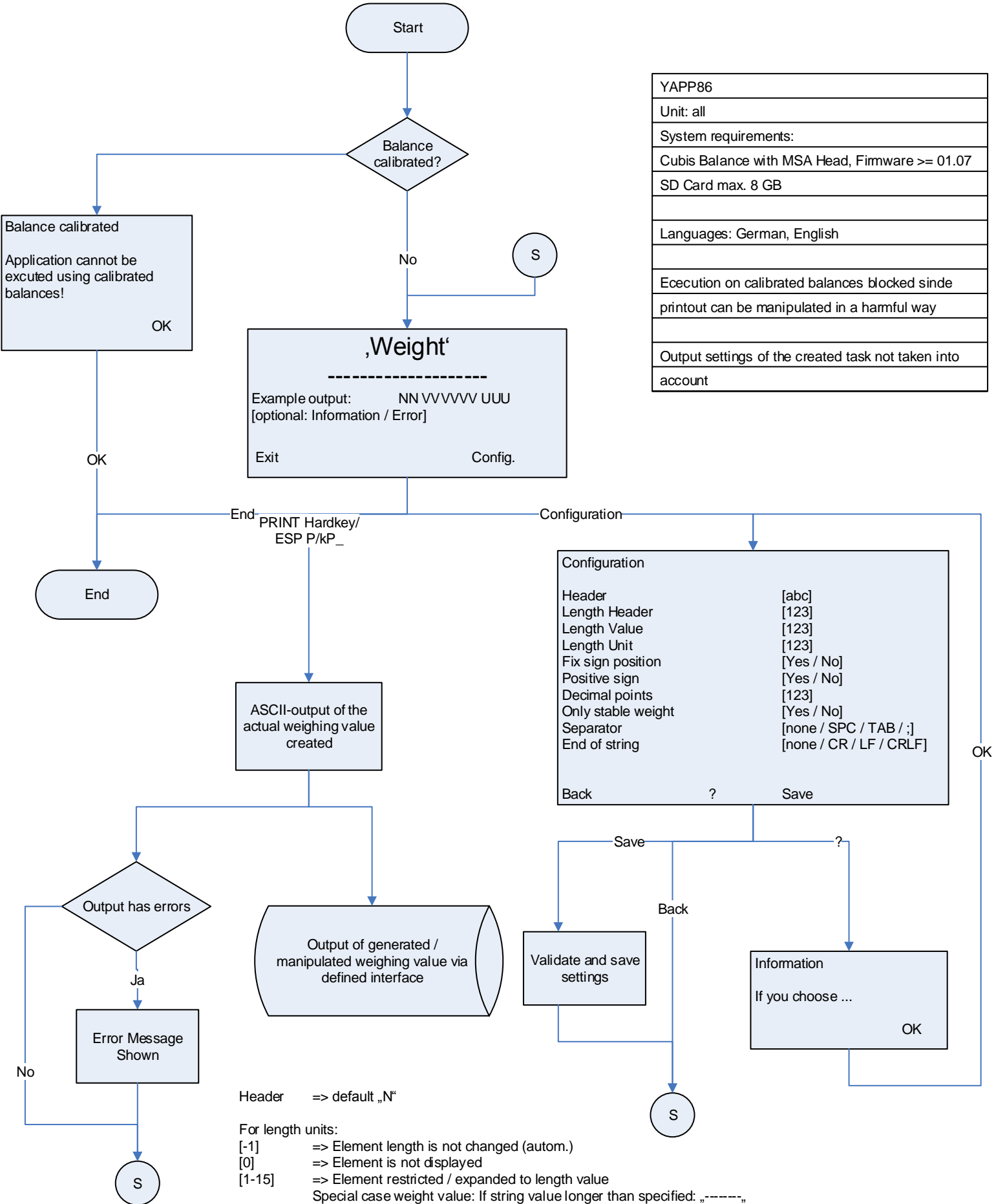


YAPP86
Unit: all
System requirements:
Cubis Balance with MSA Head, Firmware >= 01.07
SD Card max. 8 GB
Languages: German, English
Execution on calibrated balances blocked since printout can be manipulated in a harmful way
Output settings of the created task not taken into account



Header => default „N“

For length units:
 [-1] => Element length is not changed (autom.)
 [0] => Element is not displayed
 [1-15] => Element restricted / expanded to length value
 Special case weight value: If string value longer than specified: „-----“

Positive Sign: „+“ is shown if value is positive

Decimal points can only be reduced (min. [0])