S  $w_2$   $u_1$ <CR><LF>

S +<CR><LF>

S -<CR><LF>

or

S I<CR><LF>

$w_1$ : Deviation

$w_2$ : Current weight value is taken as target value. The monitoring limits [weight value – deviation, weight value + deviation] are calculated

$u_1$ : Unit of weight

+: Balance overload

-: Balance underload

I: Command cannot currently be executed

Example 1:

Command: SR 100.00

Response: S S 199.528 g

S D 362.359 g

S S 362.358 g

The target weight value is 199.528 g. As long as the weight values remain within the range of  $\pm 100.00$  g from the target value [99.528 g, 299.528 g], no further weight value will be sent in the response

The current weight value without stability is 362.359 g. The range limit was therefore exceeded and the weight value was sent in the response

Once stability is reached again, the corresponding weight value is sent in the response and taken as the target value for new measurements. The range limits are now set at 362.358 g  $\pm 100.00$  g [262.358 g, 462.358 g]

Example 2:

Command: SR

Response: S S 199.528 g

S D 232.359 g

S S 234.247 g

The target weight value is 199.528 g. As long as the weight values remain within the range of  $\pm 12.5\%$  from the target value [174.587 g, 224.469 g], no further weight value will be sent in the response

The current weight value without stability is 232.359 g. The range limit was therefore exceeded and the weight value was sent in the response

Once stability is reached again, the corresponding weight value is sent in the response and taken as the target value for new measurements. The range limits are now set at 232.247 g  $\pm 12.5\%$  [204.966 g, 263.528 g]

## Level 2

### SU - Send weight value at stability with current weight unit (with motorized draft shield)

If the balance has a motorized draft shield and is set to automatic draft shield, the draft shield is shut first and then the weight value is sent at stability. The draft shield may open after this command is executed, depending on the motorized draft shield setting (see command M07).

Syntax:

Command: SU<CR><LF>

Response: S S w<sub>1</sub> u<sub>1</sub><CR><LF>

S: Weight value at stability

w<sub>1</sub>: Weight value

u<sub>1</sub>: Unit of weight

+: Balance overload

-: Balance underload

S +<CR><LF>

S -<CR><LF>

or

S I<CR><LF>

I: Command cannot currently be executed

Example:

Command: S

Response: S S 28 pcs Current value at stability is 28 pcs

### WS - Query door position of motorized draft shield; open or close door

Command WS queries and sets the position of the left, right, and upper doors of the draft shield.

This command only functions if the balance has a motorized draft shield and is set to automatic draft shield.

Syntax:

Command: WS n<sub>1</sub><CR><LF>

n<sub>1</sub>: 0 to 11: Door combination (optional parameter)

0 = Close all doors

1 = Open right door

2 = Open left door

3 = Open top door

4 = Open left and right doors

5 = Open left and upper doors

6 = Open right and upper doors

7 = Open all doors

10 = Left draft shield button

11 = Right draft shield button

Response: WS n<sub>2</sub><CR><LF>

n<sub>2</sub>: 0 to 9: Door position

0 = All doors closed

1 = Right door open

2 = Left door open

3 = Upper door open

4 = Left and right doors open

5 = Left and upper doors open

6 = Right and upper doors open

7 = All doors open

8 = An error has occurred

9 = Doors partially open

or

WS A<CR><LF>

A: Command executed



Example:	WS I<CR><LF>	I: Command cannot currently be executed
Command:	WS 2<CR><LF>	Open right draft shield door
Response:	WS A<CR><LF>	Right draft shield door opened
Command:	WS<CR><LF>	Query draft shield door position
Response:	WS 2<CR><LF>	Right draft shield door open

## PWR – Turn balance on/off (standby)

This command puts the balance in standby mode or turns it back on.

Syntax:		
Command:	PWR 1<CR><LF>	1: Turn balance on
	PWR 0<CR><LF>	0: Turn balance off (standby)
Response:	PWR A<CR><LF>	A: Command executed
	or	
	PWR I<CR><LF>	I: Command cannot currently be executed
Example:		
Command:	PWR 1	
Response:	PWR A	Balance is on again

## I10 – Query/set the balance ID

Command I10 queries and sets the balance ID.

This ID is retained even after the balance is switched off or after command @ has been executed.

Syntax:		
Command:	I10 "Text <sub>1</sub> "<CR><LF>	Text <sub>1</sub> : Text for ID (optional)
Response:	I10 A<CR><LF>	A: Command executed; ID saved
	or	
	I10 A "Text <sub>1</sub> "<CR><LF>	A: Command executed; ID queried
	I10 I<CR><LF>	I: Command cannot currently be executed
	I10 L<CR><LF>	L: Syntax error; command cannot be executed

Example 1:

Command:	I10 "My new ID"<CR><LF>	
Response:	I10 A<CR><LF>	"My new ID" saved to balance

Example 2:

Command:	I10<CR><LF>	Query ID
Response:	I10 A "My new ID"<CR><LF>	ID displayed

## I11 - Query balance type

This command is used to query the balance type.

Syntax:

Command: I11<CR><LF>

Response: I11 A "MSA623S"<CR><LF> A: Command executed; balance type queried

I11 I<CR><LF> I: Command cannot currently be executed

I11 L<CR><LF> L: Syntax error; command cannot be executed

## I14 - Query balance components

This command is used to query detailed information about balance components (weighing platform, display, optional data output, and draft shield).

Syntax:

Command: I14 n<sub>1</sub><CR><LF>

n<sub>1</sub>: Desired information

0 = Components

1 = Description of components

2 = Software ID number

3 = Software version

4 = Serial number

5 = IP number

Response: I14 A n<sub>1</sub> Index Info<CR><LF>

A: Command executed; information queried

I14 I<CR><LF>

I: Command cannot currently be executed

I14 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: I14 0<CR><LF>

Response: I14 B 1 "Display"<CR><LF>

I14 B 3 "Optional"<CR><LF>

I14 A 4 "Draft shield"<CR><LF>

Example 2:

Command: I14 1<CR><LF>

Response: I14 B 1 1 "MSA"<CR><LF>

I14 A 1 2 "623S"<CR><LF>

Example 3:

Command: I14 2<CR><LF>

Response: I14 B 2 1 "1C26 4482"<CR><LF>

I14 A 2 2 "00-39-74"<CR><LF>

Example 4:

Command: I14 3<CR><LF>

Response: I14 B 3 1 "01-60-06"<CR><LF>

I14 B 3 2 "00-39-74"<CR><LF>

I14 A 3 3 "04-10-03 DO\_DB25"<CR><LF>

Example 5:

Command: I14 4<CR><LF>

Response: I14 B 4 1 "327925844"<CR><LF>  
I14 A 4 2 "12345678"<CR><LF>

Example 5:

Command: I14 5<CR><LF>

Response: I14 A 5 1 "175.16.253.177"<CR><LF>

## M01 - Query/set application filter

This command is used to query or set the application filter.

Syntax:

Command: M01 n<sub>1</sub><CR><LF>

n<sub>1</sub>: Application filter (optional)

0 = Final readout

1 = Filling mode

2 = Without filtering

3 = Low filtering

Response: M01 A<CR><LF>

A: Command executed; value saved

or

M01 A n<sub>1</sub><CR><LF>

A: Command executed; application filter queried

M01 I<CR><LF>

I: Command cannot currently be executed

M01 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: M01 2<CR><LF>

Set application filter to "Low filtering"

Response: M01 A<CR><LF>

Application filter parameter set

Example 2:

Command: M01<CR><LF>

Query application filter

Response: M01 A 2<CR><LF>

Value of application filter parameter displayed

## M02 - Query/set filter adjustment

This command is used to query and adjust the filter for standard weighing on the balance. This setting is retained even after the balance is switched off or after command @ has been executed.

Syntax:

Command: M02 n<sub>1</sub><CR><LF>

n<sub>1</sub>: Filter adjustment (optional)

0 = Very stable

1 or 2 = Stable

3 = Unstable

4 = Very unstable

Response: M02 A<CR><LF>

A: Command executed; value saved

or

M02 A n<sub>1</sub><CR><LF>

A: Command executed; filter adjustment queried

M02 I&lt;CR&gt;&lt;LF&gt;

I: Command cannot currently be executed

M02 L&lt;CR&gt;&lt;LF&gt;

L: Syntax error; command cannot be executed

Example 1:

Command: M02 3&lt;CR&gt;&lt;LF&gt;

Set filter adjustment to "Unstable"

Response: M02 A&lt;CR&gt;&lt;LF&gt;

Application filter parameter set

Example 2:

Command: M02&lt;CR&gt;&lt;LF&gt;

Query filter adjustment

Response: M02 A 3&lt;CR&gt;&lt;LF&gt;

Value of filter adjustment parameter displayed

### M03 - Query/set automatic zeroing

This command is used to query and adjust automatic zeroing for standard weighing on the balance. This setting is retained even after the balance is switched off or after command @ has been executed.

Syntax:

Command: M03 n<sub>1</sub><CR><LF>n<sub>1</sub>: Automatic zeroing (optional)

0 = Deactivate

1 = Activate

Response: M03 A&lt;CR&gt;&lt;LF&gt;

A: Command executed; value saved

or

M03 A n<sub>1</sub><CR><LF>

A: Command executed; value queried

M03 I&lt;CR&gt;&lt;LF&gt;

I: Command cannot currently be executed

M03 L&lt;CR&gt;&lt;LF&gt;

L: Syntax error; command cannot be executed

Example 1:

Command: M03 1&lt;CR&gt;&lt;LF&gt;

Activate automatic zeroing

Response: M03 A&lt;CR&gt;&lt;LF&gt;

Automatic zeroing parameter set

Example 2:

Command: M03&lt;CR&gt;&lt;LF&gt;

Query automatic zeroing

Response: M03 A 1&lt;CR&gt;&lt;LF&gt;

Parameter value displayed

## M04 - Query/set I/O inputs

This command is used to allocate I/O inputs with specific functions. When parameters  $n_1$  and  $n_2$  are omitted from the command, the assignment is queried. To use the I/O inputs, first configure the appropriate settings in Menu > Configure devices > Configure interfaces > Control inputs/outputs > Peripheral port or Available ports (for example, choose "Control inputs" to allocate all five inputs for a port). This setting is retained even after the balance is switched off or after command @ has been executed.

Syntax:

Command:	M04 $n_1$ $n_2$ <CR><LF>	$n_1$ : Input 0, 1 = Not integrated in Cubis 2 = Standard data input 1 3 = Standard data input 2 4 = Standard data input 3 5 = Standard data input 4 6 = Standard data input 5 7 = Optional data input 1 8 = Optional data input 2 9 = Optional data input 3 10 = Optional data input 4 11 = Optional data input 5 $n_2$ : Function 0 = Off 1,3 = Right door of draft shield 2 = Left door of draft shield 4 = Zero 5 = Tare 6 = Print 27 = Ionizer 31 = Calibrate/adjust
Response:	M04<CR><LF> or M04 A $n_1$ $n_2$ <CR><LF> M04 I<CR><LF> M04 L<CR><LF>	A: Command executed; setting saved  A: Command executed; setting queried I: Command cannot currently be executed L: Syntax error; command cannot be executed
Example 1:		
Command:	M04<CR><LF>	Query settings
Response:	M04 B 0 0<CR><LF> M04 B 1 0<CR><LF> M04 B 2 2<CR><LF> M04 B 3 6<CR><LF> M04 B 4 0<CR><LF> M04 B 5 0<CR><LF> ... M04 A 11 0<CR><LF>	Settings displayed
Example 2:		
Command:	M04 2 5<CR><LF>	Set first standard data input to "Tare"
Response:	M04 A<CR><LF>	Parameter value saved on the balance

## M07 - Query/activate/deactivate automatic draft shield

If the balance has a motorized draft shield, this command can be used to query the automatic draft shield settings. The motorized draft shield can be activated or deactivated.

Command @ is used to disable this command.

Syntax:

Command: M07  $n_1$ <CR><LF>  $n_1$ : Automatic draft shield function (optional parameter)  
 0 = Deactivated  
 1 = Activated (close -> function -> open)  
 2 = Activated (close -> execute function)

Response: M07  $n_1$ <CR><LF>

or

M07 A<CR><LF> A: Command executed  
 M07 I<CR><LF> I: Command cannot currently be executed

Example 1:

Command: M07<CR><LF>  
 Response: M07 1<CR><LF> Automatic draft shield switched on with function  
 "Close -> function -> open"

Example 2:

Command: M07 0<CR><LF> Deactivate automatic draft shield  
 Response: M07 A<CR><LF> Command executed

## M12 - Acoustic signal (beep)

This command triggers an acoustic signal.

Syntax:

Command: M12  $n_1$ <CR><LF>  $n_1$ : Tone pitch  
 0 = Medium  
 1 = High  
 2 = Low

Response: M12 A<CR><LF> A: Command executed  
 or  
 M12 I<CR><LF> I: Command cannot currently be executed  
 M12 L<CR><LF> L: Syntax error; command cannot be executed

Example 1:

Command: M12 0<CR><LF> Trigger acoustic signal  
 Response: M12 A<CR><LF> Acoustic signal triggered

## M13 – Activate/deactivate touchscreen softkeys

Command M13 locks or releases the touchscreen softkeys.

Syntax:

Command:	M13 1<CR><LF>	1: Softkeys released
	M13 0<CR><LF>	0: Softkeys locked
Response:	M13 A<CR><LF>	A: Command executed
	or	
	M13 I<CR><LF>	I: Command cannot currently be executed

Example:

Command:	M13 0	
Response:	M13 A	Touchscreen softkeys locked

## M21 – Query/set unit of weight

This command sets or queries the current unit of weight.

Syntax:

Command:	M21<CR><LF>	Query unit of weight
Command:	M21 n <sub>1</sub> n <sub>2</sub> <CR><LF>	n <sub>1</sub> : Display unit 0,1,2 = Display (does not matter which of the three numbers is chosen)
		n <sub>2</sub> : Unit of weight
		0 = Gram                   g
		1 = Kilogram             kg
		2 = Not assigned
		3 = Milligram           mg
		4 = Microgram          µg
		5 = Carat               ct
		6 = Not assigned
		7 = Pound               lb
		8 = Ounce               oz
		9 = Troy ounce         ozt
		10 = Grain              GN
		11 = Pennyweight     dwt
		12 = Momme           mom
		13 = Mesghal          msg
		14 = Hong Kong tael   tlh
		15 = Singapore tael   tls
		16 = Taiwanese tael   tlt
		17 = Not assigned
		18 = Tola               tola
		19 = Baht               baht
		25 = No unit           --
		26 = Piece             pcs
		27 = Percent          %
		28 = User-defined unit 1   free1
		29 = User-defined unit 2   free2

Response:	M21 A<CR><LF>	A: Command executed
	or	
	M21 I<CR><LF>	I: Command cannot currently be executed
	M21 L<CR><LF>	L: Syntax error; command cannot be executed
Example 1:		
Command:	M21 1 3<CR><LF>	Set current unit to mg
Response:	M21 A<CR><LF>	Command executed
Example 2:		
Command:	M21<CR><LF>	Query current unit
Response:	M21 B 0 3<CR><LF>	Current unit is milligram
	M21 B 1 3<CR><LF>	
	M21 B 2 3<CR><LF>	

## M24 - Query/activate/deactivate "Print" key; print stable or unstable weight values

This command queries whether the "Print" key is deactivated and changes this setting if desired.

This command can also be used to specify whether weight values should be printed immediately or only once the balance has completed its work.

If the "Print" key is deactivated, this setting is only valid for the SICS interface.

Command @ is used to disable this command.

Syntax:		
Command:	M24 n <sub>1</sub> <CR><LF>	n <sub>1</sub> : "Print" key function 0 = Print stable weight value 1 = Print weight value immediately, even if unstable 2 = Deactivate "Print" key
Response:	M24 n <sub>1</sub> <CR><LF>	
	or	
	M24 A<CR><LF>	A: Command executed
	M24 I<CR><LF>	I: Command cannot currently be executed
Example 1:		
Command:	M24 <CR><LF>	
Response:	M24 2<CR><LF>	"Print" key deactivated
Example 2:		
Command:	M24 0<CR><LF>	Activate "Print" key and print only stable weight values
Response:	M24 A<CR><LF>	Command executed





## C1 – Execute calibration/adjustment (as set in menu)

Command C1 is used to trigger calibration and adjustment via the interface. The calibration/adjustment key must be set to "Fixed function" in Menu > Configure calibration/adjustment > Define calibration/adjustment functions (so that calibration and adjustment take place in one step without interruption). Set the "Fixed calibration/adjustment function" to the desired function in the same menu (for example, "Internal calibration/adjustment" or "External calibration/adjustment with standard weight"). The balance must be unloaded before command C1 is executed. Command @ can be used to cancel this command while it is being executed.

