



Installation and Commissioning LevMixer® Order No.: S870AINST



The reliability and the service life of your equipment significantly depend on its installation and configuration according to the specifications.

Installation and commissioning by our service professionals will ensure that your LevMix-system successfully operates right from the start.

Sartorius Service installs and configures according to regulatory requirements and quality standards.

Specifications:

- Check of scope of delivery
- Check of device for transport damage
- Manufacturer's ID label check
- Installation of system components
- Start-up test
- Integration with customer mixing tank
- Brief instructions to operating personnel (no application training)
- General maintenance discussion

Customer site prerequisites:

- System present at installation location
- Unobstructed access to system
- Connection for power supply available at installation location
- If integration with customer mixing tank is requested, tank has to be provided at installation location
- Assistant available, e.g. for lifting and mounting heavy parts
- When ordering, a person in charge at site has to be appointed and communicated

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius-stedim.com

Specifications subject to change without notice.

Date: April 2013

