

YAPR05 "Formulation"
Version: 2.0R
Q-App Version: 09-16-11
System requirements:
Cubis MSA, Firmware > 01.07.03
SD Card max. 32GB SDHC
Languages: German, English, French
Unit: mg, g, kg
Not compatible with verified balances!
Maximum 100 recipes in memory.
Maximum 10 compounds per recipe.
Back buttons usually lead to the previous screen.

- recipe\_id and comp\_id must not be empty
- recipe\_id can contain the following characters: a-z A-Z 0-9 | \_ - 0

- batch must not be empty and can contain the following characters: a-z A-Z 0-9 | \_ - 0

- Displayed weight value is overwritten with negative target weights (weighing to zero)

