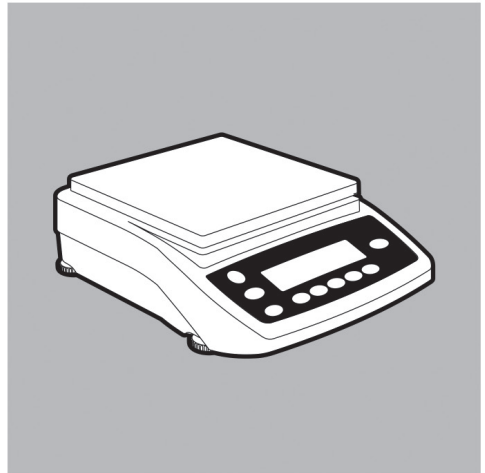
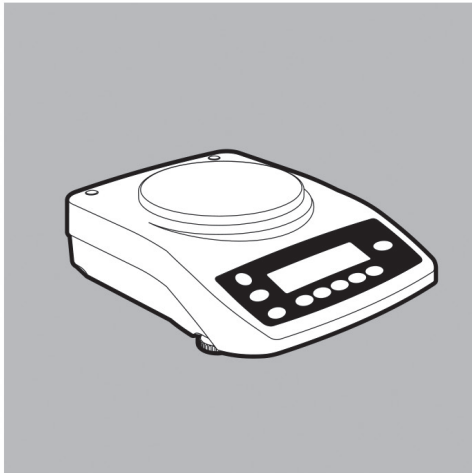
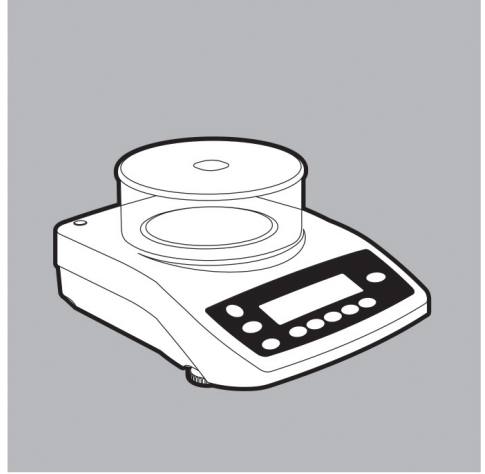
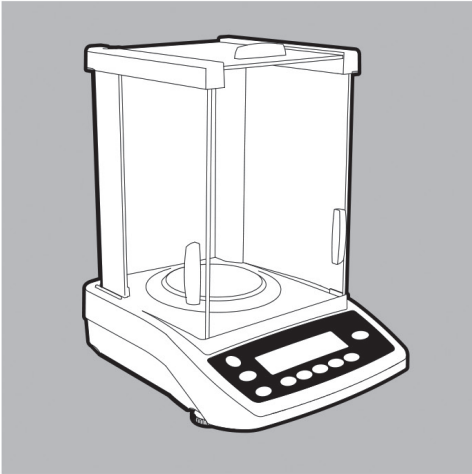


한글 사용설명서

Sartorius Entris

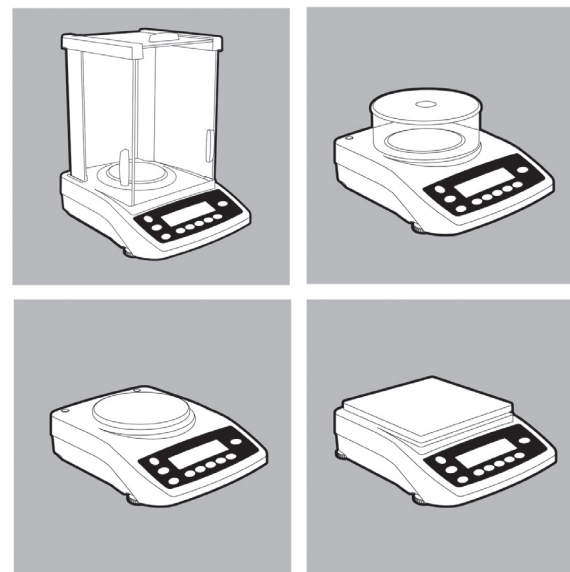
Laboratory Balances



한글 사용설명서

Sartorius Entris

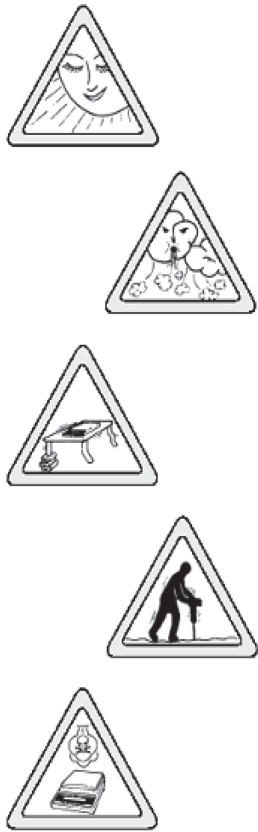
Laboratory Balances



목 차

| | |
|--------------------------------------|----|
| 설치 환경 및 보관 운송 ----- | 4 |
| 설치 순서 ----- | 5 |
| 전원 ----- | 6 |
| 주변기기 접속 / 워밍업 / 수평 조절 ----- | 7 |
| 하부계량 ----- | 8 |
| 디스플레이 설명 ----- | 9 |
| Calibration and Adjustment(보정) ----- | 10 |
| Calibration 하는 방법 ----- | 11 |
| 메뉴 설정 ----- | 12 |
| 파라미터 셋팅 ----- | 14 |
| 장치정보 ----- | 19 |
| Counting ----- | 20 |
| Weighing in percent ----- | 22 |
| Animal weighing/ Averaging ----- | 24 |
| Mass Unit Conversion ----- | 26 |
| Density Determination ----- | 28 |
| Data interface ----- | 29 |
| 문제 발생 시 ----- | 30 |
| 유지 관리 ----- | 31 |
| 제품 사양 ----- | 31 |
| 액세서리 ----- | 35 |
| 제품보증서 ----- | 37 |

설치 환경



싸토리우스의 Entris Series저울은 연구실, 공장내의 어떠한 조건 하에서도 신뢰할 수 있는 계량 결과를 제공하도록 설계되어 있습니다. 아래의 유의점을 지켜 설치 장소를 선택하면, 더욱 빠르고 정확한 계량이 이루어집니다.

- 저울은 안전한 면에 설치한다.
- 태양과 히터의 고열을 피한다.
- 열린 창문이나 문으로부터의 통풍을 피한다.
- 극심한 진동을 피한다.
- 충격, 부식성 가스로부터 저울을 보호한다.

저울을 폭발위험이 있는 장소나 지역에서는 사용하지 마십시오. 저울을 장시간 습도가 높은 곳을 놓아두지 마십시오. 저울을 저온에서 고온의 장소로 이동시킬 경우 공기 중의 수분이 차가운 저울의 표면에 내려 앉아 결로가 발생할 수 있으므로 이러한 경우에는 저울의 전원을 연결하기 전 두시간 정도 워밍업을 하여 환경에 적응시키십시오. 그 후 전원을 연결시키면 저울내외의 온도차에 의한 습도의 영향을 피할 수 있습니다.

보관과 운송에 대하여

허용 보관온도 -40°C 에서 $+70^{\circ}\text{C}$

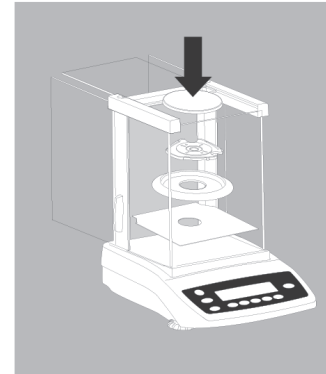
포장은 최대 1m 의 높이에서 떨어져도 내용물인 전자저울은 파손되지 않도록 설계되어 있습니다.

포장을 뜯은 후, 운송에 의한 파손이 있는지 확인해 주십시오.

파손된 경우 싸토리우스 코리아 (주)로 연락주시면 바로 조치하여 드리겠습니다.

포장재는 나중의 운송을 위해 보관해 두시면 편리하며, 재포장을 할 경우 파손방지를 위해 모든 접속케이블을 제거하여 주십시오. 저울은 극도의 고·저온, 다습, 충격, 진동 등을 피해 주십시오.

