

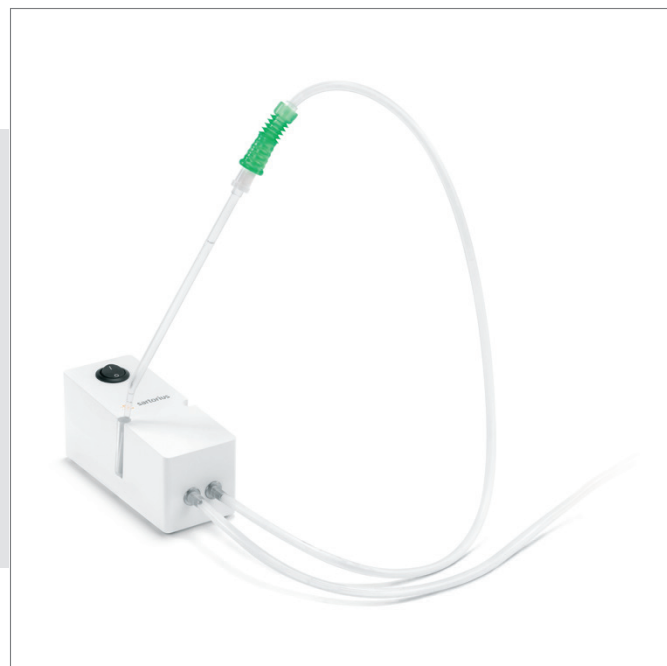


YCP05SP01

Suction Pump

Advantages

- Powerful
- Robust



Technical Specifications

Power Supply	
Type	Sartorius power supply module 6971790 with replaceable country-specific power plug adapters
Power supply, primary	Voltage V: 100 – 240 ± 10% Current A: 0.2 Frequency Hz: 50 – 60
Power supply, secondary	15 Vdc, +/- 5%, 530 mA (max.) 8 W (max.): 0 bis 40°C and 15 Vdc, +/- 5%, 330 mA (max.) 5 W (max.): 0 bis +50 °C
Installation site above sea level max. [m]	3000
Protection class according to EN/IEC 60950-1	II
Protection class according to EN/IEC 60529	IP40
Dimensions (L × W × H) [mm]	100 × 60 × 50
Weight [g]	90.0

Device	
Power supply	Only with Sartorius power supply module: 6971760
Input voltage VDC	15.0 ± 5%
Power consumption, average [W]	2
Media to be pumped, volume to be pumped Medium to be pumped	Ultrapure water
Volume to be pumped, max. ml/h	250
Dimensions Device (L × W × H) [mm]	150 × 55 × 70
Hoses (L) [mm]	600.0
Weight Device [g]	900.0
Hoses [g]	10.0

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-2011-e160701
Order No.: 98649-020-98
Ver. 07 | 2016