### Syntax:

Command:	C1<CR><LF>	
Response:	C1 B<CR><LF>	B: Command started
	or	
	C1 " 0.00 g"<CR><LF>	Prompt to unload the balance
	C1 " 500.00 g"<CR><LF>	Prompt to load the balance
	C1 A<CR><LF>	A: Command executed
	C1 I<CR><LF>	I: Command cannot currently be executed
	C1 L<CR><LF>	L: Syntax error; command cannot be executed

Example 1 ("Internal calibration/adjustment" has been set in the menu):

Command:	C1<CR><LF>	
Responses:	C1 B<CR><LF>	Calibration/adjustment started
	C1 A<CR><LF>	Calibration/adjustment complete

Example 2 ("External calibration/adjustment with standard weight" has been set in the menu):

Command:	C1<CR><LF>	
Responses:	C1 B<CR><LF>	Calibration/adjustment started
	C1 " 0.00 g"<CR><LF>	Prompt to unload the balance
	C1 " 500.00 g"<CR><LF>	Prompt to load the balance with 500 g
	C1 A<CR><LF>	Calibration/adjustment complete

## Remote Control

### P112 – Write text in selected line in display

A specific text is written in the selected line in the display. The number of lines is currently unlimited (recommendation: use max. 20 lines). If the text has more than 50 characters, it will be cut off at the end of the line. The "" parameter is used to write an empty line in the display and thus hide any existing text (from the active application).

### Syntax:

Command:	P112 n "Text"<CR><LF>	n: Line number Text: The text to appear in the display
Response:	P112 A<CR><LF>	A: Command executed
	or	
	P112 I<CR><LF>	I: Command cannot currently be executed

Example 1:

Command:	P112 3 "Tare the balance."	
Response:	P112 A	Text written in the 3rd line of the display



## P121 – Turn on bar graph in checkweigher

If the "Checkweighing" application is active, this command can change the SetP, max., and min. checkweighing limits and turn on the bar graph. The checkweighing limits are used in the unit set on the balance.

Syntax:	
Command:	P121 SetP Max Min<CR><LF>
	SetP: Target value Max: Maximum deviation Min: Minimum deviation
Response:	P121 A<CR><LF> or P121 I<CR><LF>
	A: Command executed I: Command cannot currently be executed
Example:	
Command:	P121 123.44 g 7.37 g 6.43 g
Response:	P121 A Limits for target value = 123.44, maximum = 123.44 + 7.37 = 130.81, and minimum = 123.44 – 6.43 = 117.01 set, and bar graph displayed again

## RM20 – Activate/deactivate user input

This command opens an edit box in the balance display so that the user can input data.

Syntax:	
Command:	RM20 n "Text <sub>1</sub> " "Text <sub>2</sub> " "Text <sub>3</sub> " <CR><LF>
	n = 1: Floating-point numbers (real) n = 2: Floating-point numbers (real) n = 8: Alphanumeric input field n = 13: Delete edit box and return to current display Text <sub>1</sub> : Input field title (max. 20 characters) Text <sub>2</sub> : Default for input field (for floating-point numbers, the number of decimal places is taken from the default) Text <sub>3</sub> : Weight unit or comment
Response:	RM20 A "P <sub>1</sub> " <CR><LF> or RM20 C<CR><LF> RM20 I<CR><LF> RM20 L<CR><LF>
	P <sub>1</sub> : User input A: Command executed C: "C" key pressed I: Command cannot currently be executed L: Syntax error; command cannot be executed
Example 1:	
Command:	RM20 1 "Reference weight" "15.000" "g"
Response:	RM20 A "22.250" Open edit box with numeric input field User input executed and input value sent back in the response
Example 2:	
Command:	RM20 8 "User name" "Name" "max. 50 characters"
Response:	RM20 A "Tom Smith" Open edit box with alphanumeric input field User input executed and user name sent back in the response
Example 3:	
Command:	RM20 13
Response:	RM20 A Delete edit box again Command executed; current task appears on the balance display

## RM30 – Assign new function to softkeys

The RM30 command assigns new functions to the softkeys (maximum 15 keys), which are displayed with the R39 command.

Syntax:

Command:	RM30 "Text <sub>1</sub> " " Text <sub>2</sub> " " Text <sub>3</sub> " ... "Text <sub>15</sub> "<CR><LF>	
		Text <sub>1</sub> : Text for softkey 1 (max. 8 characters)
		Text <sub>2</sub> : Text for softkey 2 (max. 8 characters)
		Text <sub>3</sub> : Text for softkey 3 (max. 8 characters)
		...
		Text <sub>15</sub> : Text for softkey 15 (max. 8 characters)
Response:	RM30 B<CR><LF>	A: Command executed
	or	
	RM30 I<CR><LF>	I: Command cannot currently be executed
	RM30 L<CR><LF>	L: Syntax error; command cannot be executed

Example:

Command:	RM30 "Result" "Min" "Max" ... "wRef"	
Response:	RM30 B	Command executed, but not displayed yet

## RM32 – Assign new order to softkeys

Command RM32 can be used to display the softkey assignment, previously defined with command RM32, in another order. Command @ is used to disable this command.

Syntax:

Command:	RM32 n <sub>1</sub> n <sub>2</sub> n <sub>3</sub> ... n <sub>15</sub> <CR><LF>	n <sub>1</sub> : First softkey to be displayed
		n <sub>2</sub> : Second softkey to be displayed
		n <sub>3</sub> : Third softkey to be displayed
		.....
		n <sub>15</sub> : 15th softkey to be displayed
Response:	RM32 A<CR><LF>	A: Command executed
	or	
	RM32 I<CR><LF>	I: Command cannot currently be executed
	RM32 L<CR><LF>	L: Syntax error; command cannot be executed

Example:

Command:	RM30 "Result" "Min" "Max" "wRef"<CR><LF>	
Command:	RM32 3 1 2 4<CR><LF>	
Response:	RM32 A	Softkeys displayed in the following order: "Max," "Result," "Min," and "wRef"

## RM34 – Create a dynamic parameter

Command RM34 is used to display a dynamic parameter in the working environment. The parameter is calculated as follows:  $\text{value} = n_3 * (\text{current weight value} + n_2)$ . Command @ is used to disable this command.

Syntax:

Command: RM34  $n_1$   $n_2$   $n_3$   $n_4$   $n_5$  "Text<sub>1</sub>" "Text<sub>2</sub>"<CR><LF>

$n_1$  = Lines in the working environment in which the parameter will be displayed (1 to 15)

If  $n_1 = 0$ : Delete all dynamic parameters from the display

If  $n_1 = -1$ : Delete dynamic parameters from the first line

...

If  $n_1 = -15$ : Delete dynamic parameters from the 15th line

$n_2$  = Constants to be added to or subtracted from the parameter

$n_3$  = Factor to be multiplied by the parameter

$n_4$  = Number of decimal places (0 to 9)

$n_5$  = Rounding factor (1, 2, 5, 10, 20, 50, 100)

Text<sub>1</sub>: Name of dynamic parameter (up to 20 characters)

Text<sub>2</sub>: Unit of weight for dynamic parameter (up to 6 characters)

Response: RM34 A<CR><LF>

A: Command executed

or

RM34 I<CR><LF>

I: Command cannot currently be executed

RM34 L<CR><LF>

L: Syntax error; command cannot be executed

Example:

Command: RM34 1 2.1 3 2 5 "text1" "g"<CR><LF> Show parameter in first line

Response: RM34 A<CR><LF> Command executed

## RM35 – Immediately change softkey designations

Command RM35 is used to change softkey designations immediately. Only softkeys which have been defined using command RM30 (and displayed using command RM39) are changed. The character “ cannot be used in the text. Command @ is used to disable this command.

Syntax:

Command: RM35  $n_1$  "Text<sub>1</sub>"  $n_2$  "Text<sub>2</sub>" ...  $n_{15}$  "Text<sub>15</sub>"<CR><LF>

$n_1$  = Position of first softkey to be changed

Text<sub>1</sub>: Text for first softkey to be changed (max. 8 characters)

$n_2$  = Position of second softkey to be changed

Text<sub>2</sub>: Text for second softkey to be changed (max. 8 characters)

.....

$n_{15}$  = Position of 15th softkey to be changed

Text<sub>15</sub>: Text for 15th softkey to be changed (max. 8 characters)

Response: RM35 A<CR><LF>

A: Command executed

or

RM35 I<CR><LF>

I: Command cannot currently be executed

RM35 L<CR><LF>

L: Syntax error; command cannot be executed

Example:

Command: RM30 "Result" "Min" "Max" "wRef"<CR><LF>

Command: RM35 2 "nRef" 4 "Set" 5 "Next"<CR><LF>

Response: RM35 A Softkeys displayed in the following order: "Result," "nRef," "Min," "Set," and "Next"

## RM36 – Assign/query function for multiple softkey lines

The command RM36 assigns or queries functions for up to 30 softkey allocations (up to a maximum of 15 keys). This is activated on the balance with the R38 command.

Syntax:

Command: RM36  $n_1$  "Text<sub>1</sub>" "Text<sub>2</sub>" "Text<sub>3</sub>" ... "Text<sub>15</sub>"<CR><LF>

$n_1 = 0$ : Display all allocations

$n_1 = 1$  to 30: Number of softkey allocation

Text<sub>1</sub>: Text for softkey 1 (max. 8 characters)

Text<sub>2</sub>: Text for softkey 2 (max. 8 characters)

Text<sub>3</sub>: Text for softkey 3 (max. 8 characters)

...

Text<sub>15</sub>: Text for softkey 15 (max. 8 characters)

Response: RM36 A<CR><LF>

A: Command executed

or

RM36 I<CR><LF>

I: Command cannot currently be executed

RM36 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: RM36 3 "Result" "Min" "Max" "wRef"

Assign 3rd softkey allocation

Response: RM36 A

Third softkey allocation assigned, but not displayed yet

Example 2:

Command: RM36 3

Query third softkey allocation

Response: RM36 3 "Result" "Min" "Max" "wRef"

Third softkey allocation sent back in the response

## RM37 – Prepare preset softkey designations for display

Command RM37 is used to copy a softkey allocation, which has previously been defined using command RM36, to command RM30. This allocation can be activated with "RM39 1" on the balance. It is even easier (without command RM37) to display allocations defined with command RM36 using command RM38. Command @ is used to disable this command.

Syntax:

Command: RM37  $n_1$ <CR><LF>

$n_1$  = Softkey allocation previously defined using command RM36

Response: RM37 A<CR><LF>

A: Command executed

or

RM37 I<CR><LF>

I: Command cannot currently be executed

RM37 L<CR><LF>

L: Syntax error; command cannot be executed

## Example 1:

Commands: RM36 3 "Result" "Min" "Max" Assign third softkey allocation  
 "wRef"  
 RM36 3 Assign third softkey allocation to RM30, but do not display yet  
 RM39 1 Display this softkey allocation

Responses: RM36 A  
 RM37 A  
 RM39 A

## Example 2: Example 1 can also be carried out as follows:

Commands: RM36 3 "Result" "Min" "Max" Assign third softkey allocation  
 "wRef"  
 RM38 3 Display third softkey allocation

Responses: RM36 A  
 RM38 A

**RM38 – Activate RM36-assigned softkey lines**

The RM38 command activates the assigned softkey allocations (assigned using the RM36 command) or assigns a new sequence to them.

Syntax:  
 Command: RM38 n1 P1<CR><LF> n1: Allocation number (1 to 30)  
 P1: Optional parameter that can be used to change the sequence (alphabetical order: ABC...MNO) for the softkeys

Example 1:  
 Where there are five softkey allocations, key 1 should be swapped with key 3 – enter "CBADE" as the parameter

Example 2:  
 With eight softkeys, swap key 2 with key 5 and key 4 with key 8 – the parameter is then "AECHBFGD"

Response: RM38 A<CR><LF> A: Command executed  
 or  
 RM38 I<CR><LF> I: Command cannot currently be executed  
 RM38 L<CR><LF> L: Syntax error; command cannot be executed

## Example 1:

Command: RM38 3 Important! RM36 command must be executed beforehand!  
 Response: RM38 A Softkeys now have new functions that were previously defined with the M36 command for the third allocation

## Example 2:

Command: RM38 2 EACBD Important! RM36 command must be executed beforehand!  
 Response: RM38 A Softkeys now have new functions that were previously defined with the M36 command for the second allocation, but with a changed sequence: "Text<sub>5</sub>" "Text<sub>1</sub>" "Text<sub>3</sub>" "Text<sub>2</sub>" "Text<sub>4</sub>"



### RM39 – Activate/deactivate RM30–assigned softkey functions

The RM39 command activates and deactivates or deletes the functions assigned to softkeys (with the RM30 command).

