

The New Gold and Carat Balance Series II

The Highest Precision "Made in Germany"

Simplifying Progress



New GL Gold and GCL Carat Balances

- Quality and Technology "Made in Germany"
- Highly accurate results, guaranteed via monolithic weigh cell
- Draft shield with a weighing chamber height of 160 mm
- isoCAL: Automatic, built-in calibration and adjustment function
- LED and touch technology display
- "PC-direct feature" for easily & directly data transfer
- Automatically detects Sartorius accessories (e.g. printer, second display)



High Capacity Balances: Cubis® II Series

- Quality and Technology "Made in Germany"
- Highly accurate results, guaranteed via monolithic weigh cell
- Large TFT touch display for easy operation
- Support samples with large and high weight
- Suitable for gold shops for weighing to check gold stock in and out



Printer YDP20-0CE

- Dot matrix printer (RS232)
- A touch-control keypad for easy operation
- Design to receive weighing data from a connected balance and print it on plain paper
- Equip with the real-time clock to ensure the printed actual date and time



Second display YSD01

- Directly connect to Sartorius balances
- Display data or status data simultaneously to the balance
- Display the weight unit symbol in "g" (gram), "ct" (carat), etc.