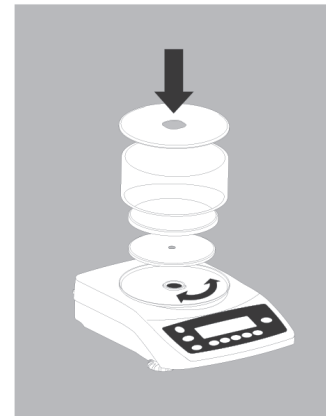
설치 순서



1. 0.1 mg / 사각유리챔버 설치

아래에 순서대로 내용물을 설치하여 주십시오

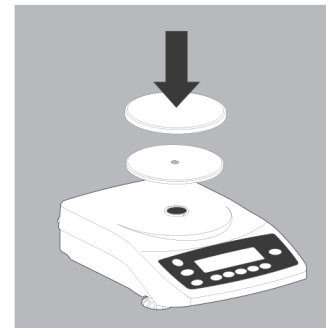
- Shield plate (바닥 판)
- Shield ring (실드 링)
- Pan support (팬 서포트)
- Weighing pan (계량 접시)



2. 1mg / 원형 유리챔버의 설치

아래에 순서대로 내용물을 설치하여 주십시오.

- 바닥뚜껑을 정위치에 놓는다.
- Pan support (팬 서포트)
- Weighing pan (계량 팬)
- Glass draft shield (원형 유리챔버)
- The upper lid (상판 뚜껑을 닫는다)



3. 삼각형 팬 (원형 계량접시의 경우)

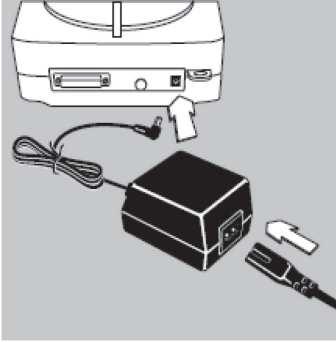
아래에 순서대로 구성품들을 설치한다.

- Pan support (팬서 포트)
- Weighing pan (계량 접시)

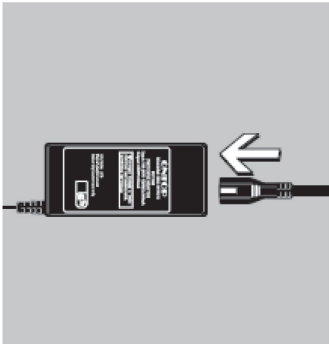
4. 사각형의 접시

- Weighing pan (계량 접시)

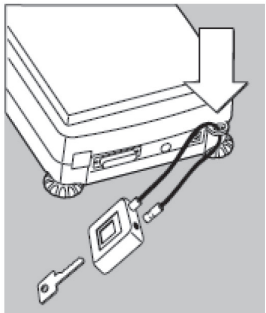
전원



- 싸토리우스에서 공급 되어진 정품 아답터만을 사용하셔야 안전하게 제품을 이용하실 수 있습니다.
- 뒷면 Angle plug에 AC 아답터를 연결한다.
- 파워코드를 연결한다.
- 전원을 넣는다.

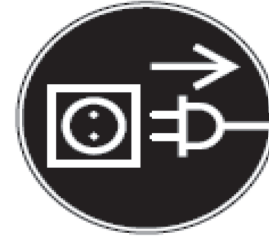


- 저울은 AC 아답터의 전원으로 작동합니다.
- AC 아답터의 기재되어 있는 전압과 사용지역의 전원이 일치하는지 확인해 주십시오.
- 최대용량이 10kg인 저울의 아답터는 110V or 220V 겸용입니다.
(단, 반드시 싸토리우스에서 공급된 정품 아답터를 사용하셔야 합니다.)



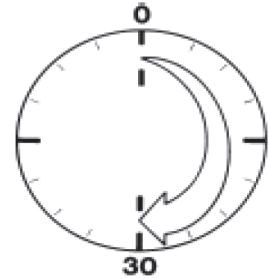
- 필요 시 도난방지용 잠금장치를 체인이나 자물쇠로 잠그면 어디에서도 안전하게 이용하실 수 있습니다.

주변기기 접속 / 워밍업 / 수평 조절



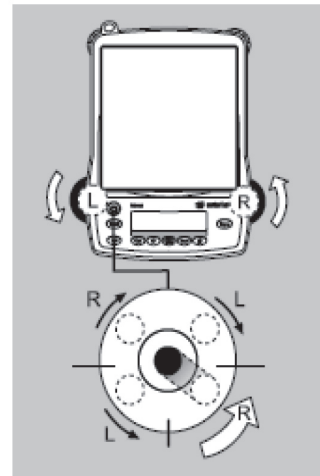
▶ 주변기기 접속

- 본체연결포트(프린터, 컴퓨터, 인터페이스 연결포트에 연결 또는 분리하기 전에 저울의 전원장치를 꼭 분리시키십시오.



▶ 워밍업

- 작동 전 30분간 워밍업을 해야 기기의 작동온도에 도달하며, 전자저울의 정확한 성능을 보증할 수 있습니다.

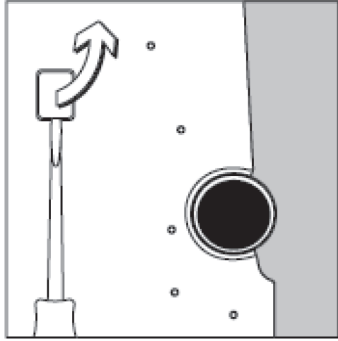


▶ 수평 조절

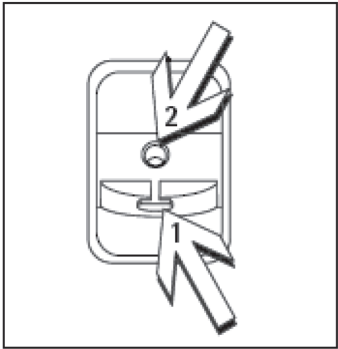
사용지점에 따라 수평조절기 다리를 적당히 돌려서 수평기의 물방울이 원의 중앙에 오도록 조정해야, 저울의 수평을 맞출 수 있습니다.

- 물방울이 상단에 위치 시 : 시계방향
- 물방울이 하단에 위치 시 : 반시계방향

하부계량



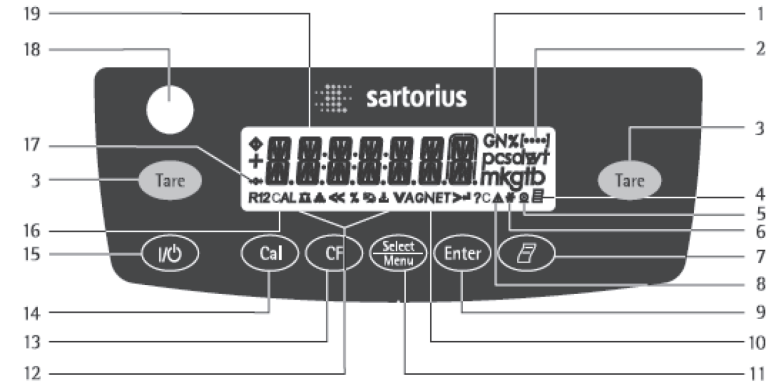
- 하부계량
전자저울의 바닥면에 하부계량을 할 수 있는 고리가 장착되어 있습니다. 계량법상으로는 비공식적인 방법입니다.
- 아래쪽 커버를 열어 주십시오.
(이때, 저울을 뒤집지 마십시오.)



- 내장된 고리(1)의 사용: 고리에 샘플을 부착할 수 있습니다.
(예: 고리선 사용)

디스플레이 설명

디스플레이 전체보기



- | | |
|---|--|
| 1. Weighing unit (단위) | 14. Start calibration/Routine (보정키) |
| 2. Menu level indicator (메뉴 표시) | 15. On/Off (전원키) |
| 3. Taring (0점, 용기소거키) | 16. Symbol : Calibration function |
| 4. Symbol "GLP printout" | 17. Symbols for zero range (Verified models only) (0점 표시) |
| 5. Symbol "Printing mode" (출력 표시) | 18. Level indicator (물방울) |
| 6. Symbol "Application program active" | 19. Weight value displayed in selected weight unit (액정) |
| 7. Data output (프린터키) | |
| 8. Calculated-value indicator | |
| 9. Start an application program (시작키) | |
| 10. Symbol (Gross or net value) (총계 표시) | |
| 11. Select an application program / Open the operating menu (메뉴 선택/메뉴 진입) | |
| 12. Symbols for active application (\square , \triangle , $\%$, \square , \square , \square , \square) (application 심볼) | |
| 13. Delete(Clear function) (취소키) this key is generally used to cancel functions – Quit application program – Cancel calibration routine/Exit the operating menu. | |

Symbols

- << 저장 및 ESC
- < 상위메뉴로 이동
- √ 아래메뉴로 이동
- > 하위메뉴로 이동
- ↩ 선택키

Calibration and Adjustment (보정)

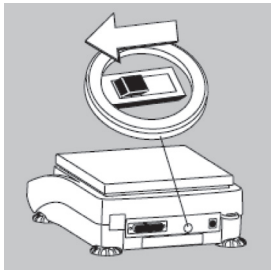
▶ 목 적

Calibration은 표시된 값과 실제 값 사이의 차이를 결정하는 것이며, Adjustment는 차이 값의 보정을 의미한다.

▶ 특 징

보정 시 필요 조건 들

- 저울 짐판 위에는 아무것도 올려 놓지 않는다.
- Tare(영점)를 잡는다.
- 0.000g이 화면에 표시되어야 한다.
- 오차 2% 내에서만 보정 가능. (예 : 200g 보정시 2% 내 오차범위 내 질량 필요)
- ※만약 위의 조건을 충족 못할 시 "Error 02"가 나타남.




▶ 외부교정

도량형법에 의해 외부 교정만으로 사용되어질 경우

- 메뉴에서 CAL- ADJ : Block 설정한다.
- 저울뒷면에 어세스 스위치를 왼쪽 그림처럼 닫는다.

▶ 내부교정 (I 가 포함된 모델 만 가능- 예, Entris224i-1S)

- Cal 키를 이용하여 내부분동으로 교정한다.

Press  Cal 키를 누른다.
 내장된 분동이 자동적으로 실행된다.
 저울이 보정된다.
 자동적으로 weighing mode로 돌아온다.

※ -I 가 포함된 모델은 Internal calibration이 가능한 모델.



Calibration 방법

▶ 내부교정

[단 계]

1. 0점을 잡는다. (Tare 키 이용)
2. Calibration키를 누른다.
(내부분동을 이용하여 자동적으로 동작한다.)
3. Calibration이 실행된다.
4. 내부분동이 제거 된후 차이 값이 보정된다.

[메커니즘]

 0.0 g
 CAL.INT.
 CAL.RUN.
 CAL.END
 0.0 g



▶ 외부교정

Parameters (changes in factory settings):
 SETUP: BAL.SCAL.: CAL.-ADJ.: CAL.EXT. (menu code 1.1.9.1)

위와 같이 설정을 변경한다.






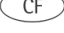
[단 계]

1. 0점을 잡는다. (Tare 키 이용)
2. Cal 버튼을 누른다.
(요구하는 무게를 올린다.)
3. 표시된 분동을 저울짐판위에 올린다.
4. Calibration이 실행된다.
5. 분동을 제거한다.

 0.0 g
 CAL.EXT.
 - 5000.0 g
 5000.0 g
 CAL.END
 + 5000.0 g
 0.0 g








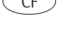




메뉴 설정

▶ 메뉴 설정을 위한 버튼 설명


| Symbol | Key | 버튼 설명 |
|--------|--|--------------------|
| V |  | 메뉴 화면이동 |
| > |  | 현재 메뉴의 아랫단계로 이동 |
| ↵ |  | 메뉴설정 |
| |  (Press and hold) | 저장 후 현재 위치에서 빠져나옴. |
| << |  | 현재 메뉴의 윗단계로 이동 |
| < |  | 메뉴단계의 표시 |
| [....] | | |

메뉴 네비게이션 (예시)

▶ 언어선택

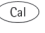

| Key (or instruction) | Display | |
|---|---|--|
|  hold | APPLIC. | 1. 메뉴화면을 연다. |
| Repeatedly  | INPUT ... | 2. 메뉴화면에 진입 후 ↓ 버튼을 이용하여 Language를 선택한다. |
| Repeatedly  | LANGUAG. | |
|  | ENGLISH ° | 3. 원하는 언어 선택한다. (English in factory) |
|  | GERMAN | 4. 독일어를 선택한다. |
|  | GERMAN ° | 5. Enter키를 누르면 오른쪽 상단에 "°" 이 표시된다 |
|  | LANGUAG. | 6. CF 버튼을 이용하여 빠져나온다. |
|  ,  | | |
| Repeatedly:  |  | 7. CF 버튼을 반복적으로 눌러서 빠져나온다. |
|  | | 8. 저장없이 빠져나올때는 전원버튼을 이용한다. |
| | 0.0 g | |


Parameter Settings (Overview)







| Level 1 [•] | Level 2 [••] | Level 3 [•••] | Menu code |
|------------------------|---|---|--------------|
| SETUP | BAL.SCAL. Balance scale parameters | AMBIENT conditions | 1. 1. 1. |
| | | APP.FILT. Application filter | 1. 1. 2. |
| | | STAB.RNG.Stability range | 1. 1. 3. |
| | | TARING Taring 1) | 1. 1. 5 |
| | | AUTZER. Auto zero | 1. 1. 6 |
| | | WT.UNIT Basic weight unit | 1. 1. 7. |
| | | DISP. Acc Display accuracy | 1. 1. 8. |
| | | CAL./ADJ. Function of the  key | 1. 1. 9. |
| | | CAL.UNIT. weight unit for calibration | 1. 1.11. |
| | | BAUrate | 1. 5. 1. |
| | INTERF. Interface | PARITY Parity | 1. 5. 2. |
| | | STOPBIT Number of stop bits | 1. 5. 3. |
| | | HANDSHK. Handshake mode | 1. 5. 4. |
| | | DATABIT Number of data bits | 1. 5. 5. |
| | | DAT.REC. Output: SBI (ASCII) or printout | 1. 5. 6. |
| | PRNT.OUT Settings for print function | PRINT (manual automatic) | 1. 6. 1. |
| | | STOPAUT. Stop automatic printing | 1. 6. 2. |
| | | AUT.CYCL. Time-dependent autom. printing | 1. 6. 3. |
| | | TAR./PRT. Tare bal./balance after ind. print | 1. 6. 4. |
| | | PRT.INIT. Printout of appl. parameters | 1. 6. 5. |
| | EXTRAS (Additional functions) | FORMAT Line format for printout | 1. 6. 6. |
| | | MENU Menu read only can edit | 1. 8. 1. |
| | | SIGNAL Acoustic signal (beep) | 1. 8. 2. |
| | | KEYS (Keypad) | 1. 8. 3. |
| | | EXT.KEY External switch function | 1. 8. 4. |
| | RESET | ONMODE Power-on mode | 1. 8. 5. |
| | | BACKLIT Display backlighting | 1. 8. 6. |
| | | MENU Factory settings | 1. 9. 1. |
| APPLIcation programs | WEIGH | | 2. 1. |
| | UNIT Toggle wt. unit | DISP.DIG. Display accuracy | 2. 2. 2. |
| | COUNT.ing | RESOLUT. Resolution | 2. 3. 1. |
| | | REF.UPDT. Autom. ref. sample updating | 2. 3. 2. |
| | PERCENT Weighing in percent | DEC.PLCS Decimal places | 2. 4. 1. |
| | ANIMALW.eighing | ACTIVITY. Animal activity | 2. 7. 1. |
| | | START | 2. 7. 2. |
| | CALC.ulation | METHOD (Operator) | 2. 8. 1. |
| | | DEC.PLCS Decimal places | 2. 8. 2. |
| | DENSITY determination | DEC.PLCS Decimal places | 2. 9. 1. |
| INPUT Input | ID NO. | ID input; max. 7 characters, e.g. as Inventory no. | 3. 1. |
| | | | |
| INFORMATION | VERSION, SER.NO., MODEL | Display software ver., serial no., model | 4. 1. .2 .3. |
| LANGUAG. (LANGUAG.) | ENGLISH (factory setting) | | 5. 1. |
| | DEUTSCH (German) | | 5. 2. |
| | FRANçais (French) | | 5. 3. |
| | ITAL.iano (Italian) | | 5. 4. |
| | ESPAÑOL (Spanish) | | 5. 5. |
| | РУССКОЕ (Russian) | | 5. 6. |
| | POLSKI (Polish) | | 5. 7. |
| | CODES Menu shows codes (not texts) | | 5. 8. |

Parameter settings: Overview

o = Factory setting; ✓ = User-defined setting



| Level 1 [•] | Level 2 [••] | Level 3 [•••] | Level 4 [••••] | Menu code |
|--------------|--|--|---|------------|
| SETUP | BAL.SCAL. Balance parameters | AMBIENT conditions (Filter adaptation) | o VERY STABLE | 1. 1. 1. 1 |
| | | | o STABLE | 1. 1. 1. 2 |
| | | | o UNSTABLE | 1. 1. 1. 3 |
| | | | o VERY UNSTABLE | 1. 1. 1. 4 |
| | | APP.FILT. Application filter | o FINAL.RD | 1. 1. 2. 1 |
| | | | FILLING | 1. 1. 2. 2 |
| | | STABILITY range | 1/4 DIG. (digit) | 1. 1. 3. 1 |
| | | | 1/2 DIG. (digit) | 1. 1. 3. 2 |
| | | | 1 - DIGIT (digit) | 1. 1. 3. 3 |
| | | | o 2 - DIGIT (digit) | 1. 1. 3. 4 |
| | | | 4 - DIGIT (digit) | 1. 1. 3. 5 |
| | | | 8 - DIGIT (digit) | 1. 1. 3. 6 |
| | | TARING Taring | o W/O STAB (W/o stability) | 1. 1. 5. 1 |
| | | | o W/ STAB (After stability) | 1. 1. 5. 2 |
| | | AUT.ZERO Auto zero | o OFF | 1. 1. 6. 1 |
| | | | o ON | 1. 1. 6. 2 |
| | WT.UNIT Basic weight through unit | | For list of units, see Chapter | 1. 1. 7. 1 |
| | | | Toggling between weight units" | 1. 1. 7.23 |
| | DISP.DIG. Display accuracy | | o ALL | 1. 1. 8. 1 |
| | | | o MINUS 1 | 1. 1. 8. 2 |
| | | | o DIVIS. 11 interval | 1. 1. 8. 6 |
| | CAL./ADJ. Function of the of the  key | | CAL.EXT. External cal./adj. | 1. 1. 9. 1 |
| | | | o CAL.INT Internal cal./adj. | 1. 1. 9. 2 |
| | | | KEY BLOCKED  blocked | 1. 1. 3. 3 |
| | CAL.UNIT Unit for calibration weight | | o GRAMS | 1. 1.11. 1 |
| | | | o KILOGRAMS | 1. 1.11. 2 |
| | | | o POUNDS | 1. 1.11. 3 |

| Level 1 [•] | Level 2 [••] | Level 3 [•••] | Level 4 [••••] | Menu code |
|-----------------|---------------------------|---|---|------------|
| SETUP | INTERF. Interface | BAUDRATE | 600 | 1. 5. 1. 3 |
| | | | o 1200 | 1. 5. 1. 4 |
| | | | 2400 | 1. 5. 1. 5 |
| | | | 4800 | 1. 5. 1. 6 |
| | | | 9600 | 1. 5. 1. 7 |
| | | | 19200 | 1. 5. 1. 8 |
| | | PARITY Parity | o ODD | 1. 5. 2. 3 |
| | | | EVEN | 1. 5. 2. 4 |
| | | | NONE | 1. 5. 2. 5 |
| | | STOPBIT No. of stop bits | o 1 STOPBIT | 1. 5. 3. 1 |
| | | | 2 BITS | 1. 5. 3. 2 |
| | | HANDSHK. Handshake mode | SOFTWARE | 1. 5. 4. 1 |
| | | | o HARDWARE. | 1. 5. 4. 2 |
| | | | NONE | 1. 5. 4. 3 |
| | | DATABIT No. of data bits | o 7 BITS | 1. 5. 5. 1 |
| | | | 8 BITS | 1. 5. 5. 2 |
| | | DAT.REC. Com- munication mode | 5BI (ASCII) | 1. 5. 6. 1 |
| | | | o PRINTER | 1. 5. 6. 2 |
| | PRNT.OUT Printing fct. | PRINT (manual automatic) | MAN.W/O stability | 1. 6. 1. 1 |
| | | | o MAN.WITH. stability | 1. 6. 1. 2 |
| | | | AUT.W/O stability | 1. 6. 1. 3 |
| | | | AUT.WITH. stability | 1. 6. 1. 4 |
| | | STOPAUT. Stop automatic printing | o OFF Not possible | 1. 6. 2. 1 |
| | | | ON Use print key  | 1. 6. 2. 2 |
| | | AUT.CYCL. Time-dependent autom. printing | o EACHVAL (1 display update) | 1. 6. 3. 1 |
| | | | AFTER 2 (2 display updates) | 1. 6. 3. 2 |
| | | TAR./PRT. Tare the balance after individual printout | o OFF | 1. 6. 4. 1 |
| | | | ON | 1. 6. 4. 2 |

| Level 1 [•] | Level 2 [••] | Level 3 [•••] | Level 4 [••••] | Menu code |
|-----------------|-------------------------------------|---|--|------------|
| SETUP | PRNT.OUT Printing fct. | PRT.INIT. Printing applica- tion parameters | OFF | 1. 6. 5. 1 |
| | | | o ALL parameters | 1. 6. 5. 2 |
| | | | MAINPAR.ameters | 1. 6. 5. 2 |
| | | | FORMAT Line for- mat for printout | |
| | | | o 16.CHAR.(digit not identified) | 1. 6. 6. 1 |
| | | | 22.CHAR.acters (w/ 1D) | 1. 6. 6. 2 |
| | EXTRAS (Additional functions) | MENU | CANCELIT.le | 1. 8. 1. 1 |
| | | | RDONLY read only | 1. 8. 1. 2 |
| | | SIGNAL Acoustic signal | OFF | 1. 8. 2. 1 |
| | | | o ON | 1. 8. 2. 2 |
| | | KEYS (keypad) | o FREE | 1. 8. 3. 1 |
| | | | LOCKED | 1. 8. 3. 2 |
| | | EXT.KEY External switch function | o PRINT Key  | 1. 8. 4. 1 |
| | | | Z I TARE Key  | 1. 8. 4. 2 |
| | | | CAL. Key  | 1. 8. 4. 3 |
| | | | SELECT Key  | 1. 8. 4. 4 |
| | | | CF Key  | 1. 8. 4. 5 |
| | | | ENTER Key  | 1. 8. 4. 6 |
| | | ON-MODE Power-on mode | o OFF/ON Off on stand-by | 1. 8. 5. 1 |
| | | | STANDBY On Stand-by | 1. 8. 5. 2 |
| | | | AUTO-ON Automatic on mode | 1. 8. 5. 3 |
| | | BACKLIT Display backlighting | OFF | 1. 8. 6. 1 |
| | | | o ON | 1. 8. 6. 2 |
| | RESET Menu reset | MENU factory settings | YES restore factory settings | 1. 9. 1. 1 |
| | | | o NO Do not restore factory settings | 1. 9. 1. 2 |

| Level 1 [●] | Level 2 [●●] | Level 3 [●●●] | Level 4 [●●●●] | Menu code |
|------------------------------|---|--------------------------------|---|------------|
| APPL. ication programs | WEIGHING toggling between weight UNITS | CONFIG. display accuracy | o ALL | 2. 1. |
| | | | MINUS 1 | 2. 2. 2. 1 |
| | | | DIV. 1 1 interval | 2. 2. 2. 2 |
| | COUNTING | RESOLUT. | o DISP.DIG. Display accuracy | 2. 2. 2. 6 |
| | | | 10 FOL 10 times > display | 2. 3. 1. 1 |
| | | | | 2. 3. 1. 2 |
| | REF.UPDT. Auto Reference- updating | o OFF | 2. 3. 2. 1 | |
| | | AUTOM.atic | 2. 3. 2. 2 | |
| | | | | |
| | PERCENT Weighing in percent | DEC.PLCS. Decimal places | NONE no decimal places | 2. 4. 1. 1 |
| | | | o 1 DEC.PL. 1 decimal place | 2. 4. 1. 2 |
| | | | 2 DEC.PL. 2 decimal places | 2. 4. 1. 3 |
| | ANIMALW. Animal weighing | ACTIVITY Animal activity | 3 DEC.PL. 3 decimal places | 2. 4. 1. 4 |
| | | | CALM (fluct.: 2% of test obj.) | 2. 7. 1. 1 |
| | | | o ACTIVE (fluct.: 5% of test obj.) | 2. 7. 1. 2 |
| | START | | V. ACTIVE (fluct.: 20% of test obj.) | 2. 7. 1. 3 |
| | | | MANUAL | 2. 7. 2. 1 |
| | | | o AUTOM.atic | 2. 7. 2. 2 |
| | DENSITY determination | DEC.PLCS decimal places | NONE no decimal places | 2. 9. 1. 1 |
| | | | o 1 DEC.PL. 1 decimal place | 2. 9. 1. 2 |
| | | | 2 DEC.PL. 2 decimal places | 2. 9. 1. 3 |
| | | 3 DEC.PL. 3 decimal places | 2. 9. 1. 4 | |


