

Masterflex® Easy-Load® Pump Heads

Reliable Fluid Transfer for
Vivaflow® Cassettes

Benefits

- Quick and simple tubing changes
- Adjustment-free automatic tubing retention
- Compatible with multiple tubing sizes
- Highly durable, low maintenance design



Product Information

Masterflex® Easy-Load® pump heads perfectly complement Vivaflow® crossflow filtration cassettes, whilst also offering the flexibility, durability and performance needed in multiple other fluid-transfer applications.

Tubing changes are fast and easy, thanks to the single-thumb operated lever and automatic tubing retention. Two versions are offered, both compatible with multiple tubing sizes, extending their suitability to further applications, and the hard-wearing materials ensure long-term durability.

Technical Specifications

Order Number	VFA012	VFA013
Description	Easy-Load® pump head (size 16)	Easy-Load® pump head (size 15)
Dimensions (L W H)	64 105 105 mm	64 105 105 mm
Minimum flow rate	0.0012 mL/min	0.034 mL/min
Maximum flow rate	2,300 mL/min	1,700 mL/min
Housing material	Polyarylamide (PARA)	Polyarylamide (PARA)
Rotor material	SS* with PTFE coated bearings	SS* with PTFE coated bearings
Rollers	4x SS*	4x SS*
Tubing sizes	13, 14, 16, 25, 17, 18	15, 24
Optimum speed range	0.02 – 600 rpm	0.02 – 600 rpm
Volume per revolution		
Size 16 tubing	0.8 mL	-
Size 15 tubing	-	1.7 mL
Maximum vacuum, size 16 or 15 tubing	660 mm Hg	660 mm Hg
Suction lift, size 16 or 15 tubing	8.8 m H ₂ O	8.8 m H ₂ O
Maximum pressure, size 16 or 15 tubing	2.7 bar	2.7 bar

*SS= Stainless Steel

Ordering Information

Description	Pack size	Order no.
Masterflex® Easy-Load® pump head (size 16)	1	VFA012
Masterflex® Easy-Load® pump head (size 15)	1	VFA013

Additional Equipment and Accessories


Description	Pack size	Order no.
Masterflex® variable-speed peristaltic pump (240 V)	1	VFP001
Masterflex® variable-speed peristaltic pump (115 V)	1	VFP002
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