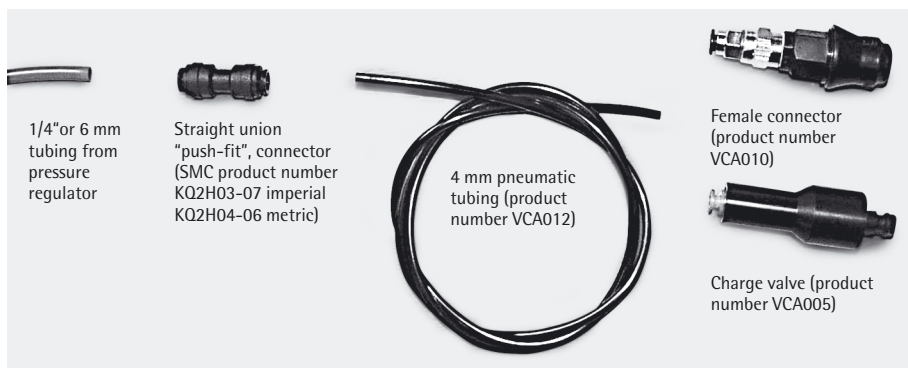


Directions for Use

Vivacell pressure products

Connection to alternative pressure sources





The majority of stirred-cell devices use 1/4" or 6 mm OD, pneumatic tubing as their inlet tube. Pressure regulators with this size of tubing can be used to pressurize Vivacell concentrators, with the addition of tubing adapters readily available from many pneumatic equipment suppliers.

Fitting instructions

Fit the 1/4" or 6 mm OD tubing from the stirred-cell pressure regulator into one side of a 1/4" - 5/32" or 6 mm - 4 mm straight union ,push-fit", connector.

Note

For 1/4" regulator tubing use SMC product number KQ2H03-07. For 6 mm regulator tubing use SMC product number KQ2H04-06. For distributors in your area please go to www.smcworld.com.

In the other side of this connector, insert standard 4 mm pneumatic tubing (product number VCA012), which is compatible with either of the recommended SMC straight union connectors. Cut this tube to the desired length; to the free end, attach the terminal connector appropriate to the Vivacell model you wish to pressurize. For the Vivacell 100 pressure head this will be the Female Connector (product number VCA010). For the Vivaspin® 20 pressure-fuge heads, you should use the Charge Valve (product number VCA005).

Operation

With all tube fittings in place, set the pressure regulator to the required pressure. Connect the tube fitting to the Vivacell or Vivaspin® concentrator. Continue with your concentration, following the operating instructions that accompany concentrator device. Here you can also find advice on suggested operating pressures, typical concentration times, and maximum pressure ratings.

Note

While the use of alternative pressure regulators with Vivacell products is possible, the Air Pressure Controller (product number VCA002) has been specially developed to compliment our range of concentrators. It includes safety features such as locking regulator controls and an integral overpressure relief valve to ensure easy and safe operation of Vivacell products.

Sartorius Stedim Biotech pressure products at a glance

Ordering information

Product	Pack size	Prod. no.
Air pressure controller (APC) complete with pressure gauge, regulator, over-pressure safety valve, female connector to Sartorius Stedim Biotech products, 1 m extension line (4 mm pneumatic tubing) with male and female connectors, and 1 m of 6 mm inlet tubing	1	VCA002
Female connector for Vivaspin® 20 pressure heads	1	VCA010
Male connector	1	VCA011
Charge valve for pressure heads	1	VCA005
4 mm pneumatic tubing (3 m)	1	VCA012
Vivacell 100 pressure head with replacement seals (5)	1	VCA800
Replacement seals for Vivacell 100 pressure head (VCA800)	10	VCA014
Vivaspin® 20 pressure head	1	VCA200

Sartorius Stedim Lab Ltd
Sperry Way
Stonehouse Park
Gloucestershire
GL10 3UT, UK

www.sartorius.com

Copyright by
Sartorius Lab Instruments
GmbH & Co. KG, Goettingen,
Germany.
All rights reserved. No part
of this publication may
be reprinted or translated in
any form or by any means
without the prior written
permission of Sartorius Lab
Instruments GmbH & Co. KG.
The status of the information,
specifications and illustrations
in this manual is indicated
by the date given below.
Sartorius Lab Instruments
GmbH & Co. KG reserves the right
to make changes to the technology,
features, specifications and design
of the equipment without notice.

Status:
December 2018,
Sartorius Lab Instruments
GmbH & Co. KG,
Goettingen, Germany

Printed in Germany on paper that has been
bleached without any use of chlorine
Vivacell Pressure Products
Publication No.: SLU6121-e181203