

Laboratory Balances for Every Application, in Any Industry.

Regulated Application

Pharmaceutical, Biotech, Academic, R&C...

Cubis® II Balance

- Configured Balances have custom modularity with thousands of hardware/software applications
- 21 CFR Part 11 and EU Annex 11 models available
- Made to stock balances available on demand



Minimal Regulated Application

Chemical, Biopharma, Environmental...

Quintix® Pro Balance

- Highest weighing accuracy with the internal motorized levelling (optional) and automatic internal adjustment isoCAL
- Energy efficiency, reduced carbon footprint, and sustainability, minimizing environmental impact from the manufacturing plant to the bench.
- Seamless connectivity features include interface options, easy data transfer, and GLP | GMP compliant printout



Basic Weigh Application

Academic, Biotech, Industrial...

Entris® II Basic Balance

- Chemical resistance
- Fastest stabilization time in its class
- PC-direct, multiple data ports
- 12 built-in applications with GLP |
- GMP compliant printout | data output

Entris® II Advance Balance

- Real-time level support
- Large graphic touch display
- 13 built-in applications with
- GLP | GMP compliant printout | data output

