



Practum[®]

Laboratory Balances

Simplifying Progress

SARTORIUS

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.



What I Like



Plug & Play
Standard printer YDP40

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.

Supervisor Lock

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

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®.

Changes Readily Documented by Interval-Controlled Printout of Weights

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



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.

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.

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.

Start Weighing Right

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



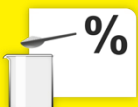
Weighing

You name it, Practum® 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® provides incomparable weighing reliability.



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® will conveniently determine and "hold" the peak value for you.



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.



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®.



Checkweighing

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



Counting

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



Density Determination

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



Practum® – Technical Specifications


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
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
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										
Dimensions D×W×H	mm	360×216×320						360×216×95									
	inch	14.1×8.5×12.6						14.1×8.5×3.75									
Data interface	Mini USB - Automatic recognition of Sartorius printer models YDP40 and YDP20-OCE - 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																
Technology and construction	Made in Germany																

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com