

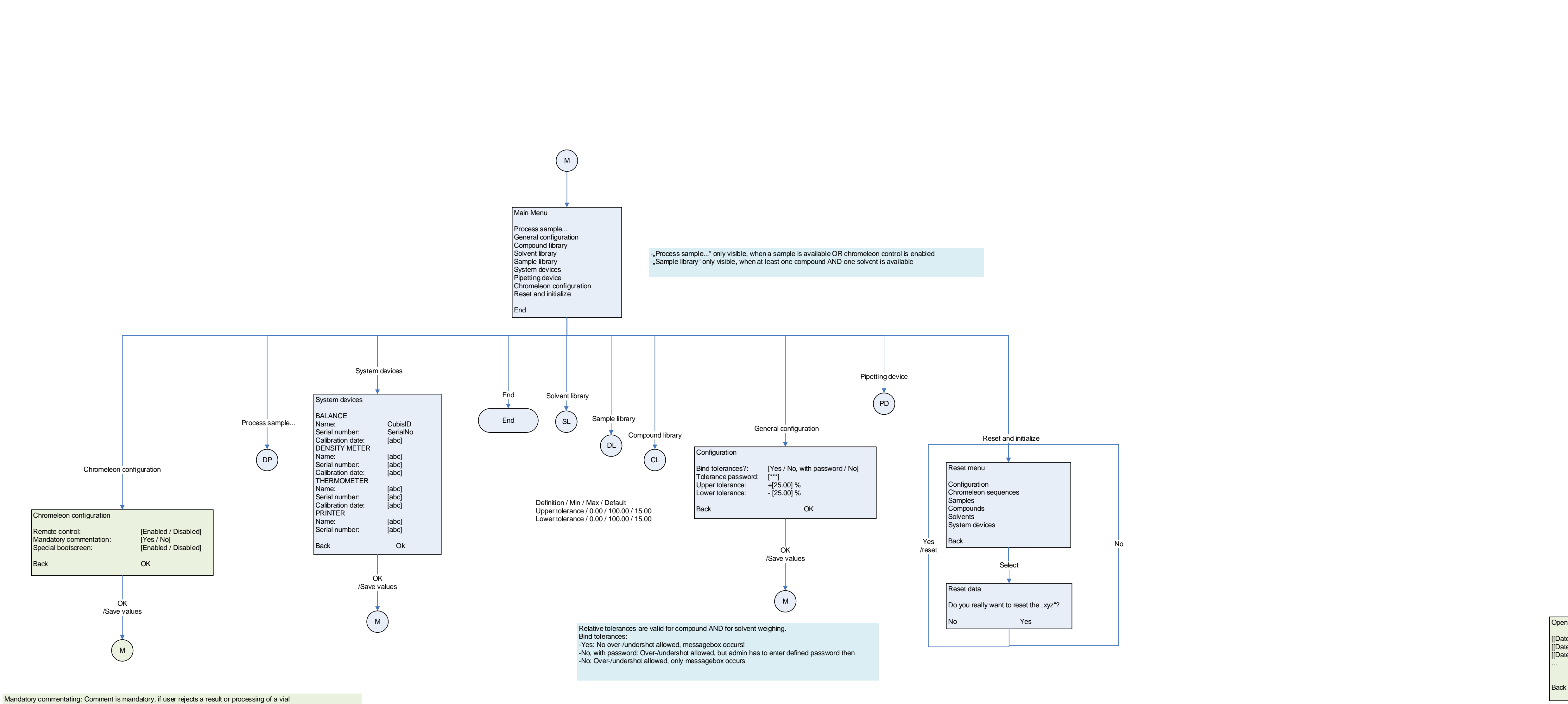
BASE PROCESS	VAPP10: Glass Solution Density
CHRONILETTE	Unit: g/ml; g/L
MANUAL/PFETTE	System measurements Custom labels with MSA (Unit: Amount x 10 ³ g/L) g/L (Unit: mol/L)
ORIGI TRIMATOR	Applicable with all methods Back: Custom labels ready to process screen
VICKIOLAR DISPENSER	Maximum 100 elements per compound/substance
CALCULATIONS	

Notes:

- Page 10 x volume of glass solution VAPP10 (unit: amount)
- Custom label print with three xk G/L
- Unit: g/ml; g/L; mol/L (Unit: amount x 10³ g/L)
- The main part of the Name is CHRONI, created by Chronometer
- In the report of the Name "T" appears, indicating the correct state of the sequence
- "N" = unprocessed compound from Chronometer
- "C" = processed compound from Chronometer
- "S" = processed (at least) by VAPP10, but not by Chronometer
- "D" = processed (at least) by Chronometer, but not by VAPP10
- "T" = processed (at least) by VAPP10, but not by Chronometer
- "U" = processed (at least) by Chronometer, but not by VAPP10
- "V" = processed (at least) by VAPP10, but not by Chronometer
- "W" = processed (at least) by VAPP10, but not by Chronometer

Data structure:

- 1 Chronometer sequence consists of 9 units
- 1 custom label consists of 1 unit x 3 positions
- 1 unit consists of 1 sample
- 1 sample consists of 100 measurements (100 measurements)
- 1 compound consists of 1 solvent and 1 or 2 substances
- 1 element consists of 1 solvent and 1 or 2 substances



Method comments: Comment is mandatory. User rights limit or processing of not selected parameter. To fully test the tool and identify screen the labels must be adapted to current ones