**Syntax:**

Command:	RM39 P <sub>1</sub> <CR><LF>	P <sub>1</sub> = 0: Previous allocation of keys is deleted. P <sub>1</sub> = 1: Softkeys are overwritten with new functions P <sub>1</sub> = 2: Softkeys are overwritten with functions for the current task
Response:	RM39 A<CR><LF> or RM39 I<CR><LF>	A: Command executed I: Command cannot currently be executed

**Example:**

Command:	RM39 1	Important! RM30 command must be executed beforehand!
Response:	RM39 A	The softkeys overwritten from left to right with the assigned functions (based on the RM30 command). If more than five functions were entered, the fifth softkey is labeled "More." Pressing the fifth softkey displays the other softkeys (5 up to a maximum of 15) in a pop-up window.

### RM44 – Query/set input with barcode scanner

Barcode scanners (or keypad input) cannot be locked in Cubis.

**Syntax:**

Command:	RM44 n <sub>1</sub> <CR><LF>	Change input n <sub>1</sub> = Index for locking barcode scanner input 0: Barcode scanner input only activated with command RM20 1: Barcode scanner always active
	RM44<CR><LF>	Query
Response:	RM44 A n <sub>1</sub> <CR><LF> or RM44 I<CR><LF> RM44 L<CR><LF>	A: Command executed I: Command cannot currently be executed L: Syntax error; command cannot be executed

**Example 1:**

Command: RM44 1<CR><LF>

Response: RM44 A<CR><LF>

**Example 1:**

Command: RM44<CR><LF>

Response: RM44 A 1<CR><LF>

**RM48 – Change order of standard keys**

Command RM48 is used to display several standard softkeys in any order. Command @ is used to disable this command.

Syntax:

Command:	RM48 n <sub>1</sub> n <sub>2</sub> n <sub>3</sub> n <sub>4</sub> n <sub>5</sub> n <sub>6</sub> n <sub>7</sub> n <sub>8</sub> <CR><LF>	n <sub>1</sub> : Allocation number n <sub>2</sub> : Position of "Internal adjustment" key n <sub>3</sub> : Position of "External adjustment" key n <sub>4</sub> : Position of "Internal calibration" key n <sub>5</sub> : Position of "External calibration" key n <sub>6</sub> : Position of key 1/10d n <sub>7</sub> : Not implemented n <sub>8</sub> : Not implemented
Response:	RM48 A<CR><LF> or RM48 I<CR><LF> RM48 L<CR><LF>	A: Command executed  I: Command cannot currently be executed L: Syntax error; command cannot be executed

Example 1:

Command:	RM48 1 1 5 2 4 3 0 0<CR><LF>	Display standard softkeys
----------	------------------------------	---------------------------

Response:	RM48 A<CR><LF>
-----------	----------------

Example 2:

Command:	RM48 0 0 0 0 0 0 0 0<CR><LF>	Hide standard softkeys
----------	------------------------------	------------------------

Response:	RM48 A<CR><LF>
-----------	----------------

## RM49 – Activate/deactivate info text

Command RM49 displays informational text. When the text is empty ("") the info text is not shown. Use "\x09" to add a tab to the text, and "\x0D" to add a line break. Command @ is used to disable this command.

Syntax:

Command: RM49  $n_1$   $n_2$  "Text"<CR><LF>

Display info text

$n_1$  = Displayed keys in window

1: Window without keys

2: Window with OK key

3: Window with C key

4: Window with OK and C key

$n_2$  = Additional information about the text

1: Info

2: Alert

3: Stop

4: Question

5: X-mark

6: Date/time

7: Sound

8: Alphabetic

9: Numeric

10: Hourglass

Text = Info text to be displayed in the window

or

RM49 0<CR><LF>

Hide info text

Response: RM49 B<CR><LF>

B: Info text displayed

or

RM49 A 1<CR><LF>

A 1: "OK" key pressed. Info text no longer displayed

RM49 A 2<CR><LF>

A: "C" key pressed. Info text no longer displayed

RM49 I<CR><LF>

I: Command cannot currently be executed

RM49 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: RM49 2 3 "Weight is too heavy!"<CR><LF>

Display info text and "OK" key

Response: RM49 B<CR><LF>

Example 2:

Command: RM49 0<CR><LF>

Hide info text

Response: RM49 A<CR><LF>

Example 3:

Command: RM49 2 3 ""<CR><LF>

Hide info text

Response: RM49 A<CR><LF>

## RM51 – Activate/deactivate selection window

Command RM51 is used to define a list to be displayed together with the desired keys and additional information on the display as a selection. Use "\x09" to add a tab to the text. Command @ is used to disable this command.

Syntax:

Command: RM51  $n_1$   $n_2$   $n_3$   $n_4$   $n_5$  "Text<sub>1</sub>" "Text<sub>2</sub>" ... "Text<sub>17</sub>"<CR><LF>

$n_1$  = Number of selected entry ( $n_1 = 1$  to 15)

If  $n_1 = 0$  or  $n_1 = 20$ : No entry is selected

If  $n_1 = 21$  to 25: Entry  $n_1 = 20$  is selected

$n_2$  = Additional information

0: Hidden

1: Displayed

$n_3$  and  $n_4$  were not implemented in Cubis, because scrolling with the slider is possible

$n_5$  = How additional information (Text<sub>2</sub>) should be displayed

0: Additional information in text form

1: Additional information in softkey form

Text<sub>1</sub>: Info

Text<sub>2</sub>: Additional information in text or function key form

Text<sub>3</sub>: First entry on the list

Text<sub>4</sub>: Second entry on the list

...

Text<sub>17</sub>: 15th entry on the list

Command: RM51<CR><LF>

Close window

Command: RM51 0<CR><LF>

Close window

Response: RM51 B<CR><LF>

B: Command executed. System is waiting for user selection

or

RM51 F  $n$ <CR><LF>

Line  $n$  from the list selected by user

RM51 A C<CR><LF>

"Cancel" key pressed

RM51 A K<CR><LF>

Function key pressed

RM51 A P<CR><LF>

P: "Prev" key pressed

RM51 A N<CR><LF>

N: "Next" key pressed

RM51 I<CR><LF>

I: Command cannot currently be executed

RM51 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: RM51 3 1 1 1 1 "Fruit selection" "All" "Apple" "Orange" "Kiwi" "Banana" "Strawberry" "Peach"<CR><LF>

Response 1: RM51 B<CR><LF>

Command executed. System is waiting for user selection

Response 2: RM51 F 3<CR><LF>

User has selected the third entry ("Kiwi")

Example 2:

Command: RM51<CR><LF>

Response: RM51 B<CR><LF>

Command executed. Window closed



**RM54 – Activate/deactivate window with info**

Command RM54 is used to define and activate/deactivate windows using an info list. Use "\x09" to add a tab to the text, and "\x0D" to add a line break. Command @ is used to disable this command.

## Syntax:

Command: RM54 n<sub>1</sub> n<sub>2</sub> "Text<sub>1</sub>" "Text<sub>2</sub>"<CR><LF>    n<sub>1</sub> = Keys

- 1: No keys in window
- 2: "OK" key only
- 3: "C" key only
- 4: User-defined key only
- 5: "OK" and "C" keys
- 6: "OK" and user-defined keys
- 7: "C" and user-defined keys
- 8: "OK," "C," and user-defined keys

n<sub>2</sub> = Number of additional text (defined using command RM49)  
Text<sub>1</sub>: Label for user-defined key  
Text<sub>2</sub>: Text for info window

Command:	RM54<CR><LF>	Close window
Response:	RM54 B<CR><LF>	Command executed. Window displayed
	RM54 A 1<CR><LF>	"OK" key pressed. Window closed
	RM54 A 2<CR><LF>	"C" key pressed. Window closed
	RM54 A 3<CR><LF>	User-defined key pressed. Window closed
	RM54 I<CR><LF>	Command cannot currently be executed
	RM54 L<CR><LF>	Syntax error; command cannot be executed

## Example:

Command:	RM54 8 1 "Next" "Gross \x0923.4 g\x0DSerial number is 1234567."<CR><LF>	Add two lines of text: "Gross 23.4 g" and "Serial number is 1234567." Display all three keys and info text "Info"
Response:	RM54 A<CR><LF>	Command executed

## Additional Sartorius Commands

### SA – Send weight value at stability and store in Alibi memory

The SA command calls up a weight value at stability and stores it in the Alibi memory. As an option, a label can be assigned when the weight value is stored in the Alibi memory.

If the balance has a motorized draft shield and is set to automatic draft shield, the draft shield is shut first and then the weight value is sent at stability. The draft shield may open after this command is executed, depending on the motorized draft shield setting (see command M07).

Syntax:

Command:	SA "Text"<CR><LF>	Text:	Label (optional)
Response:	SA A "w <sub>1</sub> " "w <sub>2</sub> " "w <sub>3</sub> " "w <sub>4</sub> " "w <sub>5</sub> " "n <sub>1</sub> " "n <sub>2</sub> " "Text"<CR><LF>	w <sub>1</sub> :	Net weight value
		w <sub>2</sub> :	Balance tare memory
		w <sub>3</sub> :	Appl. tare memory 1
		w <sub>4</sub> :	Appl. tare memory 2
		w <sub>5</sub> :	Gross weight value
		n <sub>1</sub> :	Balance serial number
		n <sub>2</sub> :	Consecutive number of data record in Alibi memory
		Text:	Label (optional), if entered
		A:	Command executed
	or		
	S I<CR><LF>	I:	Command cannot currently be executed

Example 1:

Command: SA "Art. 23"

Response: SA A "N2 228.86[6] g" "T 0.00[0] g" "T1 0.00[0] g"  
"T2 99.50[5] g" "G# 328.37[1] g" "Ser No. 23201202" "Mem No. 503" "Mem ID Art. 23"

Example 2:

Command: SA

Response: SA A "N1 173.51[1] g" "T 0.00[0] g" "PT1 125.00[0] g"  
"T2 0.00[0] g" "G# 298.51[1] g" "Ser No. 23201202" "Mem No. 504" "Mem ID"

### CM – Execute application command

This command executes application commands.

Syntax:

Command:	CMD "m <sub>1</sub> .c <sub>1</sub> " P <sub>1</sub> <CR><LF>	m <sub>1</sub> :	Application module
		.	Separator for application module and command
		c <sub>1</sub> :	Application command
		P <sub>1</sub> :	Optional parameter
Response:	CMD "m <sub>1</sub> .c <sub>1</sub> " P <sub>1</sub> A<CR><LF>	A:	Command executed
	or		
	CMD I<CR><LF>	I:	Command cannot currently be executed

Example 1:

Command: CMD WEIGH.DO\_TARE1 1

Response: CMD WEIGH.DO\_TARE1 1 A Balance tared

Example 2:		
Command:	CMD MESSAGE.SHOW_ERROR	"Weight is too low!"
Response:	CMD MESSAGE.SHOW_ERROR	"Weight is too low!" A Error message displayed
Example 3:		
Command:	CMD MESSAGE.HIDE_ERROR	
Response:	CMD MESSAGE.HIDE_ERROR A	Error message deleted from display
Example 4:		
Command:	CMD RECIPE.START	
Response:	CMD RECIPE.START A	Formulation application started

## PAR – Query parameter

This command queries the values of a current valid parameter.

Syntax:		
Command:	PAR "m <sub>1</sub> .P <sub>1</sub> "<CR><LF>	m <sub>1</sub> : Application module . : Separator for application module and parameter P <sub>1</sub> : Parameter from application module
Response:	PAR A h <sub>1</sub> v <sub>1</sub> <CR><LF>	h <sub>1</sub> : Header of queried parameter v <sub>1</sub> : Value of queried parameter A: Command executed
	or	
	PAR I<CR><LF>	I: Command cannot currently be executed
Example 1:		
Command:	PAR USER.TITLE	Query name of active user
Response:	PAR A User Tom Smith	Header and name of active user sent back in the response
Example 2:		
Command:	PAR CHECK.MIN	Query minimum limit for checkweighing application
Response:	PAR A 12.230 g	Minimum limit for active application is sent back in the response
Example 3:		
Command:	PAR DENSITY.RHO_SAM	Query density of sample being weighed
Response:	PAR A 1.4 g/cm <sup>3</sup>	Density of current sample is sent back in the response
Example 4:		
Command:	PAR TASK.TITLE	Query name of active task
Response:	PAR A Task Determine density	Header and name of active task sent back in the response
Example 5:		
Command:	PAR COUNT.WREF	Query average weight for the piece count application
Response:	PAR A 9.95010 g	Current average piece weight sent back in the response



## MN36 – Assign a function to several menus

Command MN36 assigns functions for up to 30 menu allocations (up to a maximum of 30 entries). This is activated on the balance with the MN38 command.

Syntax:

Command: MN36 n<sub>1</sub> "Text<sub>1</sub>" "Text<sub>2</sub>" "Text<sub>3</sub>" ... "Text<sub>15</sub>" <CR><LF>

n<sub>1</sub>: 1 to 30: Number of menu allocation  
 Text<sub>1</sub>: Text for menu entry 1 (max. 30 characters)  
 Text<sub>2</sub>: Text for menu entry 2 (max. 30 characters)  
 Text<sub>3</sub>: Text for menu entry 3 (max. 30 characters)  
 ...  
 Text<sub>30</sub>: Text for menu entry 30 (max. 30 characters)

To stop text from being displayed, enter \~  
 before the text (e.g., "\~Next")

To use the character "\" enter "\\"

Command: MN36 A<CR><LF>                   A:       Command executed  
 or  
 MN36 I<CR><LF>                        I:       Command cannot currently be executed  
 MN36 L<CR><LF>                        L:       Syntax error; command cannot be executed

Example:

Command: MN36 3 "Select" "Next" "Previous"

Response: MN36 A                       Assign third menu allocation  
   Third menu allocation assigned, but not displayed yet



## TX36 – Assign text to several text pages

Command TX36 is used to assign text to up to 30 text pages (with 15 lines of text each). This text is activated on the balance using command TX38 and can be changed using command TX37.

Syntax:

Command: TX36  $n_1$  "Text<sub>1</sub>" "Text<sub>2</sub>" "Text<sub>3</sub>" ... "Text<sub>15</sub>"<CR><LF>

$n_1$  = 1 to 30: Number of text page

Text1: Text for line 1 (max. 30 characters)

Text2: Text for line 2 (max. 30 characters)

Text3: Text for line 3 (max. 30 characters)

...

Text15: Text for line 15 (max. 30 characters)

Response: TX36 A<CR><LF>

A: Command executed

or

TX36 I<CR><LF>

I: Command cannot currently be executed

TX36 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: TX36 3 "Tare balance" "Place sample on balance" "Press [Next]"

Assign text to third text page

Response: TX36 A

Third text page assigned, but not displayed yet

## TX37 – Overwrite a line on a text page

Command TX37 is used to overwrite a selected line on a selected text page. Command TX36 must be executed for this text page beforehand. Command TX38 displays this page.

Syntax:

Response: TX37  $n_1$   $n_2$  "Text"<CR><LF>

$n_1$  = 1 to 30: Number of text page

$n_2$  = 1 to 15: Number of text line

Text: Text for line  $n_2$  (max. 30 characters)

Response: TX37 A<CR><LF>

A: Command executed

or

TX37 I<CR><LF>

I: Command cannot currently be executed

TX37 L<CR><LF>

L: Syntax error; command cannot be executed

Example 1:

Command: TX37 3 2 "Place third sample on the balance"

Important! TX36 command must be executed beforehand!

Response: TX37 A

Second line overwritten on the third page

## TX38 – Activate/deactivate TX36–assigned text pages

The TX38 command activates/deactivates the assigned text pages (assigned using the TX36 command) or assigns a new sequence to them. This is a faster alternative to command P112, which writes individual text lines to the display.

Syntax:

Command:	TX38 n <sub>1</sub> <CR><LF>	n <sub>1</sub> :	1 to 30: Number of text page
Response:	TX38 A<CR><LF>	A:	Command executed
or	TX38 I<CR><LF>	I:	Command cannot currently be executed
	TX38 L<CR><LF>	L:	Syntax error; command cannot be executed

Example 1:

Command:	TX38 3	Important! TX36 command must be executed beforehand!
Response:	TX38 A	Working environment now has new text that was previously defined with the TX36 command for the third allocation

Example 2:

Command:	TX38 0	Deactivate text page
Response:	TX38 A	



Sartorius Weighing Technology GmbH  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289  
[www.sartorius-mechatronics.com](http://www.sartorius-mechatronics.com)

Copyright by Sartorius,  
Goettingen, Germany.  
No part of this publication may be  
reprinted or translated in any form or  
by any means without the prior written  
permission of Sartorius.  
All rights reserved. The status of  
the information, specifications and  
illustrations in this manual is indicated  
by the date given below. Sartorius  
reserves the right to make changes to  
the technology, features, specifications,  
and design of the equipment without  
notice.

Date:  
September 2011  
Sartorius Weighing Technology GmbH  
Goettingen, Germany

Specifications subject to change  
without notice. RS · KT  
Publication no.: WMS6008-e11093