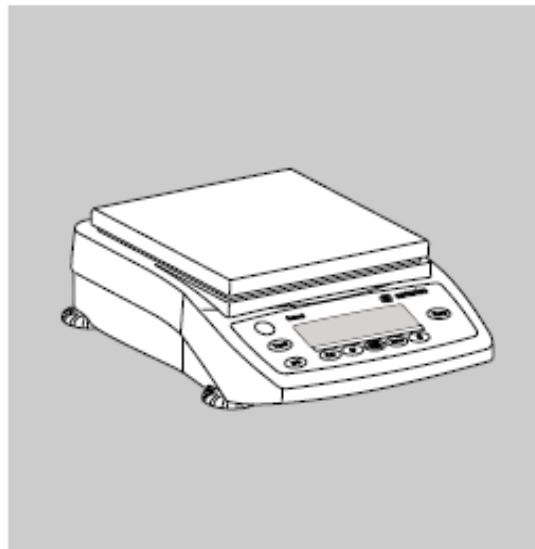
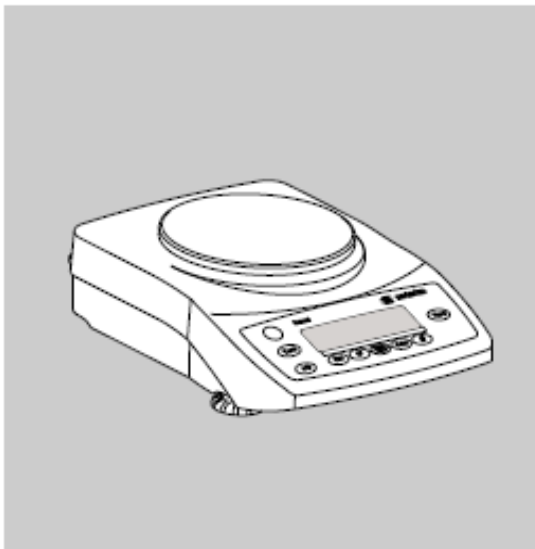
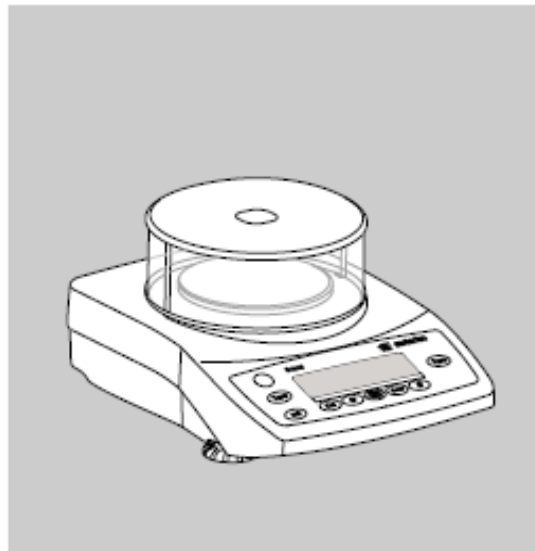
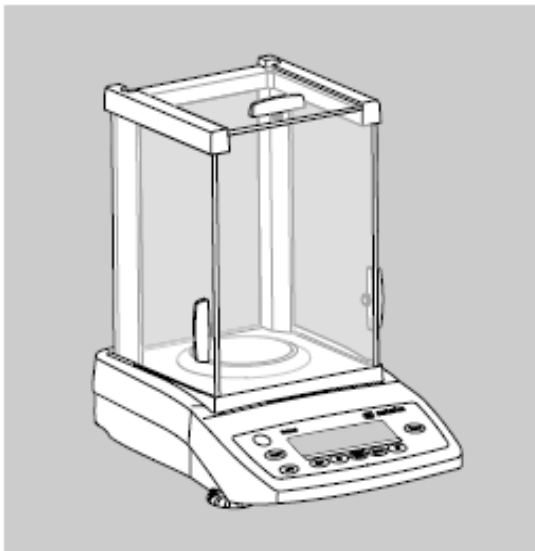


한글 설명서

Sartorius Extend Series Sartorius Gem and Gold Extend Series

Electronic Analytical and Precision Balances and Precious Metal Scales




목 차

- 설치환경 및 보관 운송
- 설치 순서
- 전원
- 워밍업및 수평조절
- 하부계량
- 디스플레이 설명
- Calibration and Adjustment (보정)
- Calibration 하는 방법
- 메뉴 설정
- 파라미터 셋팅
- ID 입력
- Counting
- Weighing in percent
- Calculation
 - Animal weighing/ Averaging
- Net-total Formulation
- Totalizing
- Mass Unit Conversion
- Density Determination
- ISO/GLP 프린팅
- Data interface
- 문제 발생
- 유지 관리
- 제품 사양
- 액세서리
- 제품보증서

Symbols

The following symbols are used in these instructions:

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

설치환경



싸토리우스의 ED Series 저울은 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 저울을 전원에 접속하기 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.

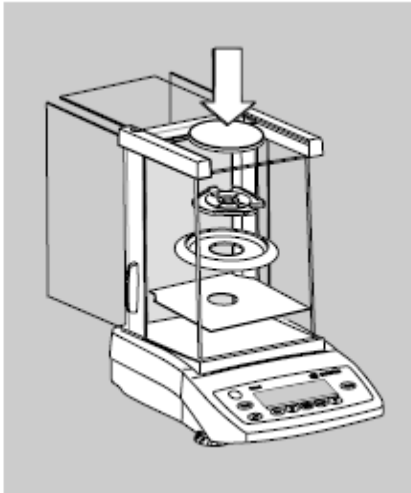
포장을 뜯은 후, 난폭한 운송에 의한 파손이 있는지 확인해 주십시오.

파손된 경우 (주)한국싸토리우스 메카트로닉스로 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,진동등을 피해 주십시오.

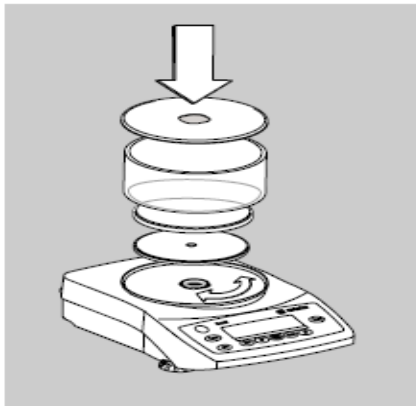
설치순서



1. 0.1 mg / 사각유리챔버 설치

아래에 순서대로 내용물을 설치하여 주십시오

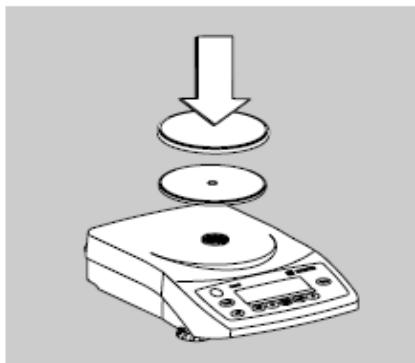
- Shield plate (바닥 판)
- Shield ring (월드링)
:ED423S-DS,ED323S-DS
ED153-DS 는 제외
- Pan support (팬 서포트)
- Weighing pan (계량 접시)



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)



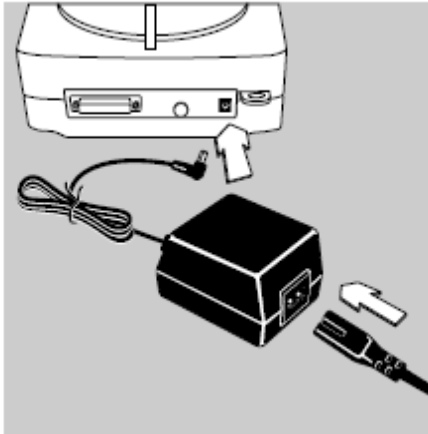
3. 삼각형팬. (원형 계량접시의 경우)

아래에 순서대로 구성품들을 설치한다.

- Pan support (팬서포트)
- Weighing pan (계량접시)

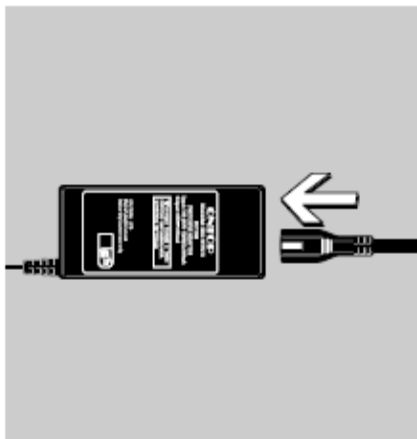
4. 사각형의 접시

- Weighing pan (계량접시)

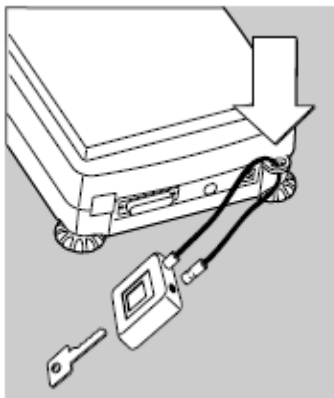


- (주)한국싸토리우스 메카트로닉스에서 공급
되어진 정품아답터만을 사용하셔야 안전하게
제품을 이용하실 수 있습니다.

