



sartorius

Cubis® High Capacity Micro Balance

Proven to deliver better reproducibility and faster measurement



40%

Cubis® vs. Main Competitor

Micro Balances are the perfect choice for drug eluting stent manufacturers in the Medical Devices industry; and results from a recent comparison study between the Cubis® High Capacity Micro Balance and an equivalent main competitor balance have shown that using the Cubis® delivers real benefits*.

Better Reproducibility

The Cubis® delivers outstanding reproducibility, delivering better weighing performance, even when measuring at less than minimum weights.

Repeatability performance of
Cubis® vs. Competitor X:

- Cubis® performed 40% better using a 2 g test weight (0.0011 mg vs. 0.0019 mg)
- Cubis® performed 20% better using a test weight near the maximum range of the balances (50 g) (0.0034 mg vs. 0.0043 mg)

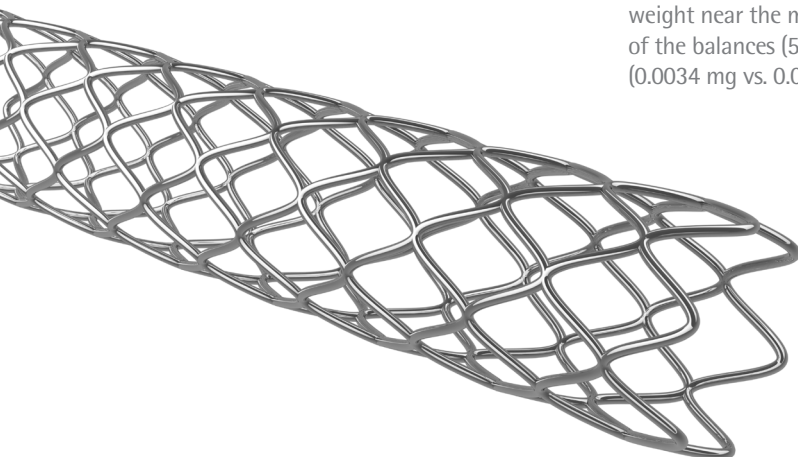
20%

Faster Measurement

The Cubis® delivers faster measurement thanks to short stabilisation times that improve efficiency when weighing larger lots of stock solutions.

Measurement stabilization time of
Cubis® vs. Competitor X:

- Cubis® performed 20% faster using a typical minimum weight (10.6 s vs. 12.1 s)



Learn more about Cubis® balances!
Get a free demo or a quote.
Visit: www.sartorius.com/md

*Source – Performance comparison between Sartorius Cubis MSA66S High Capacity Micro Balance and the main competitor balance. Comparison was conducted at Sartorius test lab in Göttingen under the same standard environmental conditions (draft, vibrations, humidity, electrostatic influences, etc.) with several repetitions using same weight.