

Cubis® II Premium Balance

Data Integrity and Connectivity

Need to comply with 21CFR11? Need the fastest stability on the market? Cubis® II – Your Balance, Your Way.

Features

- The world's highest resolution in a micro weighing instrument with up to 60 million divisions
- Integrated Software - no need for additional computer
- Full end-to-end data integrity. 21 CFR Part 11 compliance, integrated audit-trail, state-of-the-art user management
- Error-free operation. Individual QApp workflows, motorized auto-leveling*
- Integrated climate sensors
- Individual sample holders
- Outstanding service support. Integrated status center, integrated service functions preventive maintenance on accredited standards

* For all models up to a maximum capacity of 8.2 kg



Discover more about
Cubis® II Premium Balance



Please scan!