장치정보

| Level 1 [•] | Level 2 [••] | Level 3 [•••] | Example | Code |
|-----------------|------------------|--|---------------|-------|
| INFO rmation | VER.NO. | Show software version | REL.32.09 | 4. 1. |
| | SER.NR. | Show serial number, e.g.: (To toggle focus between upper and lower display sections: Press ) | 29712345 | 4. 2. |
| | MODEL | Show model designation (to change focus from upper to middle to lower display section and back: Press ) | ENTRIS 124-15 | 4. 3. |

Display of Menu Items: Selecting Languages or Codes

| | | |
|------------|------------------------------------|-------|
| LANGUAG. | ENGLISH (factory setting) | 5. 1. |
| (LANGUAG.) | DEUTSCH (German) | 5. 2. |
| | FRAN.çais (French) | 5. 3. |
| | ITAL.iano (Italian) | 5. 4. |
| | ESPAÑOL (Spanish) | 5. 5. |
| | РУССКИЙ (Russian) | 5. 6. |
| | POLSKI (Polish) | 5. 7. |
| | CODES Menu shows codes (not texts) | 5. 8. |




Counting

Display symbol: 

▶ 목 적

같은 무게를 가진 샘플의 수량이 많을 시 수량을 정확하게 파악할 수 있습니다. 처음 기준 샘플들의 무게를 이용하여 기준값을 만든 후 일정 수의 샘플들을 전자저울에 올려 놓으면 올려놓은 개수를 알 수 있는 기능이다.

▶ 기준 샘플수량 변화 방법

| [키버튼] | [내 용] |
|---|---|
|  key | 1. 키를 누른다. |
| | 2. 기준 샘플의 개수를 입력한다. (1에서 100까지) (정확한 개수 측량을 위해서는 수량을 많이 입력한다.) |
|  key | 3. 키를 이용하여 원하는 개수를 선택한다. |
|  key | 4. 선택 후 길게 키를 길게 누르고 있다. (입력 완료) |

▶ 기준 샘플수량 업데이트

자동 비교 샘플의 업그레이드는 카운팅기능의 정확도를 위해 활용된다. 활성화 또는 비활성화가 가능하며, 자동 비교 샘플 업데이트 기능은 충족되어 질 때 수행된다. (디스플레이상에 OPT 가 표시됨)

▶ 준비단계

메뉴 안 Counting application 선택한다. (아래의 그림은 factory setting이 된 모습)

APPL IC.ation programs

| | |
|-------------|-----------------------------|
| COUNT. | |
| RESOLUT.ion | |
| o DISP.ACC. | Display accuracy |
| 10-FOLD | 10-fold higher |
| REF.UPDT. | Autom. ref. sample updating |
| o OFF | Off |
| AUTOM. | Automatic |

o = Factory setting

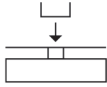

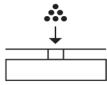




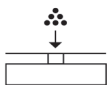


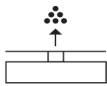

프린팅 된 모습

Printout: Counting
 nRef 10 : Ref. sample quantity
 wRef 21.14 g : Reference weight
 for 1 unit
 Qnt + 500 pcs : Calculated quantity

(기준 1개의 무게가 21.14 g)

예제

같은 샘플의 개수 측정 – 파라미터 셋팅 : application : count (menu code 2,3)

| Step | Key (or instruction) | Display/Data output |
|-------------------------------------|--|--|
| 1. 빈 용기를 저울위에 올려놓는다. |  | + 22.6 g |
| 2. 0점을 잡는다. (Tare 키를 이용) |  | 0.0 g |
| 3. 기준 샘플을 올린다. (예 : 20개 정도) |  | |
| 4. 비교 샘플의 양을 변화시키기 원하면 다음키를 누른다. |  | REF 10 pcs |
| 5. 원하는 기준 샘플 개수를 선택한다. | Repeatedly:  Press briefly  press and hold | REF 20 pcs |
| 6. 비교샘플의 개수와 비교 무게를 확인한다. |  | + 20 pcs * nRef 20 pcs wRef 1.07 g |
| 7. 측정하고자 하는 개수 만큼 샘플을 저울에 올려 놓는다. |  | + 500 pcs |
| 8. 프린터 출력을 원하면, 프린트 버튼을 누른다. |  | Qnt + 500 pcs + 1.07 g Δ* + 535.0 g * + 500 pcs * |
| 9. 총 무게와 개수를 버튼을 이용 하여 볼 수 있다. | Repeatedly:  | |
| 10. 샘플 제거 |  | - 2 pcs * |
| 11. 반복적으로 측정을 원하면 step7 부터 다시 측정한다. | | |
| 12. 기준값을 삭제한다. |  | 0.0 g |




Weighing in Percent

Display symbol: %

▶ 목 적

이 기능은 기준 중량을 비율로 변환하여 확인하고자 할 때 사용되는 기능입니다.

▶ 기준 샘플수량 변화 방법

| [키버튼] | [내 용] |
|---|---|
|  key | 1. 키를 누른다. |
| | 2. 기준 샘플의 개수를 입력한다. (1에서 100까지) (정확한 개수 측량을 위해서는 많은 수량을 입력한다.) |
|  key | 3. 키를 이용하여 원하는 개수를 선택한다. |
|  key | 4. 선택 후 키를 길게 누르고 있다. (입력 완료) |

▶ 준비단계

메뉴의 Weighing in percent application을 선택한다. (아래의 그림은 factory setting된 모습)

| APPL I.C.ation programs | 프린팅 된 모습 |
|-----------------------------|----------|
| PERCENT Weighing in percent | |
| DEC.PLCS Decimal places | |
| o NONE Decimal places | |
| 1 DEC.PL. 1 decimal place | |
| 2 DEC.PL. 2 decimal places | |
| 3 DEC.PL. 3 decimal places | |

o = Factory setting

Printout: Weighing in percent



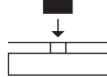

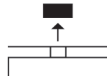
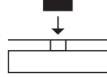




| | | |
|------|----------|--|
| pRef | 100 | : Reference percentage |
| Wxx% | 111.6 g | : Reference weight for selected reference percentage xx% |
| Prc | + 94.9 % | : Calculated reference percentage |

(111.6g 을 100%로 설정된 모습을 나타냄)


예제

퍼센트 중량의 결정 – 파라미터 셋팅 : application : Percent (menu code 2.4)

기준 퍼센트 : REF 100%

| Step | Key (or instruction) | Display/Data output |
|-----------------------------------|---|---|
| 1. 0점을 잡는다. |  | 0.0 g |
| 2. 설정된 기준 퍼센트 확인 |  | REF 100 % |
| 3. 100%에 해당하는 무게의 비교샘플을 올려 놓는다. |  | |
| 4. Enter키를 이용하여 확인한다. |  | + 100.0 % * pRef 100 % Wxx% + 111.6 g |
| 5. 기준 샘플을 제거한다. |  | |
| 6. 모르는 무게의 샘플을 올려놓는다. |  | + 94.9 % * |
| 7. 원하는 퍼센트가 나타난다. |  | Prc + 94.9 % |
| 8. 버튼을 이용하여 실제 중량 퍼센트 중량을 볼 수 있다. | Repeatedly:  | + 105.9 g * + 94.9 % * |
| 9. 기준샘플의 입력값을 제거하려면 CF 이용하여 실시한다. |  | + 105.9 g |
| 10. 필요 시 프린터 출력 |  | N + 105.9 g |

Animal weighing / Averaging




Display symbol: 

▶ 목 적

불안정한 샘플이나 불안정한 환경에서 측정시 사용합니다.

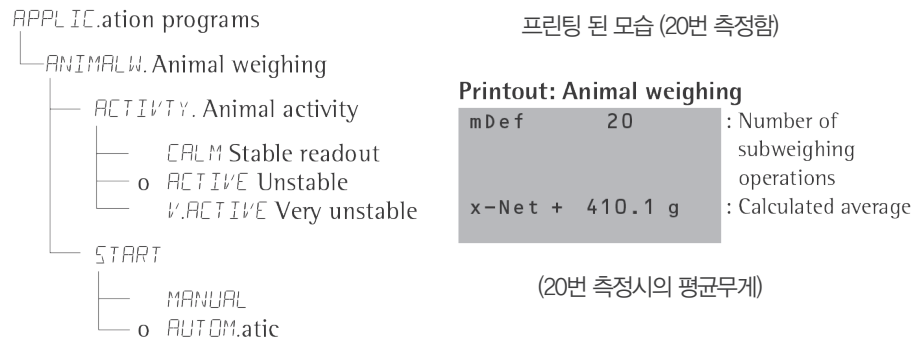
또, 여러가지의 샘플의 평균값을 이 프로그램을 통해 계산할 수 있습니다.

▶ 기준 샘플수량 변화 방법

| [키버튼] | [내 용] |
|---|-----------------------------------|
|  key | 1. 키를 누른다. (원하는 수를 선택) (1...:100) |
|  key | 2. 키를 이용하여 원하는 개수를 선택한다. |
|  key | 3. 선택 후 길게 키를 길게 누르고 있다. (입력 완료) |

▶ 준비단계

Animal weighing application을 setting한다. (아래의 그림은 factory setting된 모습)

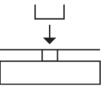




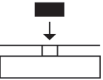
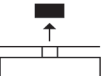


o = Factory setting

예제

자동 시작이 설정된 20번 측정의 동물계량

파라미터 셋팅 : Application : animal weighing (menu code 2.7.)

| Step | Key (or instruction) | Display/Data output |
|-------------------------------------|--|--|
| 1. 저울 위에 동물계량용 접시를 놓는다. |  | 22.6 g |
| 2. 0점을 잡는다. |  | 0.0 g |
| 3. 버튼을 눌러 횟수 확인 |  | REF 30 |
| 4. 원하는 횟수를 입력한다. (20 입력) | Repeatedly: Press briefly press and hold  | REF 20 |
| 5. Enter키를 이용 저장. |  | + 0.0 g * |
| 6. 동물계량용 접시를 올려 놓은 후 움직이는 샘플을 측정한다. |  | 888 19 20 ... 1 |
| 7. 횟수만큼 계산된 값이 나타난다. | | + 410.1 g Δ* mDef 20 x-Net + 410.1 g |
| 8. 샘플제거 후 원하면 다음 샘플을 측정한다. |  | + 0.0 g * |

Mass Unit Conversion

▶ 목 적








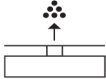

이 기능을 이용해서 디스플레이 상에 다른 무게 단위를 나타낼 수 있습니다.

▶ 특 징

- 기본적인 단위를 셋팅할 수있다. (표 참조)
- 셋팅 된 단위는 저장되어 있다.
- 전원을 넣으면 항상 기본으로 설정되어진 단위가 나타난다.

예제

pound (lb), ounces(ozt)로 변환 – 파라미터 셋팅

| Step | Press key | Display Printout |
|--|--|-------------------------------------|
| 1. Select / Menu 버튼을 누른다. |  | NONE ⁰ [•] |
| 2. pound(lb)를 찾는다. | Repeatedly:  | POUNDS |
| 3. Enter 키를 이용 입력한다. |  | POUNDS ⁰ |
| 4. Select / Menu 버튼을 눌러 ounces(ozt)를 찾는다.(TROY OZ) |  , Repeatedly:  | NONE ⁰ [••] TROY.OZ. |
| 5. Enter키를 입력한다. |  | TROY.OZ. ⁰ |
| 6. 또 다른 단위의 설정을 원하면 위와 같이 입력한다. | | [•••] |
| 7. CF키를 눌러 빠져나온다. |  | 0.00 g |
| 8. 샘플을 올려 놓는다. |  | + 100.00 g |
| 9. Toggling 을 이용 단위변환 가능. (Enter키 이용) | Repeatedly:  | + 0.22046 lb + 3.5275 ozt |

※ 다음의 단위들은 세계에서 공통적으로 사용되어지는 단위입니다.

그러나 한국에서는 모든 단위를 사용 할 수 없으며, 법정계량 단위 만을 이용해야 합니다.

(상업적 목적 사용 시 mg, g, kg, cat만 사용 가능. 연구용 목적 시 제외)

The following weight units are available in your balance: :

| Menu item | Unit | Conversion factor | Display symbol |
|----------------------------|------------|-------------------|----------------|
| 1) USERDEF. 1) | Grams | 1.000000000000 | o |
| 2) GRAMS (Factory setting) | Grams | 1.000000000000 | g |
| 3) KILOGR. | Kilograms | 0.001000000000 | kg |
| 4) MILLIGR. | Milligrams | 1000.0000000000 | mg |

Density Determination

Display symbol: $\Delta\Delta$




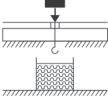

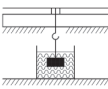



▶ 목 적

이 기능을 부력법을 이용하여 고체의 밀도를 측정할 수 있습니다. 측정을 위해서는 홀더와 매달림 와이어를 사용하여야 측정할 수가 있으며, 저울에는 포함되어 있지 않습니다.

예제

고체 밀도의 결정.

파라미터 셋팅- Application : DENSITY : DEC,PLCS : 1DEC,PL (menu code 2.9.1.2)

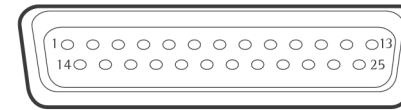
| Step | Key (or instruction) | Display/Data output |
|---------------------------|---|---------------------|
| 1. 밀도 키트를 장착 한다. | | |
| 2. 0점을 잡는다. |  | 0.0 g |
| 3. application을 시작 |  | |
| 4. "AIR" 확인한다. |  | AIR ? |
| 5. 공기중의 샘플무게를 측정한다. |  | + 20.0 g ?* |
| 6. 공기중의 무게 저장한다. |  | |
| 7. 저울에서 무게 제거한다. |  | WATER ? |
| 8. 하부계량용 홀더 장착한다. | | |
| 9. "water"를 확인한다. |  | 0.0 g ?* |
| 10. 액체에 샘플을 넣는다. |  | + 15.0 g ?* |
| 11. 결과값을 확인한다. (Enter 확인) | | + 4.0 g ?* |
| | | RhoFL 0.6237 o |
| | | Wa + 20.0 g |
| | | Wfl + 15.0 g |
| 12. 결과를 삭제 한다. |  | Rho 4.0 o |

Data interface

▶ 목 적

싸토리우스 전자저울은 PC 및 다른 장비와 연결을 할 수 있게 RS-232C 인터페이스가 장착되어 있습니다. 컴퓨터를 통해서 저울의 기능 및 응용 프로그램의 변환, 시작 또는 모니터링을 할 수 있다.

Female interface connector



Pin Assignment Chart, 25-pin, RS-232:

| | |
|---------|--|
| Pin 1: | Shield |
| Pin 2: | Data output (Tx/D) |
| Pin 3: | Data input (Rx/D) |
| Pin 4: | Internal ground (GND) |
| Pin 5: | Clear to Send (CTS) |
| Pin 6: | Not connected |
| Pin 7: | Internal ground (GND) |
| Pin 8: | Internal ground (GND) |
| Pin 9: | Not connected |
| Pin 10: | Not assigned |
| Pin 11: | + 12 V (operating voltage for Sartorius printer) |
| Pin 12: | Reset _ Out |
| Pin 13: | + 5 V |
| Pin 14: | Internal ground (GND) |
| Pin 15: | Universal remote switch |
| Pin 16: | Not connected |
| Pin 17: | Not connected |
| Pin 18: | Not connected |
| Pin 19: | Not connected |
| Pin 20: | Data Terminal Ready (DTR) |
| Pin 21: | Not connected |
| Pin 22: | Not connected |
| Pin 23: | Not connected |
| Pin 24: | Not connected |
| Pin 25: | + 5 V |

※ 더 자세한 정보를 원하시면,

"Data interface description for Entris Models"

www.Sartorius.com 에서

다운받으실 수 있습니다.