- 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. 최대용량
10 kg 까지의 저울의 아답터는 Free voltage
이므로 110v or 220V에서 사용가능합니다.
단, 반드시 (주)한국싸토리우스에서 공급된 정품
아답터를 사용하셔야 합니다.



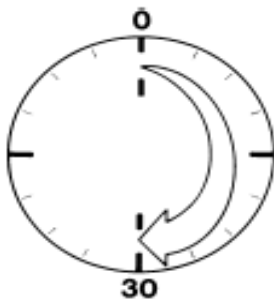
- 필요시 도난방지용 잠금장치를 체인이나
자물쇠로 잠그면 어디서라도 안전하게 사용
하실 수 있습니다.

워밍업 및 수평조절



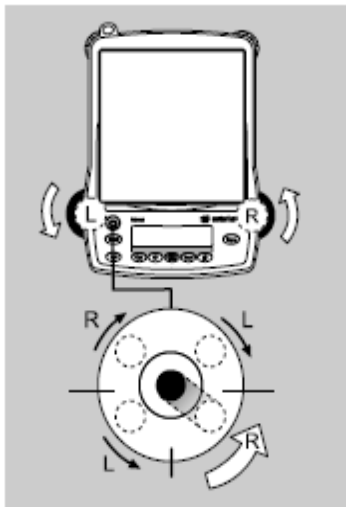
주변기기와의 접속

- Peripheral port (프린터, 컴퓨터), interface port 연결 또는 분리 전에 저울에 전원장치를 꼭 분리시키것을 확인하십시오.



워밍업

- 작동전 30분간의 워밍업은 전자저울의 정확한 값을 보증할 것이며, 30분간의 워밍업은 기기의 작동온도에 도착하게되 사용 시 전자저울에 성능을 보증할 것입니다.

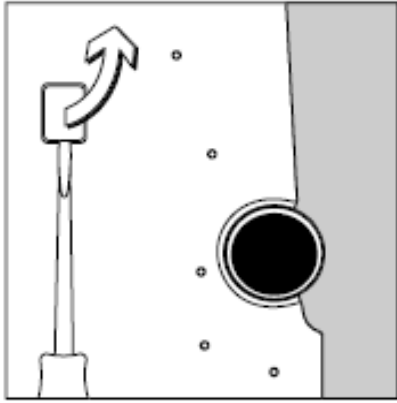


수평의 조절

사용지점에 따라 수평조절기 다리를 적당히 돌려서 수평기의 물방울이 원의 중앙에오도록 조정하면, 저울의 수평을 맞출수 있습니다.

- 물방울이 상단에 위치시: 시계방향
- 물방울이 하단에 위치시: 반시계방향

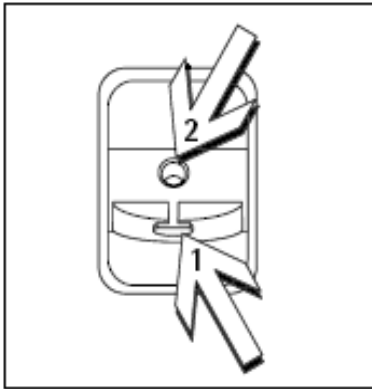
하부계량



- 하부계량
전자저울의 바닥면에 하부계량을 할 수 있는
고리가 장착되어 있다.

계량법상으로는 비공식적인 방법입니다.

- 아래쪽 커버를 열어 주십시오.
(이때 저울을 뒤집지 마십시오.)

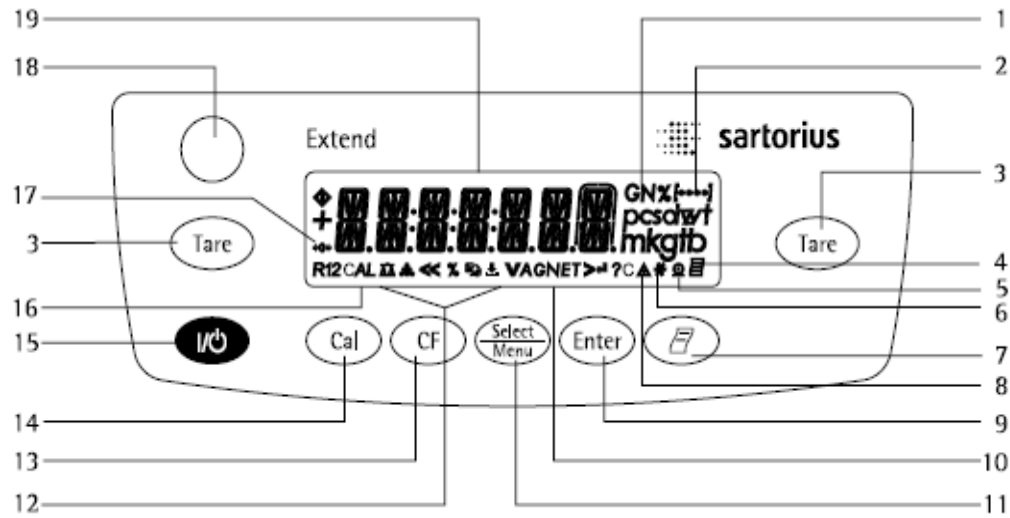


- 내장된 고리(1)의 사용:고리에 샘플을
부착할 수 있습니다. (예: 고리선 사용)

- 빈구멍(2):특별한 고리의 사용시
(ED153,ED822,ED8201,ED5201,ED2201)
제외 고리(2)를 이용시 한국싸토리우스에
문의하십시오 (T)02-575-6945

디스플레이 설명

디스플레이 전체보기.



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

Calibration and Adjustment (보정)

목 적

Calibration은 표시된 값과 실제값사이의 차이를 결정하는 것이며, Adjustments 차이값의 보정을 의미한다.

특 징

보정시 필요조건들

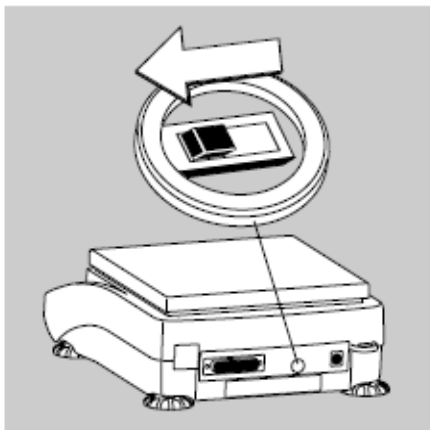
- 저울 짐판위에는 아무것도 올려 놓지 않는다.
- Tare(영점)를 잡는다.
- 0.000 g이 화면에 표시되어야 한다.
- 오차 2% 내에서만 보정가능. (예, 200g 보정시 2% 내 오차범위내 질량필요)

* 만약 위의 조건을 충족 못할시 " Error 02 " 가 나타남.

외부교정

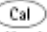
도량형법에 의해 외부 교정만으로 사용되어질 경우

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



내부교정 (0.1 mg analytical balance 와 -CW 모델만 가능)

- CAL 키를 이용하여 내부분동으로 교정한다.

Press 

CAL 키를 누른다.

내장된 분동이 자동적으로 실행된다.

저울이 보정된다.

자동적으로 weighing mode로 돌아온다.

* -CW 모델은 Internal calibration 이 가능한 모델.

Calibration 하는 방법.

내부교정.

단 계

메커니즘

1. 0점을 잡는다. (Tare 키 이용)
2. Calibration 키를 누른다.
(내부분동을 이용하여 자동적으로 동작한다.)
3. Calibration 이 실행된다.
4. 내부분동이 제거 된후 차이값이 보정된다.



외부교정

Parameters (changes in factory settings):

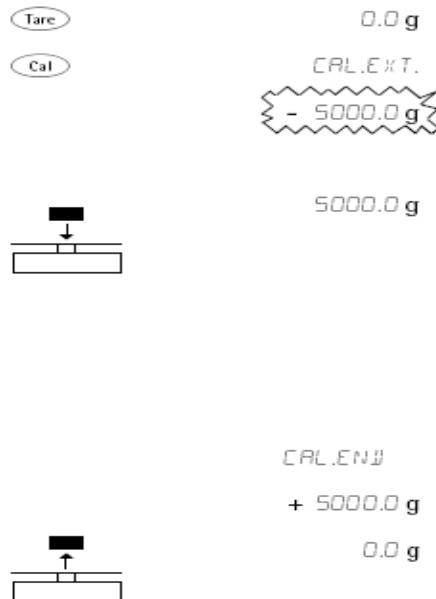
SETUP: BAL.SCAL.: CAL.-ADJ.: CAL.EXT. (menu code 1.1.9.1)

위와 같이 설정을 변경한다.

메커니즘





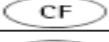

단 계

1. 0점을 잡는다. (Tare 키 이용)
2. CAL 버튼을 누른다.
(요구되어진 무게를 올린다.)
3. 표시된 분동을 저울집판위에 올린다.
4. Calibration이 실행된다.
5. 분동을 제거한다.



메뉴설정

메뉴설정을 위한 버튼 설명









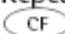


Symbol	Key
V	
>	
↵	
	 (press and hold)
<<	
<	
[....]	

위의 심볼을 순서대로 설명.


- ; 메뉴로 화면이동
- ; 하나에 메뉴의 아랫단계로 이동
- ; 메뉴설정
저장 후 그 위치에서 빠져나옴.
- ; 하나에 메뉴의 윗단계로 이동
- ; 메뉴단계의 표시.

