

System requirements:
Cubis MSA, Firmware > 01.07.00
SD Card max. 32GB SDHC
Languages: English, German, French, Spanish
Unit: mg
A user must be logged in!
Compatible with verified balances (only positive weighings possible).

[illegible]