S/RICH SILS

Cubis[®] II

The Only 100% configurable balance

- Customizable Modularity thousands of configuration possibilities
- Adaptable Software Integrated QApps
- Best-in-Class Performance monolithical weighing system
- Full End-to-End Data Integrity
- Intelligence by Design automatic motorized leveling

Simplifying Progress