메뉴 네비게이션 (예시)

언어선택

Key (or instruction)	Display	단 계
 (hold)	APPLIC.	1. 메뉴화면을 연다.
Repeatedly: 	INPUT *** LANGUAG.	2. 메뉴화면에 진입 후 ↓버튼을 이용하여 language를 선택한다.
	ENGLISH °	3. 원하는 언어 선택한다. (English in factory)
	ESPAÑOL	4. 스페인어 선택한다.
	ESPAÑOL °	5. Enter 키를 누르면 오른쪽 상단에 "0"이 표시된다
	LENGUA	6. CF 버튼을 이용하여 빠져나온다.
 ,  Repeatedly: 		7. CF 버튼을 반복적으로 눌러서 빠져나온다.
	0.0 g	8. 저장없이 빠져나올때는 전원버튼을 이용한다.

Parameter Settings: Menu

Level 1 [●]	Level 2 [●●]	Level 3 [●●●]	Menu code
SETUP	BAL.SCAL. Balance/scale parameters	AMBIEN Ambient conditions	1. 1. 1.
		APPFIL Application filter	1. 1. 2.
		STABRNG Stability range	1. 1. 3.
		TARNG Taring ¹⁾	1. 1. 5.
		AUTOZER Auto zero	1. 1. 6.
		WTUNIT Basic weight unit	1. 1. 7.
		DISPLA Display accuracy ¹⁾	1. 1. 8.
		CALADJ Function of the  key	1. 1. 9.
		CALUNIT Weight unit for calibration ¹⁾	1. 1. 11.
	INTERF. Interface	BAUD Baud rate	1. 5. 1.
		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.
	PRNTOUT Settings for print function	DATA REC. Output: SBI (ASCII) or printout	1. 5. 6.
		PRNT (manual/automatic)	1. 6. 1.
		STOPAUT Stop automatic printing	1. 6. 2.
		AUTOZCL Time-dependent autom. printing	1. 6. 3.
		TAR/PR Tare bal/scale after ind. print	1. 6. 4.
		PRTUNIT Printout of appl. parameters	1. 6. 5.
		FORMAT Line format for printout	1. 6. 6.
		GLP ISO/GLP-compliant printout	1. 6. 7.
		TIME: 12/24 h	1. 6. 8.
	EXTRAS Additional functions	DATE: Format	1. 6. 9.
		MENU	1. 8. 1.
		SIGNAL Acoustic signal (beep)	1. 8. 2.
		KEYS Keypad	1. 8. 3.
		EXTKEY External switch function	1. 8. 4.
		ONMODE Power-on mode	1. 8. 5.
	RESET	BACKLIT Display backlighting	1. 8. 6.
		MENU Factory settings	1. 9. 1.
APPLIC. Application programs	WEIGH		2. 1.
	UNIT Toggle wt. unit	DISP.WIG. Display accuracy ¹⁾	2. 2. 2.
	COUNT Counting	RESOLU Resolution	2. 3. 1.
		REFUPBIT Autom. ref. sample updating	2. 3. 2.
	PERCENT Weighing in percent	DEC.PLCS Decimal places	2. 4. 1.
	NET-TOT Net-total formulation	COMP.PRT Printout of components	2. 5. 1.
	TOTAL Totalizing	COMP.PRT Printout of components	2. 6. 1.
	ANIMALW Animal weighing	ACTIVITY Animal activity	2. 7. 1.
		STAR	2. 7. 2.
	CALC. Calculation	METHOD (operator)	2. 8. 1.
WEIGHT Density determination		DEC.PLCS Decimal places	2. 8. 2.
		DEC.PLCS Decimal places	2. 9. 1.
INPUT Input	IDNO	ID input; max. 7 characters	3. 1.
INFO Information	VERSION, SER.NO., MODEL	Display software ver., serial no., model	4. 1./2./3.
LANGUAG.	ENGLISH (factory setting)		5. 1.
	DEUTSCH (German)		5. 2.
	FRANÇ. (French)		5. 3.
	ITAL. (Italian)		5. 4.
	ESPAÑOL (Spanish)		5. 5.
	РУССКИЙ (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.	AMBIENT	✓ STABLE Very stable	1. 1. 1. 1
			o STABLE	1. 1. 1. 2
			UNSTABLE	1. 1. 1. 3
			✓ UNSTABLE Very unstable	1. 1. 1. 4
		APP.FILT.	o FINAL R. Final readout mode	1. 1. 2. 1
			FILLING Filling mode	1. 1. 2. 2
		STAB.RNG.	1/4 SIG. (digit)	1. 1. 3. 1
			1/2 SIG.	1. 1. 3. 2
			1 - SIGIT	1. 1. 3. 3
			o 2 - SIGIT	1. 1. 3. 4
			4 - SIGIT	1. 1. 3. 5
			8 - SIGIT)	1. 1. 3. 6
		TARING ¹⁾	W/O STAB (w/o stability)	1. 1. 5. 1
			o W/ STAB After stability)	1. 1. 5. 2
		AUTOZER.	OFF	1. 1. 6. 1
			o ON	1. 1. 6. 2
		WT.UNIT	For list of units, see	1. 1. 7. 1
			"Toggling between Weight Units"	1. 1. 7. 23
		DISP.SIG.)	o ALL	1. 1. 8. 1
			MINUS 1	1. 1. 8. 2
			SIG. 1 1 interval	1. 1. 8. 6
		CAL./ADJ.	o CAL. EXT. External cal./adj. ¹⁾	1. 1. 9. 1
			CAL. INT. Internal cal./adj. ²⁾	1. 1. 9. 2
			LOCKED Cal key blocked	1. 1. 3. 3
		CAL.UNIT. Unit ¹⁾	o GRAMS	1. 1.11. 1
			KILOG. Kilograms	1. 1.11. 2
			POUNDS	1. 1.11. 3

¹⁾ Setting cannot be changed on verified balances/scales

²⁾ Only on models with built-in motorized calibration weight


Level 1

Level 2

Level 3


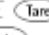




Level 4

Menu code

Level 1 [•]	Level 2 [••]	Level 3 [•••]	Level 4 [••••]	Menu code
SETUP	INTERF. Interface	BAUD rate	600 o 1200 2400 4800 9600 19200	1. 5. 1. 3 1. 5. 1. 4 1. 5. 1. 5 1. 5. 1. 6 1. 5. 1. 7 1. 5. 1. 8
		PARITY Parity	o 0 EVEN NONE	1. 5. 2. 3 1. 5. 2. 4 1. 5. 2. 5
		STOP BIT No. of stop bits	o 1 BIT 2 BITS	1. 5. 3. 1 1. 5. 3. 2
		HANDSHK. Handshake mode	SOFTWARE o HARDWARE NONE	1. 5. 4. 1 1. 5. 4. 2 1. 5. 4. 3
		DATA BIT No. of data bits	o 7 BITS 8 BITS	1. 5. 5. 1 1. 5. 5. 2
		DAT.REC. Com- munication mode	5 BIT (ASCII) ¹⁾ o PRINTER (GLP-printout)	1. 5. 6. 1 1. 5. 6. 2
	PRNT. OUT Printing fct.	PRINT (manual/ automatic)	MAN. W/O W/o stability o MAN. WITH W/ stability AUT. W/O Autom. w/o stability AUT. WITH. Autom. w/ stability	1. 6. 1. 1 1. 6. 1. 2 1. 6. 1. 3 1. 6. 1. 4
		STOPAUT. Stop automatic printing	o OFF Not possible ON Use print key 	1. 6. 2. 1 1. 6. 2. 2
		AUT.CYCL. Time-dependent autom. printing	o EACH/PL (1 display update) AFTER 2 (2 display updates)	1. 6. 3. 1 1. 6. 3. 2
		TARE/PRNT. Tare the bal./scale after individual printout	o OFF ON	1. 6. 4. 1 1. 6. 4. 2

¹⁾ Note concerning verified balances/scales as legal measuring instruments in the EU*:
 In the setting "SBI", the non-verified display digit is not automatically identified.
 Please take the corresponding measures or adjust the settings on the peripheral device.

* Including the signatories of the Agreement on the European Economic Area.

[•]	[••]	[•••]	[••••]	
SETUP	PRT. OUT Printing fct.	PRT. INIT. Printing appli- cation parameters	OFF 1. 6. 5. 1 o ALL All parameters 1. 6. 5. 2 MAINPAR. Main parameters 1. 6. 5. 2	
		FORMAT Line format for printout	16 CHAR. 16 characters (w/o ID) 1. 6. 6. 1 o 22 CHAR. 22 characters (w/ ID) 1. 6. 6. 2	
		GLP Printout as ISO/GLP- compliant	o OFF 1. 6. 7. 1 CAL.-ADJ. Only for calib./adj. 1. 6. 7. 2 ALWAYS All printouts 1. 6. 7. 3	
		TIME	24H 24-hour format 1. 6. 8. 1 12H 12-hour format "AM/PM" 1. 6. 8. 2	
		DATE	JJ.MM.YY Day/month/year 1. 6. 9. 1 MM.JJ.YY Month/day/year 1. 6. 9. 2	
	EXTRAS Additional functions	MENU	o CANE. IT Can change settings 1. 8. 1. 1 RJ. ONLY 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 LOCKEJ 1. 8. 3. 2	
		EXT. KEY Function of the external switch	o PRINT  1. 8. 4. 1 Z/TARE  1. 8. 4. 2 CAL.  1. 8. 4. 3 SELECT  1. 8. 4. 4 CF  1. 8. 4. 5 ENTER  1. 8. 4. 6	
		ONMODE Power-on mode	o OFF / ON Off/on/standby 1. 8. 5. 1 STANDBY On/standby 1. 8. 5. 2 AUTO ON Auto on 1. 8. 5. 3	
		BACKLIT Display backlighting	OFF 1. 8. 6. 1 o ON 1. 8. 6. 2	
	RESET Reset menu	MENU Factory settings	YES Restore fcty. settings 1. 9. 1. 1 o NO Do not restore settings 1. 9. 1. 2	

Level 1
[•]

Level 2
[••]

Level 3
[•••]

Level 4
[••••]

Menu code

APPLIC.	WEIGH	UNIT	o	ALL	2. 1.
Applic. programs	Toggle units	Display accuracy	o	MINUS 1	2. 2. 2. 1
				1/15. 1 interval	2. 2. 2. 2
					2. 2. 2. 6
	COUNTING	RESOLUT.	o	1/15. ACC. Display accuracy	2. 3. 1. 1
				10-FOL 10 times > disp.	2. 3. 1. 2
		REF. UPDT.	o	OFF	2. 3. 2. 1
				AUTO	2. 3. 2. 2
	PERCENT ²⁾ Weighing in percent	DEC.PLCS Decimal places	o	NONE No dec. places	2. 4. 1. 1
				1 DEC.PL. 1 decimal place	2. 4. 1. 2
				2 DEC.PL. 2 decimal places	2. 4. 1. 3
				3 DEC.PL. 3 decimal places	2. 4. 1. 4
	NET-TOT Net-total	COMP.PRT. Component printout	o	OFF	2. 5. 1. 1
				ON	2. 5. 1. 2
	TOTAL ²⁾ Totalizing	COMP.PRT. Component printout	o	OFF	2. 6. 1. 1
				ON	2. 6. 1. 2
	ANIMALW. ²⁾ Animal weighing	ACTIVITY. Animal activity	o	CALM Fluct.: 2% of test obj.)	2. 7. 1. 1
				ACTIVE (fluct.: 5% of test obj.)	2. 7. 1. 2
				2 ACTIVE (fluct.: 20% of test obj.)	2. 7. 1. 3
		START	o	MANUAL	2. 7. 2. 1
				AUTO. Automatic	2. 7. 2. 2
	CALC. ²⁾ Calculation	METHO. (operator)	o	MUL. Multiplier	2. 8. 1. 1
				1/1. Divisor	2. 8. 1. 2
		DEC.PLCS Decimal places	o	NONE No dec. places	2. 8. 2. 1
				1 DEC.PL. 1 decimal place	2. 8. 2. 2
				2 DEC.PL. 2 decimal places	2. 8. 2. 3
				3 DEC.PL. 3 decimal places	2. 8. 2. 4
	DENSITY ²⁾ Density determination	DEC.PLCS Decimal places	o	NONE No dec. places	2. 9. 1. 1
				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

¹⁾ Setting cannot be changed on verified balances/scales

²⁾ not on ED...-PCE models

ISO/GLP 프린팅을 위한 ID 입력

Level 1

Level 2

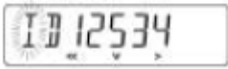



Level 2

Menu code


LEVEL 1 [.]]	LEVEL 2 [.]]	LEVEL 3 [.]]	MENU CODE
INPUT Input	ID NO.	ID input; max. 7 characters Permitted characters: 0 to 9; A to Z; (hold) ; space	3. 1.

Input – Enter – ID No. – Enter – 7가지의 영어문자와 숫자등입력가능

ID를 입력하는 방법은 아래와 같이 실시한다.

key: Press and hold to repeat			
Display	Key	Display symbol	Function
	First position:		
	Enter	>	Go to next position
	Select Menu	V	Select current position
	CF	<<	Exit without saving changes
	Middle positions:		
	Select Menu	V	Select current position
	Enter	>	Go to next position
	CF	<	Go to previous position
	Last position:		
	Select Menu	V	Select current position
	CF	<	Go to previous position
	Enter	↵	Store and exit




Counting

Display symbol: 

목 적

동등한 무게의 여러 샘플의 많은 양의 개수를 정확하게 파악할 수 있다.
 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. (1 에서 100까지) (정확한 개수 측량을 위해서는 많은 수 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 질때 수행된다. (디스플레이상에 OPT 가 표시됨)

준비단계

메뉴안 Counting application 선택한다.
 (아래의 그림은 factory setting 된 모습)

APPLIC. Application program

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

o = Factory setting

프린팅된 후

Printout: Counting

nRef	+	10	: Reference sample quantity
wRef	+	21.14 g	: Reference weight
Qnt	+	500 pcs	: Calculated quantity

(기준 1 개 의 무게가 21.14 g)

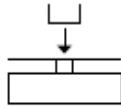
예제 : 같은 샘플의 개수 측정

파라미터 셋팅 : application : count (menu code 2.3)

Step

Key (or instruction) Display/ Data output

1. 빈 용기를 저울위에 올려놓는다.



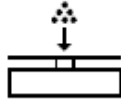
+ 22.6 g

2. 0점을 잡는다.
(Tare 키를 이용)

Tare

0.0 g

3. 기준 샘플을 올린다.
(예: 20개 정도)



4. 비교 샘플의 양을 변화시키
원하면 다음키를 누른다.

Select Menu

REF 10 pcs

5. 원하는 기준 샘플 개수를
선택한다.

Repeatedly: Select Menu
Press briefly
) Select Menu press
and hold

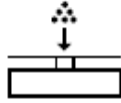
REF 20 pcs

6. 비교샘플의 개수와
비교 무게 확인한다.

Enter

+ 20 pcs *
nRef 20 pcs
wRef 1.07 g

7. 측정하고자 하는 개수 만큼
샘플을 저울에 올려 놓는다.



+ 500 pcs

8. 프린터 출력 원하면, 프린터
를 누른다.

Printer

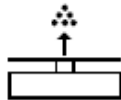
Qnt + 500 pcs

9. 총무게와 개수를 버튼을 이
하여 볼 수 있다.

Repeatedly: Select Menu

1.07 g Δ *
+ 535.0 g *
+ 500 pcs *

10. 샘플 제거



- 2 pcs *

11. 반복적으로 원하면 step 7
부터 다시 측정한다.

12. 기준값을 삭제한다.

CF

0.0 g

Weighing in Percent

Display symbol: %

목 적

이 기능은 기준 중량을 비율로 변환하여 확인하고자 할때
사용되어지는 기능

기준 샘플의 변화 방법

키버튼



key

내 용

1. 키를 누른다.
2. 기준 샘플의 개수를 입력한다. (1 에서 100까지)
(정확한 개수 측량을 위해서는 많은 수 입력한다.)
3. 키를 이용하여 원하는 개수를 선택한다.
4. 선택 후 길게 키를 길게 누르고 있다. (입력 완료)



key



key

준비단계

메뉴의 Weighing in percent application 선택한다.
(아래의 그림은 factory setting 된 모습)

```
APPLIC. Application program
├── PERCENT Weighing in percent1)
│   ├── DEC.PLCS. Decimal places
│   │   ├── NONE No decimal places
│   │   ├── o 1 DEC.PL. 1 decimal place
│   │   ├── 2 DEC.PL. 2 decimal places
│   │   └── 3 DEC.PL. 3 decimal places
```

o = Factory setting

¹⁾ not on ED...-PCE models

Printout: Weighing in percent

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



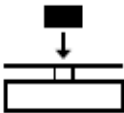

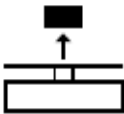
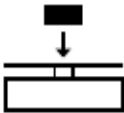




프린팅된 모습

(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. 0 점을 잡는다.		0.0 g
2. 설정된 기준 퍼센트 확인		REF 100 %
3. 100% 에 해당하는 무게의 비교샘플을 올려놓는다.		
4. enter 키를 이용하여 확인한다.		+ 100.0 % * pRef 100 % Wxx% 111.6 g
5. 기준 샘플을 제거한다.		
6. 모르는 무게의 샘플을 올려놓는다.		+ 94.9 % *
7. 원하는 퍼센트가 나타남.		Pr c + 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. 필요시 프린터 출력		N + 105.9 g



Calculation

Display symbol: C

목 적

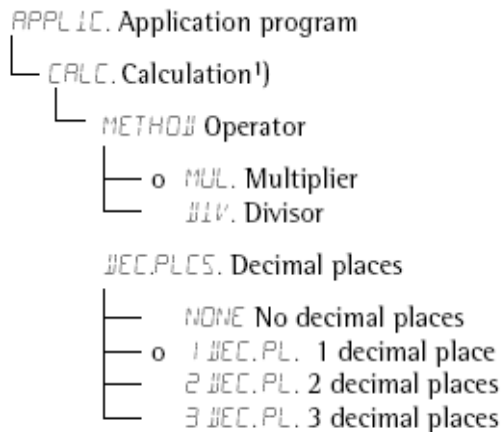
이 프로그램을 가지고 무게 값을 계산할 수 있다 (곱셈 나눗셈을 이용)
 단위 면적당 무게를 계산 할 수 있으며, 종이의 무게도 (grams per square meter)
 계산 할 수있다.

계산 인자의 셋팅.

- | 키버튼 | 내 용 |
|---|--|
|  key | 1. 키를 누른다. (7 digits 까지의 숫자 선택)
2. 필요시 소수점 자리 입력 (0.000001 에서 9999999) |
|  key | 3. 선택 후 길게 키를 길게 누르고 있다. (입력 완료) |

준비단계

메뉴안에 Calculation application 선택 한다.
 (아래의 그림은 factory setting 된 모습)



☒ = Factory setting

¹⁾ not on ED...-PCE models

Printout: Calculation









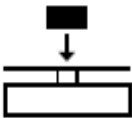


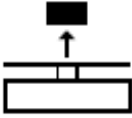
Mul	+	1.2634	: Multiplier
Div	+	0.6237	: Divisor
Res	+	79.7 <input checked="" type="radio"/>	: Result

예제


종이의 단위면적당의 무게 계산 (예제에서는 A4용지 사용)
 $0.210\text{m} \times 0.297\text{m} = 0.06237\text{m}^2$
 단위 지역당 무게 결정을 위해서는 전체무게에서 표면적을 나눈다.

파라미터 셋팅

Application : Calculation : Method : Div (2.8.1.2)

Step	Key (or instruction)	Display/Data output
1. 0점을 잡는다.		0.00 g
2. 버튼을 눌러 활성화 시킨다		-----0.
3. 나눗셈인자를 입력 (보기에서는 0.06237)):  , 5x  , 2x  , Repeatedly or press and hold;  ,  , etc.	---00000 ---06000 ---06237
4. enter 버튼을 이용 나눗셈인자 저장시킨다.		+ 0.0 ° Div 0.6237
5. 단위면적당 무게 측정 (저울위에 A4 용지 놓는다.)		+ 79.7 ° *
6. 프린터 원하면 출력		Res + 79.7 °
7. 무게와 계산된 값 볼수 있음	Repeatedly: 	+ 4.97 g * + 79.7 ° *
8. 무게 제거 9. 필요시 step 5 돌아가 다시 반복		+ 0.0 ° *




Animal weighing / Averaging

Display symbol: 

목 적

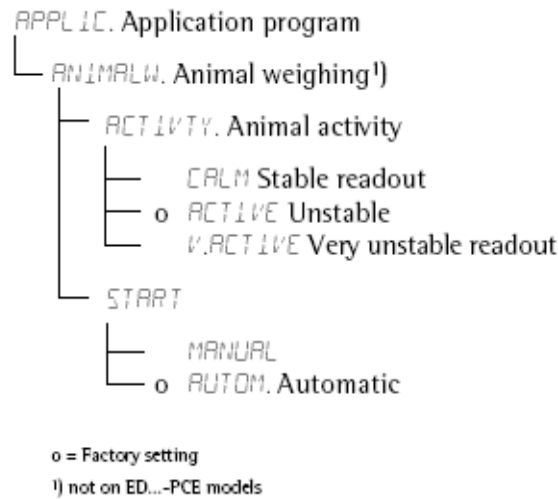
불안정한 샘플 측정시 이용 및 불안정한 환경에서 측정시 사용한다.
또, 여러가지의 샘플의 평균값을 이 프로그램을 통해 계산할 수 있다.

인자의 변환

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 application setting 한다.
(아래의 그림은 factory setting 된 모습)



Printout: Animal weighing

```

mDef      20      : Number of sub-
                  weighing operations
x-Net + 410.1 g   : Calculated average

```

프린팅 된 모습
(20 번 측정함)

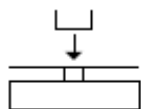
(20번 측정시의 평균무게)

예제

자동 시작이 설정된 20번 측정의 동물계량
파라미터 셋팅 : Application : animal weighing (menu code 2.7.)

Step	Key (or instruction)	Display/Data output
------	----------------------	---------------------

1. 저울위에 동물계량용
접시 놓는다.



22.6 g

2. 0점을 잡는다.

Tare

0.0 g

3. 버튼을 눌러 횟수 확인

Select Menu

REF 30

4. 원하는 횟수로 입력시킨다.
(20 입력)

Repeatedly:
Press briefly
Select Menu press
and hold

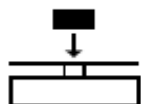
REF 20

5. enter 키를 이용 저장.

Enter

+ 0.0 g *

6. 동물계량용 접시를 올려
놓은 후 움직이는 샘플을
측정한다.



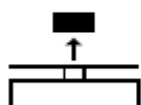
888
20
19
....
1

7. 횟수되로 계산된 값이
나타난다.

+ 410.1 g Δ*

mDef 20
x-Net + 410.1 g

8. 샘플제거 후
원하면 다음 샘플을 측정한다.



+ 0.0 g *

Net-total Formulation

Display symbol: ↓

목 적

정해진 양 까지 다른 구성성분의 값을 측정 할 수있다.
프린트시 전체값과 구성요소의 각각의 값을 둘다 출력할 수있다.

특 징

- 0에서 99 개의 구성 성분 까지 측정 가능.
- 구성 무게들 저장 (값이 저장후에 자동적으로 Tare , 자동 출력가능)
- CF 버튼을 누르면 구성성분의 무게가 지워지며, 전체의 무게만 남음.
- Toggling 기능 이용시

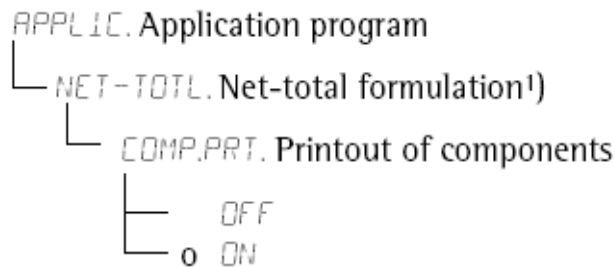


(< 2 sec).

전체값과 구성성분의 값 서로 변환 가능.

준비 단계

Net- total application 기능 설정.
(아래의 그림은 factory setting 된 모습)



o = Factory setting

¹⁾ Factory setting on ED...-PCE models

Printout: Net-total formulation

Comp 2+ 278.1 g	: Second component
T-Comp+ 2117.5 g	: Sum of components

프린팅시 모습

2번째 구성성분의 무게
전체의 구성성분의 무게.

예제

상자안에 구성물질의 개수

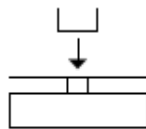
파라미터 셋팅.

application : NET-TOT (menu code 2.5)

Step

Key (or instruction) Display/Data output

1. 빈용의 무게 측정



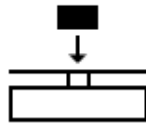
65.0 g

2. 0점을 잡는다.



0.0 g

3. 첫번째 물질을 넣는다.



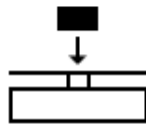
+ 120.5 g

4. enter 버튼을 이용 저장.



+ 0.0 g * NET
Comp 1+ 120.5 g

5. 두번째 물질을 넣는다.



+ 70.5 g * NET

6. enter 이용 저장.



+ 0.0 g * NET
Comp 2+ 70.5 g

7. step 5.6 을 이용 더 많은
샘플을 넣는다.

Repeat steps
5 and 6

8. 버튼을 이용하여 현재의
총 무게를 알 수 있다.




+ 191.0 g *

9. 전체무게 프린팅
(각각의 구성성분은 제거
된 상태임)



+ 2117.5 g
T-Comp+ 2117.5 g

Totalizing

Display symbol: 
목 적

이 기능을 가지고 전자저울에 최대 용량을 넘지 않는 범위 내에서
연속적이고 상호 독립적인 무게값을 더할 수 있다.

특 징

- 99 값 까지 기억할 수 있다.
- 구성 성분의 무게들을 저장할 수 있다.
- Toggle 기능을 이용하여 현재의 각각의 값과 기억된 전체의 값을 볼 수있다.
- 각각의 구성성분의 무게의 전체를 프린팅 할 수있다.
- 기능을 종료시킴으로서 전체무게를 프린팅 할 수 있다.

준비 단계

메뉴안에서 Totalizing application 을 선택한다.
(아래의 그림은 factory setting 된 모습)

```
APPLIC. Application program
├─ TOTAL Totalizing¹)
│   └─ COMP.PRT. Printout of components
│       └─ OFF
│           └─ o ON
```

o = Factory setting

¹) not on ED...-PCE models

Printout: Totalizing

```
Comp 2+ 278.1 g : Second component
S-Comp+ 2117.5 g : Totalizing memory
```

예 제

무게 값의 합계기능을 이용
파라미터 셋팅

APPLIC : TOTAL : COMP.PRT: ON(menu code 2.6.2.1)

Step

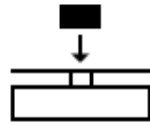
Key (or instruction) Display/Data output

1. 0점을 잡는다.

Tare

0.0 g

2. 샘플을 올려 놓는다.
(380g)



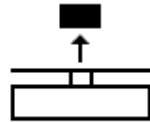
+ 380.0 g

3. 값을 저장한다. (enter 키 이용)

Enter

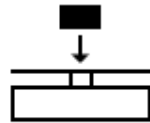
+ 380.0 g *
Comp 1+ 380.0 g

4. 샘플제거



+ 0.0 g *

5. 다른 샘플을 올려놓는다.
(575g)



+ 575.0 g *

6. 값을 저장한다 (enter 키 이용)

Enter

+ 955.0 g *
+ 575.0 g *
Comp 2+ 575.0 g

7. 저장된 합계의 값을 볼 때

Select Menu

+ 955.0 g Δ*

8. 더 많은 샘플 측량시
step 5, 6 실시

Repeat steps
5 and 6

9. 전체값 프린팅
구성성분을 지울 때 (CF 키 사용)

CF

0.0 g
S-Comp+ 2117.5 g

Mass Unit Conversion

목 적

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


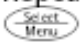


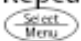


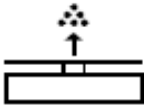
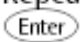
특 징

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

예 제

pound (lb), ounces (ozt) 로 변환

파라미터 셋팅

Step	Key (or instruction)	Display/Data output
1. 옆의 버튼을 누른다.		NONE 0 [•]
2. 버튼을 이용하여 원하는 것을 설정한다.	Repeatedly: 	POUNDS
3. enter 키를 이용 입력		POUNDS 0
4. 두번째 단위를 찾는다. (TROY OZ)	 , Repeatedly: 	NONE 0 [••] TROY OZ.
5. enter 키를 이용 입력한다.		TROY OZ. 0
6. 또 다른 단위의 설정을 원하면 위와 같이 입력한다.		[•••]
7. CF 키를 눌러 빠져나온다.		0.00 g
8. 샘플을 올려 놓는다.		+ 100.00 g
9. Toggling 을 이용 단위변환가 (enter 키 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 목적사용시 mg, g , kg cat만 사용가능. 연구용 목적시 제외)

Menu item	Unit	Conversion	Display
-----------	------	------------	---------

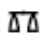
		factor	symbol
1) USER DEF. ¹⁾	Grams	1.000000000000	o
2) GRAMS (Factory setting)	Grams	1.000000000000	g
3) KILOGR.	Kilograms	0.001000000000	kg
4) CARATS	Carats	5.000000000000	o
5) POUNDS	Pounds	0.00220462260	lb
6) OUNCES	Ounces	0.03527396200	oz
7) TROY OZ.	Troy ounces	0.03215074700	ozt
8) HK TAEI	Hong Kong tael	0.02671725000	tl
9) SING.TAEI.	Singapore tael	0.02645544638	tl
10) TWN.TAEI	Taiwanese tael	0.02666666000	tl
11) GRAINS	Grains	15.4323583500	GN
12) PENNY.WT.	Pennyweights	0.64301493100	dwt
13) MILLIGR.	Milligrams	1000.000000000	mg
14) PT.P.L.B.	Parts per pound	1.12876677120	o
15) CHN.TAEI	Chinese tael	0.02645547175	tl
16) MOMMES	mommies	0.26670000000	m
17) AUSTRICT.	Austrian carats	5.00000000000	Kt
18) TOLA	Tola	0.08573333810	o
19) BAHT	Baht	0.06578947436	b
20) MESGHAL	Mesghal	0.21700000000	o
21) TONS	Tons	0.00000100000	t
22) LB / OZ ²⁾	Pounds : ounces	0.03527396200	lb oz
23) NEWTON	Newton	0.00980665000	N

¹⁾ = User-defined weight unit; can be loaded in the balance/scale over an optional RS-232 or USB interface using a computer program.

²⁾ = The format for display of pounds/ounces cannot be changed: xx.yy.yy x=lb, y=oz

 Some weight units may be blocked from use in legal metrology, depending on national verification laws.

Density Determination

Display symbol: 

목 적

이 기능을 부력법을 이용하여 고체의 밀도를 측정할 수 있다.

측정을 위해서는 홀더와 매달림 와이어를 사용하여야 측정할 수가 있으며,




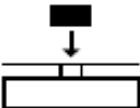

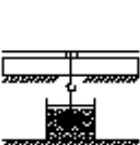



저울에는 포함되어 있지 않습니다.

파라미터 셋팅

Application : DENSITY : DEC.PLCS : 1DEC.PL (menu code 2.9.1.2)

예제

고체 밀도의 결정.

Step	Key (or instruction)	Display/Data output
1. 밀도키트 장착		
2. 0점을 잡는다.		0.0 g
3. application 시작		
4. "AIR " 확인한다.		AIR ?
5. 공기중의 샘플무게를 측정		+ 20.0 g *
6. 공기중의 무게 저장		
7. 저울에서 무게 제거		WATER ?
8. 하부계량용 홀더 장착		
9. "water " 확인한다.		0.0 g *
10. 액체에 샘플을 넣는다		+ 15.0 g *
11. 결과값을 확인. (enter 확인)		+ 4.0 ° *
		Wa + 20.0 g
		WfL + 15.0 g
		Rho 4.0 °
12. 결과 삭제.		

ISO/ GLP – Compliant Printout/ Record

특 징

제품정보, ID text, 날짜, 시간 등을 용지상에 프린팅 할 수 있다.

저울 값 사이에서 (GLP header 와 GLP footer 사이)

다음 파라미터를 포함하여.

GLP header :

- Date
- Time at beginning of measurement
- Balance/Scale manufacturer
- Balance/ Scale model
- Serial number
- Software version
- Identification number of the current sampling operation

GLP footer

- Date
- Time at end of measurement
- Field for operator signature

ISO/GLP 출력

싸토리우스 전자저울에서 ISO/GLP 문서를 출력하기 위해서는 YDP03-OCE 프린터 연결을 해야한다.

설 정

출력을 위한 메뉴코드 설정. (configuration 참조)

- ISO/GLP -compliant 출력 (보정할 때)

SETUP: PRNT.OUT:GLP:CAL.-ADJ. (menu code 1.6.7.2)

ISO/GLP -compliant 출력 (항 시)

SETUP: PRNT.OUT:GLP:ALWAYS ON (code 1.6.7.3)

- 출력 형식 (22 문자 : 출고시)

SET:PRNT.OUT :FORMAT:22CHAR. (menu code 1.6.6.2)

- 시간 설정

SET:PRNT.OUT : TIME:24H (menu code 1.6.8.1)

또는

SET:PRNT.OUT : TIME:12H (menu code 1.6.8.2)

- 날짜 설정

SET:PRNT.OUT : DATE:DD.MMM.YY (menu code 1.6.9.1)

SET:PRNT.OUT : DATE:MMM.DD.YY (menu code 1.6.9.2)

ISO/ GLP- Compliant 출력을 원하지 않을 때

SETUP:PRNT.OUT PRINT: AUT.W/O 또는 AUT.WITH

(menu code 1.6.1.3.1.6.1.4)

16 문자로 출력 원할 때 (1.6.6.1)

기능 키



출력을 하고자 할 때



application 동작할 때 자동적으로 데이터 출력 시



GLP footer 로 보내기

application 에서 빠져 나오기.

The ISO/GLP-compliant printout can contain the following lines:

-----	Dotted line
17-Aug-2005 10:15	Date/time (beginning of measurement)
SARTORIUS	Balance/scale manufacturer
Mod. ED8201	Model
Ser. no. 10105355	Balance/scale serial number
Ver. no. 00-32-02	Software version
ID 2690 923	ID.
-----	Dotted line
L ID	Measurement series no.
nRef 10 pcs	Counting: reference sample quantity
wRef 21.14 g	Counting: reference weight
Qnt + 567 pcs	Counting result
-----	Dotted line
17-Aug-2005 10:20	Date/time (end of measurement)
Name:	Field for operator signature
-----	Blank line
-----	Dotted line

ISO/GLP-compliant printout for external calibration/adjustment:

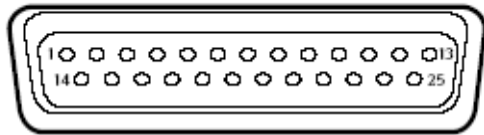
-----	Dotted line
17-Aug-2005 10:30	Date/time (beginning of measurement)
SARTORIUS	Balance/scale manufacturer
Mod. ED8201	Model
Ser. no. 10105352	Balance/scale serial number
Ver. no. 00-32-02	Software version
ID 2690 923	ID.
-----	Dotted line
Cal. Ext. Test	Calibration/adjustment mode
Set + 5000.0 g	Calibration weight
Diff. + 0.2 g	Difference determined in calibration
Cal. Ext. Complete	Confirmation of completed calibration procedure
Diff. 0.0 g	Difference from target following adjustment
-----	Dotted line
17-Aug-2005 10:32	Date/time (end of measurement)
Name:	Field for operator signature
-----	Blank line
-----	Dotted line

Data interface

목 적

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

Female interface connector



Pin Assignment Chart, 25-pin female interface connector, RS-232:

- Pin 1: Shield
- Pin 2: Data output (TxD)
- 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 connected
- Pin 11: +12 V
(operating voltage
for Sartorius printer)
- Pin 12: Reset _ Out *)
- Pin 13: +5 V output
- 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 output

*) = Hardware restart

* 더 자세한 정보를 원하시면,
"Data interface description fo
ED, GK and GW model "
www. Sartorius.com 에서의
download center 에서 다운받
실 수 있습니다.

* For the remote switch

문제 발생

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

Display	Cause	Solution
No segments appear on the display	No AC power is available	Check the AC power supply
	The power supply is not plugged in	Plug in the power supply

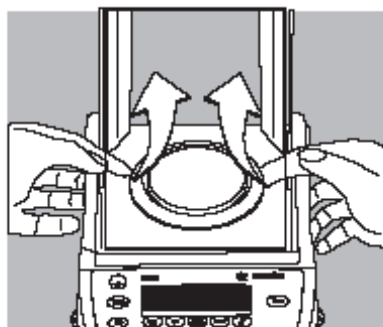
HIGH	The load exceeds the balance/ scale capacity	Unload the balance/scale
LOW or ERR 54	Something is touching the weighing pan	Move the object that is touching the weighing pan
APP.ERR.	Cannot store data: Load on weighing pan too light or no sample on pan while application is active	Increase load
WIS.ERR.	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.: - balance/scale not tared - load on weighing pan	Calibrate only when zero is displayed - Press Tare to tare the balance/scale - Unload the balance/scale
ERR 10	The Tare key is blocked when there is data in the second tare memory (net-total); only 1 tare function can be used at a time	Press CF to clear the tare memory and release the tare key
ERR 11	Tare memory not allowed	Press Tare
The weight readout changes constantly	Unstable ambient conditions (excessive vibration or draft) at the place of installation A foreign object is caught between weighing pan and balance/scale housing	Set up the balance/scale in another area Remove the foreign object
The weight readout is obviously wrong	The balance/scale was not calibrated/adjusted Balance/scale not tared before weighing	Calibrate/adjust the balance/scale Tare or zero the balance/scale before weighing

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

Contact information: Please point your Internet browser to: <http://www.sartorius.com>

* 위 문제로 응급조치 후 해결이 되지 않을 시 에는 (주)싸토리우스 코리아
로 연락하셔서 조치를 받으시기 바랍니다.T) (02) 575- 6945

유지, 관리



서비스

-문제 발생시 수리작업은 교육을 받은 싸토리
우스의 기술자가 직접 수행을 합니다.
만약, 임의로 수리 시 발생하는 문제에 대해서
는 책임을 질 수 가 없습니다.

관 리

- 전자저울의 전원을 뽑는다.



(프린터등이 연결시 전자저울의 전원을 끈 상태에서 분리한다.)
(내부에 액체가 들어가지 않도록 주의한다.)

- 알코올등을 이용하여 젖은 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.
- 스테레스 스틸 판 (사각의 밀판) 0.1mg 전자저울에서 밀판을 완전 분리 후 청소를 실시한다.
- 크리닝용 액체를 약간 묻혀서 구석구석 닦는다. (후 마른 천이용 마무리)

Overview

Specifications

Specifications

Built-in motorized calibration weight	All models with the designation suffix ED...-CW, GK..., GW.... or with a readability of 0.0001 g and all verified models		
AC power source/power requirements, voltage, frequency	AC adapter 230 V or 115 25 V, +15% to - 20%, 48-60 Hz		
Power consumption	VA	maximum 16; typical 8 (STNG6)	
Approx. hours of operation with the YRB05Z rechargeable battery pack (backlighting on)	h	35	

Ambient Conditions

The specifications given here are ensured under the following ambient conditions:

Operating temperature range	+10 to +30°C (273 to 303 K, 50 to 86°F)
Allowable ambient operating temperature	+5 to +40°C (41 to 104°F)
Proper functioning is ensured within an ambient operating temperature range of 5 to 40°C (41 to 104°F).	

Specifications for Individual Models

Model		ED224S	ED124S	GK1403
Weighing capacity		220 g	120 g	1400 ct
Readability		0.0001 g	0.0001 g	0.001 ct
Tare range (subtractive)		220 g	120 g	1400 ct
Repeatability (std. deviation)	≤±	0.0001 g	0.0001 g	0.001 ct
Linearity	≤±	0.0002 g	0.0002 g	0.002 ct
Response time (average)	s	2.5	2.5	1.5

response time (average)	s	2.5	2.5	1.5
Sensitivity drift within +10 to +30°C	≤±/K	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)	200 (E2)
Net weight, approx.:	kg	4.8	4.8	4.7
Weighing pan size	mm	90 Ø	90 Ø	90 Ø
Whg. chamber height	mm	230	230	160
Dimensions (WxDxH)	mm	230 × 303 × 330		230 × 303 × 260

Model		GK1203	GK703/ GK703-ST	GK303
Weighing capacity		1200 ct	700 ct	300 ct
Readability		0.001 ct	0.001 ct	0.001 ct
Tare range (subtractive)		1200 ct	700 ct	300 ct
Repeatability (std. deviation)	≤±	0.001 ct	0.001 ct	0.001 ct
Linearity	≤±	0.002 ct	0.002 ct	0.002 ct
Response time (average)	s	1.5	1.5	1.5
Sensitivity drift within +10 to +30°C	≤±/K	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 (F2)	50 (F2)
Net weight, approx.:	kg	4.7	4.7	4.7
Weighing pan size	mm	90 Ø	90 Ø/35 Ø	90 Ø
Whg. chamber height	mm	160	160/38	160
Dimensions (WxDxH)	mm	230 × 303 × 260	Model GK703-ST: 230 × 303 × 138	

Specifications for Individual Models

Model		ED623S ED623S-CW	ED423S ED423S-CW ED423S-DS	ED323S ED323S-CW ED323S-DS
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 (std. 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 to +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: 0.05–0.4 (depends on filter level selected)		
External calibration weight (of at least accuracy class...)	g	500 (E2)	200 (E2)	200 (F1)
Net weight, approx:	kg	3.2 3.6	3.2 3.6 4.4	3.2 3.6 4.4
Weighing pan size	mm	115 Ø	115 Ø	115 Ø
Dimensions (WxDxH)	mm	230 × 303 × 136	230 × 303 × 136 ED...-DS: 230 × 303 × 330	

Model		ED153 ED153-CW ED153-DS	GK3102	GK2202
Weighing capacity		150 g	3 100 ct	2200 ct
Readability		0.001 g	0.005 ct	0.005 ct
Tare range (subtractive)		150 g	3 100 ct	2200 ct
Repeatability (std. deviation)	≤±	0.001 g	0.005 ct	0.005 ct
Linearity	≤±	0.002 g	0.01 ct	0.01 ct
Response time (average)	s	1.3	1	1
Sensitivity drift within +10 to +30°C	≤±/K	3.3 · 10 ⁻⁶	2 · 10 ⁻⁶	2 · 10 ⁻⁶
Adaptation to ambient conditions		By selection of 1 of 4 optimized filter levels; display update: 0.05–0.4 (depends on filter level selected)		
External calibration weight (of at least accuracy class...)	g	100 (F1)	500 (E2)	200 (E2)
Net weight, approx:	kg	2.6 3.0 3.8	4.4	4.4
Weighing pan size	mm	115 Ø	115 Ø	115 Ø
Dimensions (WxDxH)	mm	230 × 303 × 136 ED...-DS: 230 × 303 × 330	230 × 303 × 260	

Model		ED6202S GW6202 ED6202S-CW	ED4202S ED4202S -CW	ED3202S GW3202 ED3202S-CW	ED2202S ED2202S -CW	ED822 ED822 -CW
Weighing capacity	g	6200	4200	3200	2200	820
Readability	g	0.01	0.01	0.01	0.01	0.01
Tare range (subtractive)	g	6200	4200	3200	2200	820
Repeatability (std. deviation)	≤±g	0.01	0.01	0.01	0.01	0.01
Linearity	≤±g	0.02	0.02	0.02	0.02	0.02
Stabilization time (typical)	s	1.1	1.1	1.1	1.1	1.0

Stabilization time (typical)	2	1.1	1.1	1.1	1.1	1.1
Sensitivity drift within +10 to +30°C	≤±/K	2 · 10 ⁻⁶	2 · 10 ⁻⁶	2 · 10 ⁻⁶	2 · 10 ⁻⁶	5 · 10 ⁻⁶
Adaptation to ambient conditions	By selection of 1 of 4 optimized filter levels; display update: 0.05–0.4 (depends on filter level selected)					
External calibration weight (of at least accuracy class...)	g	5000 (E2)	2000 (E2)	2000 (F1)	2000 (F1)	500 (F2)
Net weight, approx:	kg	3.1 3.1 3.5	3.1 3.1 3.5	3.1 3.1 3.5	3.1 3.1 3.5	2 2 2.6
Weighing pan size	mm	180 × 180	180 × 180	180 × 180	180 × 180	150 Ø
Dimensions (WxDxH)	mm	230 × 303 × 91				230×303×87

Model		ED8201 ED8201-CW	GW7201	ED5201 ED5201-CW	ED2201 ED2201-CW
Weighing capacity	g	8200	7200	5200	2200
Readability	g	0.1	0.1	0.1	0.1
Tare range (subtractive)	g	8200	7200	5200	2200
Repeatability (std. deviation) ≤±g	0.1	0.1	0.1	0.1	0.1
Linearity ≤±g	0.1	0.1	0.1	0.1	0.1
Response time (average)	s	1	1	1	1
Sensitivity drift within +10 to +30°C	≤±/K	10 · 10 ⁻⁶			
Adaptation to ambient conditions		By selection of 1 of 4 optimized filter levels; display update: 0.05–0.4 (depends on filter level selected)			
External calibration weight (of at least accuracy class...)	g	5000 (F2)	5000 (F2)	5000 (F2)	2000 (F2)
Net weight, approx.:	kg	2.7 3.5	2.7	2.7 3.5	2.7 3.5
Weighing pan size	mm	180 × 180			
Dimensions (WxDxH)	mm	230 × 303 × 91			

