## S/RICH SILS

## Entris® II Balance Line

## Best Value in its Class for Basic Weighing Tasks

- Highest quality standards engineered in Germany
- isoCAL internal calibration & adjustment feature
- High chemical resistance for maximum durability
- Plug-and-Play PC connection
- Hybrid screen for excellent readability and use
- CalAuditTrail Provides gapless documentation (only for Entris® II BCA)

## Simplifying Progress

