

Entris® II Balances

Best Value in Its Class for Basic Weighing Tasks

Features Entris® Basic Balance

- High chemical resistance – ensured using parts made from hard-wearing Polybutylene terephthalate (PBT), stainless steel and glass
- isoCAL – automatic internal calibration and adjustment
- Fastest stabilization time in its class
- Real „PC-direct“ feature
- 12 built-in applications with GLP | GMP compliant printout | data output

Features Entris® Advanced Balance

- Simplifies balance leveling with built-in real-time level support
- Increases reliability of weighing results using integrated protection systems
- Controlled access to balance settings with user management
- Gapless documentation thanks to CalAuditTrail
- Even easier usability with graphic touch display
- Provides gapless documentation with automatic documentation of calibration and levelling processes, and date and time stamp
- One additional built-in application: Pipette Smart Test



Discover more about
Entris® II Balances



Please scan!