

Save Costs, Material, and Time with Cubis® II

Calculate Your Savings

Cost \$/mg	Min ^{USP41} at your side	Typical Min ^{USP41} new Cubis® II UHR	Number of tests per year	Savings (\$) per year

Achieve constant minimum sample weights across the entire weighing range with the new Cubis® II Ultra-High Resolution Balances

What are Your Benefits?



Save on Costs

Reduce expensive material costs



Save on Material

Improve your sustainability through less consumption of material



Save Time

Measure faster with the new updated monolithic weighing system

How is the minimum sample weight calculated and why do you need it?

The USP places strict requirements on balances that are used for exact weighing in order to reduce faulty analyses in pharmaceutical processes to an acceptable and negligible level. These requirements are binding and are set out in USP chapter 41 on the topic of "Balances".

The requirement for repeatability restricts the weighing range to a working range usable for the corresponding purposes in the form of a minimum sample weight. The minimum sample weight is calculated from the standard deviation s of a repeatability measurement with at least 10 repetitions according to:

$$m_{\min, \text{USP}} = \begin{cases} 2\,000 \cdot s & \text{if } s \geq 0.41 \cdot d \\ 2\,000 \cdot 0.41 \cdot d & \text{otherwise} \end{cases}$$



Learn More:

