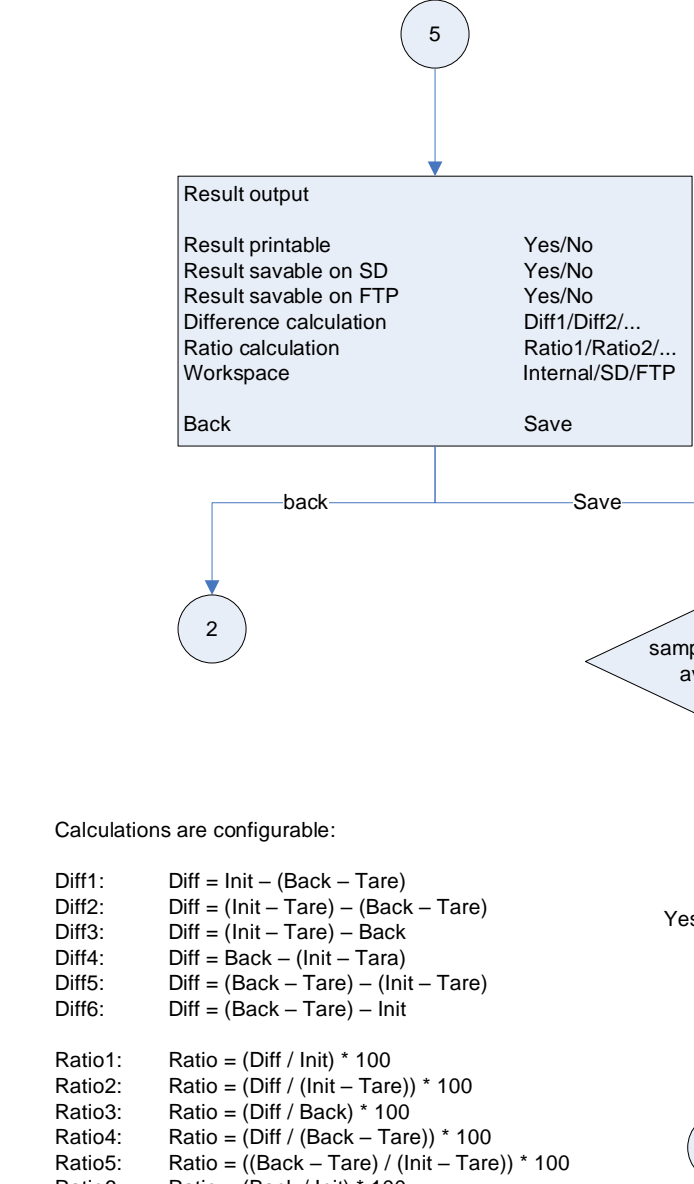


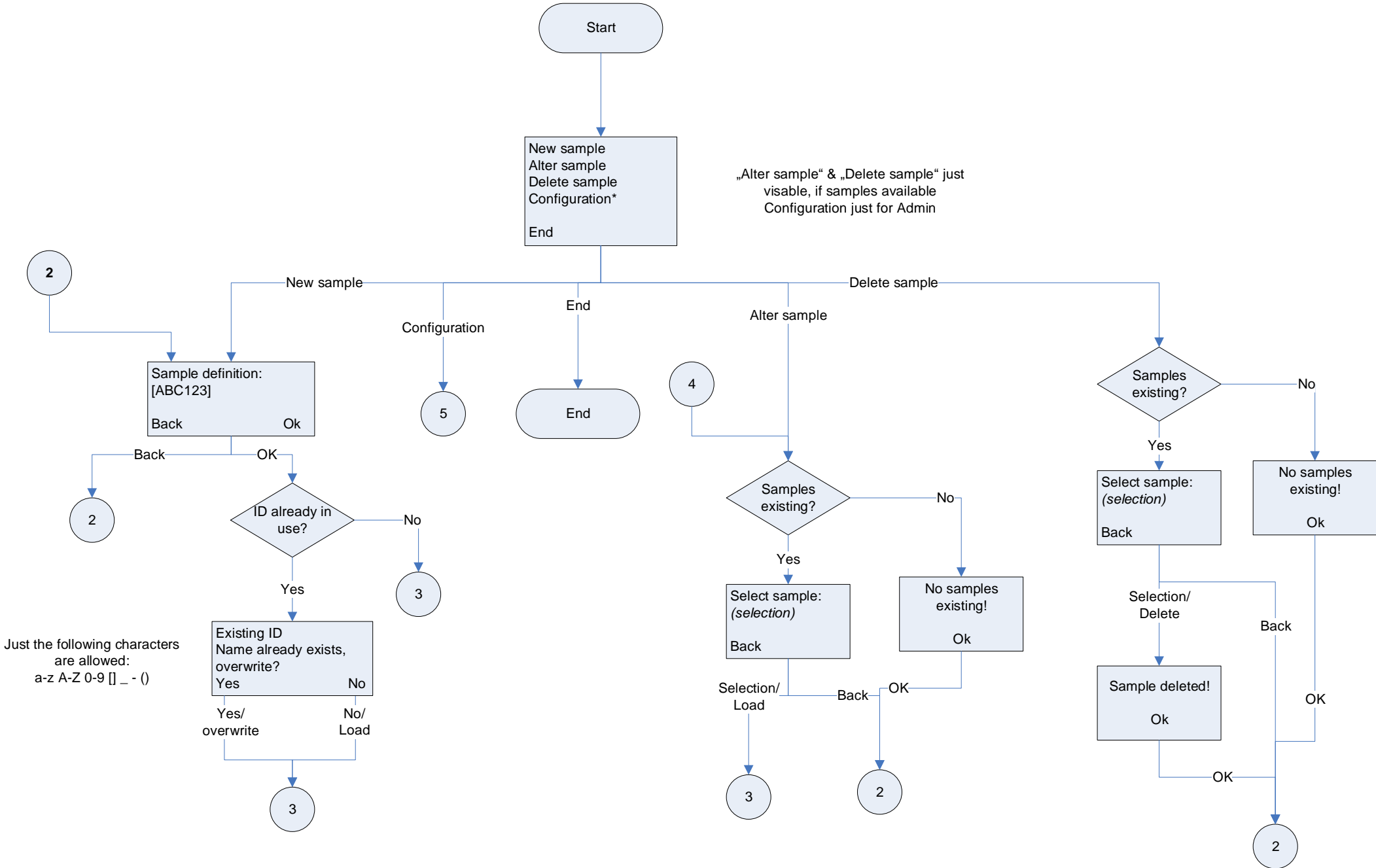
YAPP035
Unit: g, mg
System requirements:
Cubis Balance with MSA Unit, Firmware > 01.07
SD Card max. 32GB SDHC
Standard languages: German, English
Sample storage depends on configuration:
- internal file system
- SD Card
- FTP Server
Copy of files not possible!



Calculations are configurable:

Diff1: Diff = Init – (Back – Tare)  
Diff2: Diff = (Init – Tare) – (Back – Tare)  
Diff3: Diff = (Init – Tare) – Back  
Diff4: Diff = Back – (Init – Tare)  
Diff5: Diff = (Back – Tare) – (Init – Tare)  
Diff6: Diff = (Back – Tare) – Init

Ratio1: Ratio = (Diff / Init) \* 100  
Ratio2: Ratio = (Diff / (Init – Tare)) \* 100  
Ratio3: Ratio = (Diff / Back) \* 100  
Ratio4: Ratio = (Diff / (Back – Tare)) \* 100  
Ratio5: Ratio = ((Back – Tare) / (Init – Tare)) \* 100  
Ratio6: Ratio = (Back / Init) \* 100



Just the following characters  
are allowed:  
a-z A-Z 0-9 [ ] \_ - ( )

