



Quintix[®] Pro

Redefining the Standard

Simplifying Progress

SARTORIUS

Quintix® Pro

Designed to be a True
Universal Balance

Quintix® Pro Laboratory Balances deliver the performance you need, with added premium features that raise the bar in usability, flexibility, and eco-conscience design.



Useability

Enhanced user experience with new premium features including:

- Easy access to menu option with a large, 7-inch, high-resolution full-touch graphic display with modern upgraded user interface
- Ensure the highest weighing accuracy with the internal motorized levelling (optional) and automatic internal adjustment isoCAL
- Quick access to the integrated digital manual for common help topic
- Fast adaptation of balance performance to the unique needs of each application through the improved semi-automatic filter adaption (AFA)
- Seamless connectivity features include interface options, easy data transfer, and GLP | GMP compliant printout

Flexibility

Flexibility for use in a wide range of environments and applications:

- Customization with different tiers of user management access with passcode protection
- Tested for use in cleanrooms grade A and B, and special environments e.g., vacuum, argon, and nitrogen.
- Smart design and chemically-resistant surfaces made of PBT, glass and stainless steel support routine cleaning protocols
- 15 integrated application programs for common weighing tasks
- Extended language package with 15 language options

Sustainability

Reduced environmental impact from the manufacturing plant to the bench:

- Climate-conscience production process utilizes recycled materials e.g., aluminum, in addition to environment-friendly packaging
- Power-saving operation with stand-by or switch-off modes
- Automatic dim function, reduces display brightness when balance is not in use to save power
- User-defined, automatic start timer powers up the instrument in time for daily use
- Digital user manual minimizes paper consumption, while allowing quick access to information



Modern User Interface

Effortless Operation

Using the Quintix® Pro is effortless thanks to a modern usability concept. The startup assistant allows quick and easy setup of the balance, while the built-in interactive manual is ready to assist, whenever you need help. With 15 available language options, the Quintix® Pro is the perfect fit for any lab.

Status Center

The Status Center on Quintix® Pro provides a comprehensive summary of the balance's most important status information with just a single click.



User Management

The advanced user management system in Quintix® Pro enables you to assign individual roles to each user and customize them according to specific user needs. This feature allows Quintix® Pro to limit menu access strictly to authorized personnel, thereby providing comprehensive control and protection over your weighing procedures.



Correct Levelling

Level Status at a Glance

With Quintix® Pro it is easy to verify if your balance is leveled correctly. A quick glance at the conveniently positioned level indicator is all you need.

Continuous Real-Time Level Monitoring

The LevelControl function of Quintix® Pro guarantees a perfectly leveled balance. It continuously monitors the level status in real time, instantly detecting any deviations. Should any tilt be detected, it responds immediately to safeguard your data: it will alert you and temporarily block data transfer to a printer or computer. Quintix® Pro will then guide you through the leveling procedure in real time.

With the Quintix® Pro LevelControl, you can rest assured that your measurements are completely accurate.



Motorized Levelling

Leveling is even easier with the motorized feet. The Quintix® Pro levels itself fully automatically at the push of a button.



Monitoring Functions and Security

isoCAL: Reliability Guaranteed

Temperature fluctuations can compromise the precision of your weighing results. Quintix® Pro eliminates this risk from the outset with its temperature and time-controlled calibration and adjustment function, isoCAL. By customizing the time interval for isoCAL, you can ensure compliance with your Standard Operating Procedures (SOPs).



Secure Weighing Data Continuously

Take advantage of the various safety level settings available on Quintix® Pro for secure data transfer. You can tailor Quintix® Pro's automatic monitoring systems to meet your specific security requirements. Whether you choose a high-, standard-, or low-level monitoring, Quintix® Pro ensures that only accurate data is processed.



Enhanced Traceability

The Quintix® Pro's internal Device-Log function records all calibration and leveling procedures, as well as any changes to the balance and weighing settings. This creates a traceable log file with date and time stamps, enhancing your quality assurance processes.



Improved Auto Adaption to Environmental Conditions (AFA)

Whether your environment is stable or unstable, Quintix® Pro can be easily adapted to specific environmental conditions with just a single click.



