

The New Practum.<sup>®</sup>  
Start Weighing Right.



The New Practum.<sup>®</sup>  
Start Weighing Right.



Get best value for your money, without compromises in precision and reliability. Rely on consistent readings and excellent repeatability ensured by Sartorius quality and technology developed and designed in Germany.

**SPECIAL OFFER**

A calibrated balance is fundamental for getting reliable weighing results – therefore every Practum® comes with a high quality external calibration weight, including a calibration certificate, to support correct results.



Practum

# What I Like...

## **Practical and Sturdy Draft Shield**

The sturdy draft shield will stand up to the daily use even by inexperienced operators. Moreover, it can be completely disassembled to conveniently clean the glass parts in a dishwasher.

## **Rugged Weigh Cell for Reliable Results – Every Time, All the Time**

High-resolution, overload protected weigh cells, made in Germany, ensure durable and reliable performance of the Practum®.

## **Hassle-Free Below-Balance Weighing of Oversize Containers**

Extra-large volumes or magnetic substances are easy to weigh accurately using the easily accessible weighing port.





**Supervisor Lock**

This practical feature enables teachers and professors to lock the menu. As a result, students will concentrate on learning.

**USB Port and Data Transfer to Microsoft® Excel for Perfect Records**

The USB port automatically detects when a Sartorius laboratory printer is connected to the Practum®. This interface also enables you to transfer your weighing data directly into a Microsoft® Excel spreadsheet, without the need for additional software.

**Changes Readily Documented by Interval-Controlled Printout of Weights**

The Practum® will document weights automatically and reliably, such as slow filling processes.

**Extra-large Feet and Up-front Level Indicator for Easy Leveling**

Two large, easy-to-adjust feet and an outstandingly simple-to-read level indicator on the front make it a breeze for you to level the Practum®. This helps ensure accurate, repeatable weighing results.



**Plug & Play**  
Standard printer YDP40

# Practum

# The New Practum<sup>®</sup>. Start Weighing Right.

You can't go wrong with the new Practum<sup>®</sup>:  
a foolproof, exceptionally rugged and reliable lab balance.  
Practum<sup>®</sup> has everything you need for standard applications.  
So you pay only for what you really need.



## **Weighing**

You name it, Practum<sup>®</sup> will weigh it – from the smallest samples to the heaviest objects in the analytical range, and from entire chunks of solids to the finest quantities of pastes, powders and liquids, Practum<sup>®</sup> provides incomparable weighing reliability.



## **Counting**

Use Practum<sup>®</sup> to count identical materials, such as tablets, screws and similar parts. Using Practum<sup>®</sup> is easier, faster and more accurate than manual counting.



## **Density Determination**

With Practum<sup>®</sup> and our Optional Density kit, you can easily and accurately determine the density of a solid object.



## **Peak Hold Function**

Forget now you see it, now you don't. Effortlessly read off the maximum force released in milliseconds when a switch is activated, or weigh a bulky sample that covers the display. Practum<sup>®</sup> will conveniently determine and "hold" the peak value for you.



## **Checkweighing**

Practum<sup>®</sup> is ideal when you need to check whether different objects, such as packaged units, are within a specific tolerance range.



## **Animal Weighing**

Does your environment "rock" in the negative sense? Is it exposed to vibration or laminar flow? Or do you need to know the exact weight of a hyper little mouse that just won't sit still? Relax with "rocksteady" readings on Practum<sup>®</sup>.



## **Percentage Weighing**

If you need to compare different samples and determine how much they vary from a reference standard, this function will do the math for you.



# Practum



## Technical Data Practum®

Model		224	124	64	513	313	213	3102	2102	1102	612	412	6101	5101	2101	6100	5100
Weighing capacity	g	220	120	60	510	310	210	3100	2100	1100	610	410	6100	5100	2100	6100	5100
Readability	mg	0.1	0.1	0.1	1	1	1	10	10	10	10	10	100	100	100	1000	1000
Repeatability	mg	0.1	0.1	0.1	1	1	1	10	10	10	10	10	100	100	100	500	500
Linearity	mg	0.2	0.2	0.2	2	2	2	30	30	30	30	30	300	300	300	1000	1000
Typical stabilization time	s	2	2	2	1	1	1	1.5	1.5	1.5	1.5	1.5	1.5	1.5	1.5	1	1
External calibration weight	g	200 E2	100 E2	50 E2	500 F1	200 E2	200 E2	2000 F1	2000 F1	1000 F1	500 F2	200 F2	5000 F2	5000 F2	2000 F2	5000 F2	5000 F2
Weighing pan size	Ø mm	90	90	90	120	120	120	180	180	180	180	180	180	180	180	180	180
Weighing chamber height	mm	209	209	209	209	209	209										
Net weight, approx.	kg	4.5	4.5	4.5	4.9	4.9	4.9	3.1	3.1	3.1	3.1	3.1	3.1	3.1	3.1	3.1	3.1
	lb	9.9	9.9	9.9	10.8	10.8	10.8	6.8	6.8	6.8	6.8	6.8	6.8	6.8	6.8	6.8	6.8
Dimensions	mm	360×216×320						360×216×95									
D×W×H	inch	14.1×8.5×12.6						14.1×8.5×3.75									
Data interface	mini USB – Automatic recognition of Sartorius printer models YDP40 – Direct data transfer to Microsoft® application programs without any additional software – Programmable interval for data output – Choice of SBI, table format or text format data transfer protocols																
Supervisor lock	Menu lock to prevent unintended manipulation																
Housing and draft shield	Robust and easy to clean housing with easy to operate draft shield (where applicable)																
Standard application programs	Weighing, Density, Percentage, Checkweighing, Peakhold, Counting and Animal weighing																
Theft lock	Kensington lock and lockdown capability for cable or chain																
In-use cover	Standard for all models																
Underfloor weighing	Integrated																
Accessories	See Practum® accessory list																
Technology and construction	Made in Germany																



◆ [www.sartorius.com](http://www.sartorius.com)

### Germany

Sartorius Lab Instruments GmbH & Co. KG  
 Weender Landstrasse 94–108  
 37075 Goettingen

Phone +49.551.308.0  
 Fax +49.551.308.3289