

# Masterflex<sup>®</sup> Peristaltic Pumps

Reliable Fluid Transfer for  
Vivaflow<sup>®</sup> Cassettes



## Benefits

- Precise, bi-directional flow control
- No risk of sample contamination
- Suitable for multiple fluid transfer applications
- Highly durable, low maintenance design

## Product Overview

Masterflex<sup>®</sup> peristaltic pumps perfectly compliment Vivaflow<sup>®</sup> crossflow filtration cassettes, whilst also offering the flexibility, durability and performance needed in multiple laboratory fluid transfer applications.

With samples confined to the tubing, there is no contact with the pump. This means there is no danger of contamination and no cleaning of the pump mechanism required between uses.

## Product Information

Masterflex® peristaltic pumps are positive displacement pumps. This means they are self-priming and non-siphoning, with no back flow. Unlike gear or piston pumps, peristaltic pumps can be used in a broader range of applications - including the transfer of liquids, gases or mixed-phase fluids - and the hard-wearing materials ensure long-term durability.

Since there is no sample contact with the pump mechanism, there is no risk of sample contamination. To avoid cross-contamination between different samples types, simply change the tubing - with the Easy Load pump head options, this takes a matter of seconds.

A variable speed, reversible motor enables easy sample transfer in either direction and the pumps are enabled for remote control of start | stop via a DB9 connector.

## Specifications

Product Number	VFP001	VFP002
Line voltage (50 - 60 Hz)	180 - 260 V	90 - 130 V
Drive speed range	20 - 600 rpm	20 - 600 rpm
Flow rate range	16 - 480 mL/min (size 16 tubing) 34 - 1,020 mL/min (size 15 tubing)	16 - 480 mL/min (size 16 tubing) 34 - 1,020 mL/min (size 15 tubing)
Power output	37 W	37 W
Maximum current	0.8 A	1.5 A
Dimensions (L   W   H)	230   183   134 mm	230   183   134 mm
Weight	4.1 kg	4.1 kg

## Masterflex Products at a Glance Ordering Information

Product	Pack Size	Order Number
Masterflex economy drive variable speed peristaltic pump (240 V)	1	VFP001
Masterflex economy drive variable speed peristaltic pump (115 V)	1	VFP002

Additional Equipment and Accessories	Pack Size	Order Number
Masterflex Easy Load pump head (size 16)	1	VFA012
Masterflex Easy Load pump head (size 15)	1	VFA013
Size 16 PVC pump tubing and luer fittings (3 m, 3.2 mm ID, 1.6 mm WT)	1	VFA004
Size 15 PVC pump tubing and luer fittings (3 m, 4.8 mm ID, 2.6 mm WT)	1	VFA003




**Germany**

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)