For remote switch*)

*) = Hardware restart

문제 발생 안내

Error codes are shown on the main display for approx. 2 seconds.
The program then returns automatically to the previous mode.

| Display | Cause | Solution |
|---------------------------------------|---|---|
| No segments appear on the display | No AC power is available The power supply is not plugged in | Check the AC power supply Plug in the power supply |
| HIGH | The load exceeds the balance capacity | Unload the balance |
| LOW or ERR 54 | Something is touching the weighing pan | Move the object that is touching the weighing pan |
| APP.ERR. | Cannot store data: Load on weighing pan too light or no sample on pan while application is active | Increase load |
| DIS.ERR. | Display error: Data output not compatible with output format | Change the configuration in the operating menu |
| PRT.ERR. | Interface port for printer output is blocked | Reset the menu factory settings or Contact your local Sartorius Service Center |
| ERR 02 | Calibration parameter not met, e.g.: – Press Tare to tare the balance – load on weighing pan | Calibrate only when zero is displayed Unload the balance |
| ERR 10 | The Tare key is blocked for active application programs; Only 1 tare function can be used at a time | After the tare memory has been deleted using the CF key, the Tare key can be used again |
| ERR 11 | Tara memory not allowed | Press Tare |
| The weight readout changes constantly | Unstable ambient conditions (excessive vibration or draft) A foreign object is caught between weighing pan and housing | Set up balance in another area Adjust Setup configuration Remove the foreign object |
| The weight readout is obviously wrong | The balance was not calibrated/adjusted Balance not tared before weighing | Calibrated/adjust the balance |

If any other errors occur, contact your local Sartorius Service Center.

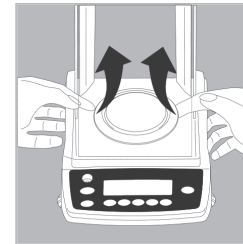
Web address: <http://www.sartorius.com>

※ 위 문제로 응급조치 후 해결 되지 않을 시 싸토리우스 코리아(주)로 연락하셔서 도움을 받으시기 바랍니다.
싸토리우스 코리아 서비스 팀 : 1600-0288

유지, 관리

▶ 서비스

문제 발생 시 수리작업은 교육을 받은 싸토리우스의 기술자가 직접 수리해야 합니다. 만약, 임의로 수리 시 발생하는 문제에 대해서는 책임을 지지 않습니다.



▶ 관리

- 전자저울의 전원을 분리한다.
(프린터 등을 연결 시 전자저울의 전원을 끈 상태에서 분리한다.)
(내부에 액체가 들어가지 않도록 주의한다.)
- 알코올등을 이용하여 적신 천과 브러쉬를 이용하여 구석구석 깔끔히 청소를 한다.
- 마른 천을 이용하여 마무리를 해준다.
※정밀저울 (0.1mg 이상)의 경우에는 팬, 팬서포트, 팬 실드를 분리 후 청소를 한다.
- 스텐레스 스틸 판 (사각의 밀판) 0.1mg
– 전자저울에서 밀판을 완전 분리 후 청소를 실시한다.
- 크리닝용 액체를 약간 묻혀서 구석구석 닦는다. (닦은 후 마른 천을 이용하여 마무리 한다.)

Overview Specifications

General Specifications

| Int. weight circuit | | All models with the designation Entris ... i-1S are equipped with an internal calibration weight. |
|---|----|---|
| Mains connection, voltage, frequency | | via Desktop power supply 6971 99, 100 – 240 Vac, ±10%, 50–60 Hz; 200 mA (max.) |
| Power consumption | VA | maximum 16; average 8 (including power supply) |
| Operating time with external battery YRB1 1Z (display backlighting on), approx. h | | 35 |

Ambient conditions

The Technical Data are valid for the following ambient conditions:

| | |
|--|------------------------------------|
| Operating temperature range | +10 ... +30 °C (+50 °F ... +86 °F) |
| Permissible ambient operating temperature | +5 ... +40 °C |
| Functionality is guaranteed up to an ambient temperature +5 °C to 40 °C. | |

| | |
|--|-----------|
| Electromagnetic Compatibility (EMC) | EN61326-1 |
| Interference emission | Class B |
| Defined immunity to interference for industrial environments | |

| Model-specific Specifications | | | | |
|---|------|--|----------------------|----------------------|
| Modele: Entris | | 224-1S 224i-1S | 124-1S 124i-1S | 64-1S 64i-1S |
| Weighing capacity | | 220 g | 120 g | 60 g |
| Readability | | 0.0001 g | 0.0001 g | 0.0001 g |
| Tare range (subtractive) | | 220 g | 120 g | 60 g |
| Repeatability (standard deviation) | < ± | 0.0001 g | 0.0001 g | 0.0001 g |
| Linearity | <± | 0.0002 g | 0.0002 g | 0.0002 g |
| Response time (average) | s | 2,5 | 2,5 | 2,5 |
| Sensitivity drift within +10 ... +30 °C | <±/K | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ |
| Adaptation to ambient conditions | | By selection of 1 of 4 optimized filter levels; display update: 0.1–0.4 (depends on filter level selected) | | |
| External calibration weight (of at least accuracy class...) | g | 200 (E2) | 100 (E2) | 50 (E2) |
| Net weight, approx. | kg | 4.4 4.8 | 4.4 4.8 | 4.4 4.8 |
| Weighing pan size | mm | 90 Ø | 90 Ø | 90 Ø |
| Whg. chamber height | mm | 230 | 230 | 230 |
| Dimensions (W×D×H) | mm | 230×303×330 | | |

| Modele: Entris | | 623-1S 623i-1S | 423-1S 423i-1S | 323-1S 323i-1S |
|---|------|---|----------------------|----------------------|
| Weighing capacity | | 620 g | 420 g | 320 g |
| Readability | | 0.001 g | 0.001 g | 0.001 g |
| Tare range (subtractive) | | 620 g | 420 g | 320 g |
| Repeatability (standard deviation) | < ± | 0.001 g | 0.001 g | 0.001 g |
| Linearity | <± | 0.002 g | 0.002 g | 0.002 g |
| Response time (average) | s | 1 | 1 | 1.1 |
| Sensitivity drift within +10 ... +30 °C | <±/K | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ |
| Adaptation to ambient conditions | | By selection of 1 of 4 optimized filter levels; display update: 0.1–0.4 (depending on the set filter level) | | |
| External calibration weight (of at least accuracy class...) | g | 500 (F1) | 200 (F1) | 200 (F1) |
| Net weight, approx. | kg | 3.2 3.6 | 3.2 3.6 | 3.2 3.6 |
| Weighing pan size | mm | 115 Ø | 115 Ø | 115 Ø |
| Dimensions (W×D×H) | mm | 230×303×136 | | |

| Modele: Entris | | 153-1S 153i-1S | 822-1S 822i-1S |
|---|------|--|----------------------|
| Weighing capacity | | 150 g | 820 g |
| Readability | | 0.01 g | 0.001 g |
| Tare range (subtractive) | | 150 g | 820 g |
| Repeatability (standard deviation) | < ± | 0.01 g | 0.001 g |
| Linearity | <± | 0.03 g | 0.002 g |
| Response time (average) | s | 1.3 | 1.3 |
| Sensitivity drift within +10 ... +30 °C | <±/K | 4 · 10 ⁻⁶ | 5 · 10 ⁻⁶ |
| Adaptation to ambient conditions | | By selection of 1 of 4 optimized filter levels; display update: 0.1–0.4 (depends on filter level selected) | |
| External calibration weight (of at least accuracy class...) | g | 100 (F1) | 500 (F2) |
| Net weight, approx. | kg | 2.6 3.0 | 2.0 2.6 |
| Weighing pan size | mm | 150 Ø | 150 Ø |
| Deviations (W×D×H) | mm | 230×303×91 | 230×303×87 |

| Modele: Entris | | 6202-1S 6202i-1S | 4202-1S 4202i-1S | 3202-1S 3202i-1S | 2202-1S 2202i-1S |
|---|------|--|----------------------|----------------------|----------------------|
| Weighing capacity | | 6200 g | 4200 g | 3200 g | 2200 g |
| Readability | | 0.01 g | 0.01 g | 0.01 g | 0.01 g |
| Tare range (subtractive) | | 6200 g | 4200 g | 3200 g | 2200 g |
| Repeatability (standard deviation) | < ± | 0.01 g | 0.01 g | 0.01 g | 0.01 g |
| Linearity | <± | 0.03 g | 0.03 g | 0.03 g | 0.03 g |
| Response time (average) | s | 1.5 | 1.5 | 1.5 | 1.5 |
| Sensitivity drift within +10 ... +30 °C | <±/K | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ | 2 · 10 ⁻⁶ |
| Adaptation to ambient conditions | | By selection of 1 of 4 optimized filter levels; display update: 0.1–0.4 (depends on filter level selected) | | | |
| External calibration weight (of at least accuracy class...) | g | 5000 (F1) | 2000 (F1) | 2000 (F1) | 2000 (F1) |
| Net weight, approx. | kg | 3.1 3.5 | 3.1 3.5 | 3.1 3.5 | 3.1 3.5 |
| Weighing pan size | mm | 180×180 | 180×180 | 180×180 | 180×180 |
| Dimensions (W×D×H) | mm | 230×303×91 | | | |


| Model: Entris | | 8201-1S 8201i-1S | 5201-1S 5201i-1S | 2201-1S 2201i-1S |
|--|------|---|-----------------------|-----------------------|
| Weighing capacity | | 8200 g | 5200 g | 2200 g |
| Readability | | 0.1 g | 0.1 g | 0.1 g |
| Tare range (subtractive) | | 8200 g | 5200 g | 2200 g |
| Repeatability (standard deviation) | < ± | 0.1 g | 0.1 g | 0.1 g |
| Linearity | <± | 0.1 g | 0.1 g | 0.1 g |
| Response time (average) | s | 1.5 | 1.5 | 1.5 |
| Sensitivity drift within +10 ... +30 °C | <±/K | 10 · 10 ⁻⁶ | 10 · 10 ⁻⁶ | 10 · 10 ⁻⁶ |
| Adaptation to ambient conditions | | By selection of 1 of 4 optimized filter levels; display update: 0.1–0.4 (depends on filter level selected) | | |
| External calibration weight (of at least accuracy class...) | g | 5000 (F2) | 5000 (F2) | 2000 (F2) |
| Net weight, approx. | kg | 2,7 3,5 | 2,7 3,5 | 2,7 3,5 |
| Weighing pan size | mm | 180 × 180 | 180 × 180 | 180 × 180 |
| Dimensions (W × D × H) | mm | 230 × 303 × 91 | | |

Accessories


| External calibration weights: | | | |
|-------------------------------|----------------|-----------------|--------------|
| For Entris balance models | Accuracy class | Weight in grams | Order no.: |
| 224 | E2 | 200 | YCW522-AC-02 |
| 124 | E2 | 100 | YCW512-AC-02 |
| 64 | E2 | 50 | YCW512-AC-02 |
| 423 | F1 | 200 | YCW523-AC-02 |
| 323 | F1 | 200 | YCW523-AC-02 |
| 153 | F1 | 100 | YCW513-AC-02 |
| 623 | F1 | 500 | YCW553-AC-02 |
| 4202 | F1 | 2000 | YCW623-AC-02 |
| 6202 | F1 | 5000 | YCW653-AC-02 |
| 3202 | F1 | 2000 | YCW623-AC-02 |
| 2202 | F1 | 2000 | YCW623-AC-02 |
| 822 | F2 | 500 | YCW554-AC-02 |
| 8201 | F2 | 5000 | YCW654-AC-02 |
| 5201 | F2 | 5000 | YCW654-AC-02 |
| 2201 | F2 | 2000 | YCW624-AC-02 |

| Product | Order No. | Product | Order No. |
|--|------------------|--|--|
| Data printer with date, time, statistics evaluation, transaction counter functions and LCD display | YDP20-OCE | Density determination kit – for Entris 224 – for Entris 124 – for Entris 64 | YDK01LP |
| Remote display, reflective (for connection to data interface port) | YRD03Z | Data cable – for connection to a PC mit USB port – for PC connection, 25-pin – for PC connection, 9-pin | YCC01-USBM2 7357312 7357314 |
| External rechargeable battery pack With battery-level indicator (LED); can be recharged using the AC adapter (charge time for completely discharged battery pack: 15 hours); see “Specifications” for hours of operation. To recharge the battery pack: Unplug the AC adapter from the balance and plug it into the battery pack | YRB11Z | Adapter cable from D-Sub 25-pin male connector to D-Sub 9-contact female connector; length: 0.25 m | 6965619 |

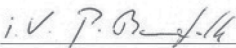
| Product | Order No. |
|---|-----------|
| Ionizing blower for eliminating static electricity | |
| – 220 V | YIB01–ODR |
| – 110 V | YIB01–OUR |
| Stat-Pen antistatic device for eliminating electrostatic charges on samples and containers (100 V bis 230 V, 50/60 Hz) | YSTP01 |
| Weighing Table | |
| – made from wood with synthetic stone | YWT09 |
| – weighing table made from synthetic stone, with vibration dampening | YWT03 |
| Wall console | YWT04 |
| Weighing pans | |
| – 1000 ml, EG 240 g, stainless steel | 641211 |
| – 500 ml | 641212 |
| – 110 ml, 90 mm Ø, aluminum | 69GP0003 |
| – 270 ml, EG 62 g, 137 mm Ø, stainless steel | YWP03G |
| – 62 mm Ø, stainless steel | 6910848 |
| – 85 ml, 70 mm Ø, aluminum | YWP06G |
| – 180 ml, 90 mm Ø, aluminum | YWP05G |
| – 174 mm d, stainless steel | YWP04G |

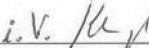


EG-/EU-Konformitätserklärung
EC / EU Declaration of Conformity



| | |
|-----------------------------------|---|
| Hersteller <i>Manufacturer</i> | Sartorius Lab Instruments GmbH & Co. KG Weender Landstrasse 94 – 108, D-37075 Goettingen, Germany |
| | erklärt in alleiniger Verantwortung, dass das Betriebsmittel <i>declares under own responsibility that the equipment</i> |
| Geräteart <i>Device type</i> | Elektronische Laborwaage <i>Electronically laboratory balance</i> |
| Baureihe <i>Type series</i> | ENTRIS.....1S |
| | in der von uns in Verkehr gebrachten Ausführung mit den grundlegenden Anforderungen der folgenden Europäischen Richtlinien übereinstimmt und die anwendbaren Anforderungen folgender harmonisierter Europäischer Normen erfüllt: <i>in the form as delivered complies with the essential requirements of the following European Directives and meets the applicable requirements of the harmonized European Standards listed below:</i> |
| 2004/108/EG 2004/108/EC | Elektromagnetische Verträglichkeit <i>Electromagnetic compatibility</i> EN 61326-1:2006 Elektrische Mess-, Steuer-, Regel- und Laborgeräte – EMV- Anforderungen – Teil 1: Allgemeine Anforderungen <i>Electrical equipment for measurement, control and laboratory use – EMC requirements – Part 1: General requirements</i> |
| 2006/95/EG 2006/95/EC | Elektrische Betriebsmittel zur Verwendung innerhalb bestimmter Spannungsgrenzen <i>Electrical equipment designed for use within certain voltage limits</i> EN 61010-1:2010 Sicherheitsbestimmungen für elektrische Mess-, Steuer-, Regel- und Laborgeräte – Teil 1: Allgemeine Anforderungen <i>Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements</i> |
| 2011/65/EU 2011/65/EU | Beschränkung der Verwendung bestimmter gefährlicher Stoffe in Elektro- und Elektronikgeräten (RoHS) <i>Restriction of the use of certain hazardous substances in electrical and electronic equipment (RoHS)</i> EN 50581:2012 Technische Dokumentation zur Beurteilung von Elektro- und Elektronikgeräten hinsichtlich der Beschränkung gefährlicher Stoffe <i>Technical documentation for the assessment of electrical and electronic products with respect to the restriction of hazardous substances</i> |
| | Jahreszahl der CE-Kennzeichenvergabe / <i>Year of the CE mark assignment:</i> 14 |
| | Sartorius Lab Instruments GmbH & Co. KG Goettingen, 2014-01-24 |


 Dr. Reinhard Baumfalk
 Vice President R&D


 Dr. Dieter Klausgrete
 Head of International Certification Management

Diese Erklärung bescheinigt die Übereinstimmung mit den genannten EG- und 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 EC and 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.

SLI13CE003-00.de,en
2011609
GOP-1.113-fo2

Sartorius Lab Instruments GmbH & Co. KG
Weender Landstrasse 94–108
37075 Goettingen, Germany
Telephone 05 51.308.0
Fax 05 51.308-32 89

www.sartorius.com

Copyright by Sartorius, Goettingen,
Germany.

No part of this publication may be reprinted
or translated in any form or by any means
without prior written from Sartorius.

All rights reserved by Sartorius in accordance
with copyright law.

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.

싸토리우스 코리아 (주)

경기도 성남시 분당구 판교역로 220 쏠리드 스페이스 8F
www.sartorius.co.kr
TEL) 031.622.5732 / Service Center) 1600-0288