Protection and Cleanability

Tested for Use in Special Environments e.g., Vacuum, Argon, and Nitrogen



Extended Protection with In-Use-Cover

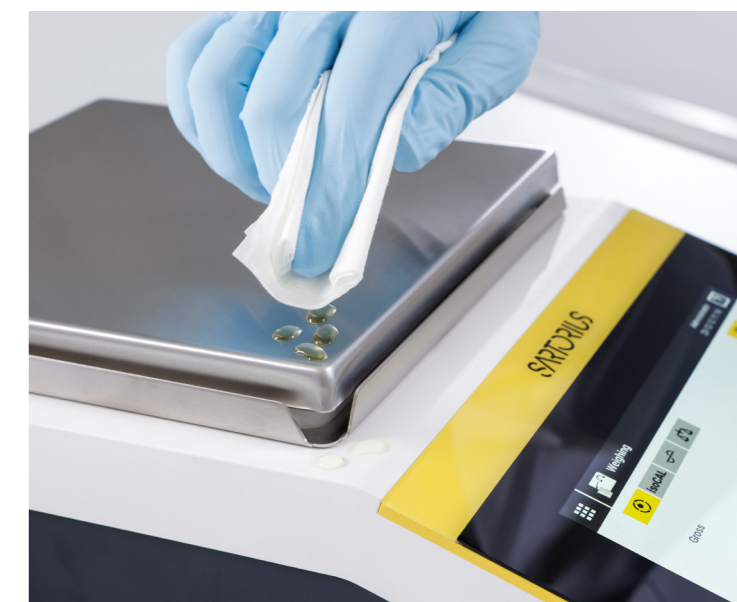
Enhance the protection of Quintix® Pro by utilizing an in-use cover. This cover envelops the entire housing of the balance, offering additional defense against contamination.

The additional protection covers for the weighing pan and draft shield base plate is particularly useful in areas where hazardous materials are handled.



Easy Cleaning

The contemporary design of Quintix® Pro facilitates easy and straightforward cleaning of the balance. The use of durable materials such as polybutylene terephthalate (PBT), stainless steel, and glass not only ensures high chemical resistance, but also prevents cross-contamination.



Documentation and Data Transfer: Easy, Fast and Reliable

Be future ready! Quintix® Pro supports multiple connection to external devices via three USB type C interfaces, RS232 9-pin interface, and an ethernet interface. Transfer weighing data or calibration reports to a USB stick or directly to a database via the ethernet interface.

Data Transfer: Direct and Correct

Effortlessly connect your Quintix® Pro to your computer or lab printer using the balance's USB Type-C interface port and Plug & Play technology. Its PC-Direct feature allows for the seamless transfer of your measurement results directly to your preferred Microsoft® Office program. This eliminates the need for additional software and prepares your data in your chosen format, ready for calculations in spreadsheets.

Data Documentation: Fast and Easy

When directly connected to a Sartorius laboratory printer, Quintix® Pro generates fast and accurate printouts of your weighing results. You can use standard printer paper or practical, continuous-length paper label rolls, or even fade-resistant paper for reliable records with unique identification of your samples.

Sample Identification: Even More Reliable and Precise with Quintix® Pro

If you're working with samples that require unique identification in regulated environments, you'll appreciate the built-in sample identification feature of Quintix® Pro. This function allows you to tailor the printout information to your specific needs. With an interactive touchscreen display and alphanumeric input capability, Quintix® Pro simplifies the process of entering complex sample IDs, batch IDs, or project IDs. Generate reliable and accurate sample records with ease using Quintix® Pro.



Universal Applications

Integrated Application Programs for Efficient Working Routines

The combination of user-friendly design and diverse integrated application programs make Quintix® Pro a truly universal balance that is suitable for any environment:

Weighing | Dosing, Mass unit conversion, Counting, Percentage weighing, Calculation, Mixing, Components, Density determination, Statistics, Check weighing, Animal weighing, Peak hold, Pipette smart check, Differential weighing, Underfloor weighing.



Service Tools

Flexible Maintenance and Reliable Backup

Define service and maintenance interval for each Quintix® Pro individually to maintain compliance with your existing SOPs. You can easily export a backup of the balance settings to external storage devices, like a USB stick. Saved records are always accessible and can also be transferred to other devices.

Each Quintix® Pro comes with an individual factory acceptance test report, confirming compliance with all specifications.



Life Cycle Services for Your Quintix® Pro

Weighing is a critical process in all laboratories, where the reliability and reproducibility of results as well as precision and accuracy are of the utmost importance. In order to consistently achieve trustful weighing results throughout the entire equipment life cycle, at Sartorius high quality products go hand in hand with professional life cycle services.

We are proud to offer an extensive range of certified and accredited services for your laboratory balances, aimed at ensuring the utmost quality of your weighing results, prolonging the equipment longevity, and minimizing downtimes. Our services ensure that your balances operate flawlessly within their tolerance limits and obtain constantly compliant, reproducible and reliable measurement data.



Our Services at a Glance



Installation and commissioning:
Reliable results right from the start

Your advantage:

- Outstanding weighing performance, steadfast results from the start
- Seamless integration of equipment into your laboratory environment
- Optimum system performance and results
- Expertly trained operating personnel



Equipment qualification (IQ | OQ):
Compliance with GMP requirements

Your advantage:

- Full compliance with regulatory requirements (GMP | GLP)
- Audit security
- All critical equipment aspects regimented due to our risk-based approach
- Service performed with all required regulatory and technical knowledge



Preventative maintenance and Service Contracts:
Long-term process safety and reliable results

Your advantage:

- Optimal operation of your system
- Extended instrument lifetime
- Compliance with regulatory requirements
- Full cost control, with service contracts tailored to your specific needs



Calibration Services:
Traceability of weighing results, regulatory compliance

Your advantage:

- Accredited calibration certificates with ISO/ IEC 17025 in most countries
- Guaranteed traceability to national standards
- Minimum weight certificates according USP <41> and Ph. Eur. 2.1.7
- Certificates for type approved balances according to OIML R76

Technical Specifications

Analytical Balances

| Model | Levelling feature | Weighing capacity [g] | Readability [mg] | Repeat-ability [mg] | Repeatability, typical [mg] | Stabilization time, typical [s] | Weighing pan size [mm] | Weighing chamber height* [mm] | Dimension (W × D × H) [mm] |
|------------------------------|---------------------------------|-----------------------|------------------|---------------------|-----------------------------|---------------------------------|------------------------|-------------------------------|----------------------------|
| QTX324IRU-1x QTX324IMU-1x | Level support Motorized feet | 320 | 0.1 | ± 0.1 | ± 0.08 | ≤1.5 | Ø 90 | 240 | 220 × 377 × 346 |
| QTX224IRU-1x QTX224IMU-1x | Level support Motorized feet | 220 | 0.1 | ± 0.1 | ± 0.08 | ≤1.5 | Ø 90 | 240 | 220 × 377 × 346 |
| QTX124IRU-1x QTX124IMU-1x | Level support Motorized feet | 120 | 0.1 | ± 0.1 | ± 0.08 | ≤1.5 | Ø 90 | 240 | 220 × 377 × 346 |

Milligram Balances

| Model | Levelling feature | Weighing capacity [g] | Readability [mg] | Repeat-ability [mg] | Repeatability, typical [mg] | Stabilization time, typical [s] | Weighing pan size [mm] | Weighing chamber height* [mm] | Dimension (W × D × H) [mm] |
|--------------------------------|---------------------------------|-----------------------|------------------|---------------------|-----------------------------|---------------------------------|------------------------|-------------------------------|----------------------------|
| QTX1503IRU-1x QTX1503IMU-1x | Level support Motorized feet | 1500 | 1 | ± 1 | ± 0.5 | ≤1 | Ø 120 | 240 | 220 × 377 × 346 |
| QTX1203IRU-1x QTX1203IMU-1x | Level support Motorized feet | 1200 | 1 | ± 1 | ± 0.5 | ≤1 | Ø 120 | 240 | 220 × 377 × 346 |
| QTX623IRU-1x QTX623IMU-1x | Level support Motorized feet | 620 | 1 | ± 1 | ± 0.5 | ≤1 | Ø 120 | 240 | 220 × 377 × 346 |
| QTX423IRU-1x QTX423IMU-1x | Level support Motorized feet | 420 | 1 | ± 1 | ± 0.5 | ≤1 | Ø 120 | 240 | 220 × 377 × 346 |
| QTX323IRU-1x QTX323IMU-1x | Level support Motorized feet | 320 | 1 | ± 1 | ± 0.5 | ≤1 | Ø 120 | 240 | 220 × 377 × 346 |

* Upper edge of the weighing pan to the lower edge of the upper draft shield panel
x = Term for country-specific models

Precision Balances

| Model | Levelling feature | Weighing capacity [g] | Readability [mg] | Repeat-ability [mg] | Repeatability, typical [mg] | Stabilization time, typical [s] | Weighing pan size [mm] | Dimension (W × D × H) [mm] |
|----------------------------------|---------------------------------|-----------------------|------------------|---------------------|-----------------------------|---------------------------------|------------------------|----------------------------|
| QTX6202IRO-1x QTX6202IMO-1x | Level support Motorized feet | 6200 | 10 | ± 10 | ± 5 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX4202IRO-1x QTX4202IMO-1x | Level support Motorized feet | 4200 | 10 | ± 10 | ± 5 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX3202IRO-1x QTX3202IMO-1x | Level support Motorized feet | 3200 | 10 | ± 10 | ± 5 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX2202IRO-1x QTX2202IMO-1x | Level support Motorized feet | 2200 | 10 | ± 10 | ± 5 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX1202IRO-1x QTX1202IMO-1x | Level support Motorized feet | 1200 | 10 | ± 10 | ± 5 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX12201IRO-1x QTX12201IMO-1x | Level support Motorized feet | 12200 | 100 | ± 100 | ± 50 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX10201IRO-1x QTX10201IMO-1x | Level support Motorized feet | 10200 | 100 | ± 100 | ± 50 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX8201IRO-1x QTX8201IMO-1x | Level support Motorized feet | 8200 | 100 | ± 100 | ± 50 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX6201IRO-1x QTX6201IMO-1x | Level support Motorized feet | 6200 | 100 | ± 50 | ± 50 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |
| QTX3201IRO-1x QTX3201IMO-1x | Level support Motorized feet | 3200 | 100 | ± 50 | ± 50 | ≤0.9 | 182 × 182 | 215 × 377 × 95 |

* Upper edge of the weighing pan to the lower edge of the upper draft shield panel
x = Term for country-specific models

Contact Details:

Contact: