



YCP05-LR01

Liquid Reservoir

Advantages

- Internal temperature control: ensures error-free calibration measurements
- Triggers red alert, if temperature is out of tolerance; ensures compliance with regulations



Technical Specifications

Dimensions (L × W × H)	
With temperature sensor [mm]	200 × 55 × 80
Without temperature sensor [mm]	135 × 55 × 35
Volume [ml]	280
Weight [g] (container fully installed)	900
Temperature sensor measuring range [°C]	+15 – +30
Temperature sensor resolution [K]	0.1
Material of container	intermediate bottom Aluminum
Material of pipette guide	cover glass

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen, Germany
Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius.com

USA Toll-free +1.800.635.2906
UK +44.1372.737159
France +33.1.70.62.50.00
Italy +39.039.4659.1
Spain +34.913.586.095
Russian Federation +7.812.327.53.27
Japan +81.3.3740.5408

Specifications subject to change without notice.
Copyright Sartorius Stedim Biotech GmbH. Printed
in the EU on paper bleached without chlorine.
Publication No.: WL-2010-e160701
Order No.: 98649-020-96
Ver. 07 | 2016