



YAPP14
Unit: n/a
System requirements:
Cubis Balance with MSA Unit, Firmware > 01.07.03
NTP Time Server
SD Card max. 32GB SDHC
Standard languages: German, English

Buttons marked with * are only available with Admin permissions.
During the workflow every hard key (except On/ Off) is blocked.