



Syntax description

Sartorius

MINI - SICS Interface



Contents

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

## 1. Intended Use

MINI-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 access to the balance (query parameters, executing commands, saving weight values to alibi memory). Automatic printing is available. Use a program such as the SartoTerminal from Sartorius to communicate with the balance.

The following commands are available:

- S – Send weight value at stability
- SI – Send weight value without stability
- SIR – Send automatic weight values at and without stability
- T – Tare the balance at stability
- TI – Tare the balance without stability

Additional Sartorius commands:

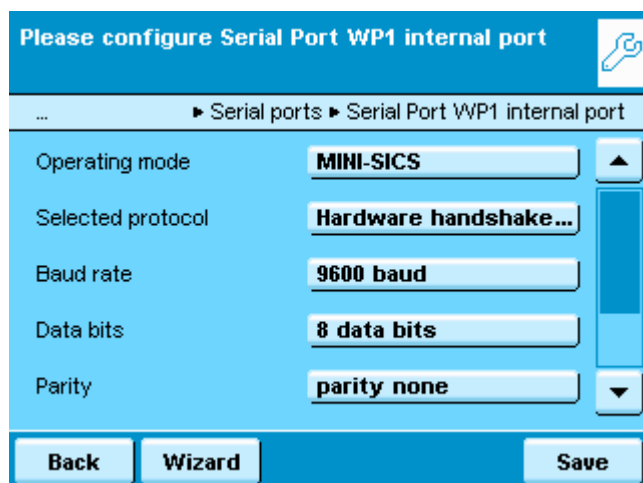
- SA – Send weight value at stability and store in alibi memory (with optional label)
- CMD – Execute application command
- PAR – Query parameter

## 2. Settings for Cubis MSA and MSU Balances

MINI-SICS commands can be used via the following interfaces:

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

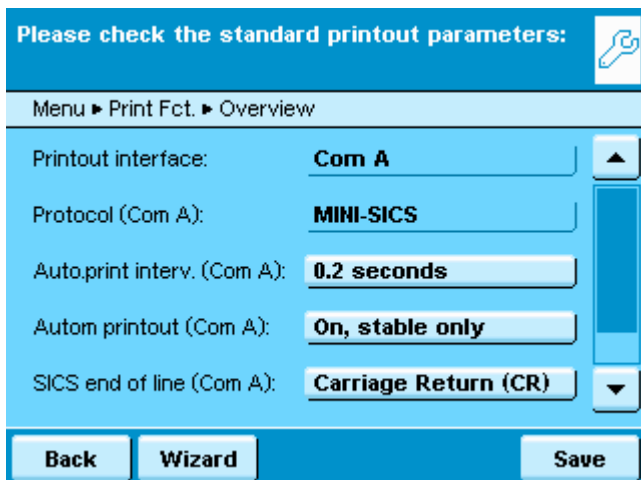
In "Menu ->Device Parameters ->Configure ports -> Configure serial ports," the interface must be set to "MINI-SICS" mode:



## Additional settings:

- ▶ Selected protocol:
  - 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
- ▶ Record data:
  - Turn off
  - Turn on

For using autom. printing in standard weighing, standard weighing shall be activated and parameter "autom. printout" in Menu -> Device parameters -> Configure Data output shall be on:



- ▶ Auto. print interval:
  - 1 time per second
  - 2 times per second (0.5 sec.)
  - 5 times per second (0.2 sec)
  - 10 times per second (0.1 sec.)
- ▶ Autom. printout:
  - Off
  - On, stable and unstable weight values
  - On, stable weight values only

For using autom. printing in task, configure parameter "autom. printout" in task configuration:

Please check the printout parameters of this task:

Task > New > Print Fct. > Overview

Printout interface: **Com A**

Protocol (Com A): **MINI-SICS**

Auto.print interv. (Com A): **0.2 seconds**

Autom printout (Com A): **On, stable only**

SICS end of line (Com A): **Carriage Return (CR)**

**Back** **Wizard** **Done** **Next**

### 3. Syntax of MINI-SICS Commands

A MINI-SICS command consists of an identifier (ID) and optional parameters.

ID [Parameter1] [Parameter2] [Parameter3] ... [Parametern]<CR><LF>

The identifiers (IDs) consist of ASCII characters and are written in capital or small letters. The parameters must be separated with a space. Each command must be concluded with carriage return and line feed (#0D#0A or <CR><LF>).

Example: Get weight value                      S<CR><LF>

### 4. Syntax of MINI-SICS Responses

The balance sends a response to each MINI-SICS command except T and TI. The responses may contain one or more weight values and/or texts. After starting MINI-SICS, balance sends response "AT".

Characters 1-2: command ID  
 Character 3: space  
 Characters 4-12: weight value  
 Character 13: space  
 Characters 14...19 unit (without spaces)  
 <CR><LF>

Response for additional Sartorius commands:

ID Status [Parameter1] [Parameter2] [Parameter3] ... [Parametern]<CR><LF>

„Status“ reproduces how the command was executed:

- A – Command was executed, no further response will be sent
- 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 texts.

## 5. Description of MINI-SICS Commands

**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.

Syntax:

Command: S<CR><LF>

Response: S w1 u1<CR><LF>

S: Weight value at stability

w1: Weight value

u1: Units of weight

S+<CR><LF>

+: Balance overload

S-<CR><LF>

-: Balance underload

or

SI <CR><LF>

I: Command cannot currently be executed

Example:

Command: S<CR><LF>

Response: S 99.528 g <CR><LF>

Current weight value at stability is 99.528 g

**SI – Send weight value without stability**

Syntax:

Command: SI<CR><LF>

Response: S w1 u1<CR><LF>

S: Weight value at stability

SD w1 u1<CR><LF>

D: Weight value without stability

w1: Weight value

u1: Units of weight

SI+<CR><LF>

+: Balance overload

SI-<CR><LF>

-: Balance underload

or

SI<CR><LF>

I: Command cannot currently be executed

Example:

Command: SI<CR><LF>

Response: S 99.528 g <CR><LF>

Current weight value at stability is 99.528 g

SD 362.359 g <CR><LF>

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 the parameter "Interval For Automatic Printout") – once, twice, 5 times, or 10 times a 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&lt;CR&gt;&lt;LF&gt;

Response: S w1 u1&lt;CR&gt;&lt;LF&gt;

SD w1 u1&lt;CR&gt;&lt;LF&gt;

SI+&lt;CR&gt;&lt;LF&gt;

SI-&lt;CR&gt;&lt;LF&gt;

or

SI &lt;CR&gt;&lt;LF&gt;

S: Weight value at stability

D: Weight value without stability

w1: Weight value

u1: Units of weight

+: Balance overload

-: Balance underload

I: Command cannot currently be executed

## Example:

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

Response: SD -94.821 g&lt;CR&gt;&lt;LF&gt;

SD 228.896 g&lt;CR&gt;&lt;LF&gt;

S 228.890 g&lt;CR&gt;&lt;LF&gt;

**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.

This command sends no response.

## Syntax:

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

## Example:

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

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

**T1 – Tare the balance 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: T1&lt;CR&gt;&lt;LF&gt;

## Example:

Command: T1

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





Example 4:

Command: CMD RECIPE.START

Response: CMD RECIPE.START A

Recipe application started

## PAR – Query parameter

This command queries the values of a current valid parameter.

Syntax:

Command: PAR "m1 P1"<CR><LF>

m1: Application module

.: Separator for application module and parameter

P1: Parameter from application module

Response: PAR A h1 v1<CR><LF>

h1: Header of queried parameter

v1: Value of queried parameter

A: Command executed

or

PAR I<CR><LF>

I: Command cannot currently be executed

Example 1:

Command: PAR

Name of active user to be queried

Response: PAR A User Tom Smith

Header and name of active user are returned

Example 2:

Command: PAR CHECK.MIN

Minimum limit for checkweighing application to be queried

Response: PAR A 12.230 g

Minimum limit for active application is returned

Example 3:

Command: PAR DENSITY.RHO\_SAM

Density of current sample to be queried

Response: PAR A 1.4 g/cm3

Density of current sample is returned

Example 4:

Command: PAR TASK.TITLE

Name of active task to be queried

Response: PAR A Task Dichtebest.

Header and name of active task are returned

Example 5:

Command: PAR COUNT.WREF

The average weight for the piece count application to be queried

Response: PAR A 9.95010 g

The current average weight is returned

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,  
Göttingen, 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.

Information status:  
März 2010, Sartorius,  
Goettingen, Germany

Specifications subject to change with-  
out notice. RS · KT  
Publication No.: WMS6010-e11031