Accessories

External calibration weights:

For model	Accuracy class	Weight in grams	Order no.:
ED423S/...-DS	E2	200	YCW5228-00
ED623S	E2	500	YCW5528-00
ED4202S	E2	2000	YCW6228-00
ED6202S, GW6202	E2	5000	YCW6528-00
ED153S/...-DS	F1	100	YCW5138-00
ED323S/...-DS	F1	200	YCW5238-00

ED3202S, ED2202S	F1	2000	YCW6238-00
ED822	F2	500	YCW5548-00
ED2201	F2	2000	YCW6248-00
ED8201, ED5201	F2	5000	YCW6548-00
or alternatively	± 25 mg	5000	YSS653-00

Product
Data printer
 with date, time, statistics
 evaluation, transaction
 counter functions and LCD

Order No.
YDP03-OCE

Remote display ¹⁾, reflective
 (for connection to
 data interface port)

YRD02Z

External rechargeable battery pack
YRB05Z

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/scale
 and plug it into the battery pack

SartoConnect ¹⁾,
 data transfer software for direct
 transmission of weight values
 to another program (e.g., MS Excel)
 – with RS-232C

connecting cable, length:

1 m (~20 in)

YSC01L

– with RS-232C

connecting cable, length:

5 m (~16 ft)

YSC01L5

– with RS-232C

connecting cable, length:

15 m (~50 ft)

YSC01L15

Product
Density determination kit¹⁾
 – for ED224S, ED124S

YDK01LP

Standard Operating Procedure

optimum use of your balance/scale
 in quality-management systems

YSL01E

Industrial AC adapter, model ING2,
 protection rating: IP65 in accordance
 with EN 60529

– for 230 V

69 71899

– for 120 V

69 71500

Data cable

– for connecting a computer
 with a USB port

YCC01-USBM2

– for computer connection,
 25-pin

73 57 312

– for computer connection, 9-pin

73 57 314

Adapter cable

6965619




from D-Sub 25-pin male
 connector to D-Sub 9-contact
 female connector; length: 0.25 m

¹⁾ Not for verified models

Universal remote control switch

Order No.

for remote control of the following functions:

, ,  or a function key

(see "Configuration" for details):

Foot switch with T-connector

YFS01

Hand switch with T-connector

YHS02

T-connector

YTC01

Note:

The T-connector is not intended for connecting multiple

The I connector is not intended for connecting multiple intelligent peripheral devices, such as PCs or YDP03-OCE printers.

Ionizing blower for eliminating static electricity	
- 220 V	YIB01-ODR
- 110 V	YIB01-OUR
Stat-Pen anti-static device for eliminating electrostatic charges on samples and containers (100 V to 230 V, 50/60 Hz)	
YSTP01	
Anti-vibration balance/scale table	
- for precise, reliable weighing operations	YWT01
- made of cast stone with shock absorbers	YWT03
Bracket for wall mounting	
YWT04	
Gem trays/Weighing bowls	
- 300 ml, weight: 86 g, stainless steel	6407
- 1000 ml, wt.: 240 g, stainless steel	641211
- 500 ml	641212
- 300 ml, wt.: 22 g, aluminum	69641304
- 110 ml, 90 mm Ø, aluminum	69GP0003
- 270 ml, wt.: 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 Ø, stainless steel	YWP04G



**Declaration of Conformity
to Council Directives 89/336/EEC and 73/23/EEC
(amended by Directive 93/68/EEC)**

The electronic precision weighing instrument of the series
ED/GK/GW/XX.....-.....

meets the applicable requirements of the test standards listed below, in conjunction
with the associated power supplies, auxiliary peripheral devices and installation

equipment listed in Annex A2 (see Annex A1 for a technical description and a list of the individual versions).

1. Electromagnetic Compatibility

1.1 Source for 89/336/EEC: Official Journal of the European Communities, No. 2004/C98/05

EN 61326 Electrical equipment for measurement, control and laboratory use
EMC requirements

Limitation of emissions: Residential areas, Class B

Defined immunity to interference: Industrial areas, continuous unmonitored operation

2. Safety of Electrical Equipment

2.1 Source for 73/23/EEC: Official Journal of the European Communities, No. 2004/C103/02

EN 61010 Safety requirements for electrical equipment for
measurement, control and laboratory use

Part 1: General requirements

EN 60950-1 Information technology equipment

Safety

Part 1: General requirements

Sartorius AG
37070 Goettingen, Germany
2005



W. Obermann
Senior Vice President, R&D
Electronic Engineering
Mechatronics Division



Dr. D. Klausgrete
Head of
International Certification Management
Mechatronics Division



Declaration of Type Conformity to Directive No. 90/384/EEC

This declaration is valid for non-automatic electromechanical weighing instruments for use in legal metrology. These weighing instruments accepted for legal metrological verification have an EC Type-approval Certificate. The model(s) concerned is/are listed below along with the associated type, accuracy class, and EC type-approval certificate number:

Model	Type	Accuracy Class	EC Type Approval No.
ED...-OCE	BD ED 100	Ⓛ	D06-09-006
ED...-PCE	BD ED 100	Ⓛ	D06-09-006

GK...-OCE	BD ED 100	①	D06-09-006
ED...-OCE	BD ED 200	②	D06-09-006
ED...-PCE	BD ED 200	③	D06-09-006
GW...-OCE	BD ED 200	④	D06-09-006

SARTORIUS AG declares that its weighing instrument types comply with the requirements of the Council Directive on non-automatic weighing instruments, no. 90/384/EEC of 20 June 1990, the associated European Standard "Metrological aspects of non-automatic weighing instruments," No. EN 45501; the amended, currently valid versions of the national laws and decrees concerning legal metrology and verification in the Member States of the European Union (EU) and the Signatories of the Agreement on the European Economic Area, which have adopted this Council Directive into their national laws; and with the requirements stipulated on the Type-Approval Certificate for verification. This Declaration of Type Conformity is valid only if the ID label on the weighing instrument has the CE mark of conformity and the green metrology sticker with the stamped

Sartorius AG
37070 Göttingen, Germany
Signed in Göttingen, 08.05.2006


Dr. G. Maaz
President of the Mechatronics Division

letter "M" (the two-digit number in large print stands for the year in which the mark has been affixed):



If these marks are not on the ID label, this Declaration of Type Conformity is not valid. Validity can be obtained, for example, by submitting the weighing instrument for final action to be taken by an authorized representative of SARTORIUS AG. The validity of this Declaration of Type Conformity shall expire upon any tampering with, repair or modification of this weighing instrument or, in some Member States, on the date of expiration. This declaration applies only to the weighing instrument without peripheral devices.

The operator of this weighing instrument shall be responsible for obtaining an authorized renewal of the verification, such as subsequent or periodic verification, of the weighing instrument for use as a legal measuring instrument.


J. Reinhold
Head of the Production Department
Mechatronics / Weighing Technology Division

L0P-3.225
PT06K01.doc

Physikalisch-Technische Bundesanstalt
Braunschweig und Berlin

PTB



EG-Bauartzulassung
EC type-approval certificate

Zulassungsinhaber: Sartorius AG

Issued to:

Weender Landstr. 94-108
37075 Göttingen

Rechtsbezug:
in accordance with:

§ 13 des Gesetzes über das Mess- und Eichwesen (verification act)
vom/dated 23. März 1992 (BGBl. I S. 711), zuletzt geändert am (last
amended on) 25.11.2003 (BGBl. I S. 2304), in Verbindung mit Richtlinie
(in connection with council directive) 90/384/EWG, geändert durch (amended
by) 93/68/EWG

Bauart:
in respect of:

Nichtselbsttätige elektromechanische Waage mit oder ohne Hebelwerk
Nonautomatic electromechanical weighing instrument with or without
lever system

Typ / Type:

BD ED 100, BD ED 200

- (I) Max 50...240 g, $e = 1...2$ mg, $n \leq 240000$
(II) Max 1...8200 g, $e = 0,01...1$ g, $n \leq 62000$

Zulassungsnummer:
Approval number:

D06-09-006 1. Revision

Gültig bis:
Valid until:

06.02.2016

Anzahl der Seiten:
Number of pages:

8

Geschäftszeichen:
Reference No.:

PTB-1.12-4023582

Benannte Stelle:
Notified Body:

0102

Im Auftrag
By order


Marcus Link



Braunschweig, 03.05.2006

Siegel
Seal

PTB-0023

Die Hauptmerkmale, Zulassungsbedingungen und Auflagen sind in der Anlage enthalten, die Bestandteil der Revision der EG-Bauartzulassung ist. Hinweise und eine Rechtsbehelfsbelehrung befinden sich auf der ersten Seite der Anlage.
The principal characteristics, approval conditions and special conditions, if any, are set out in the Annex which forms an integral part of this Revision of the EC type-approval certificate. For notes and information on legal remedies, see first page of the Annex.