

PIXUL™ 96-well Plate Washer

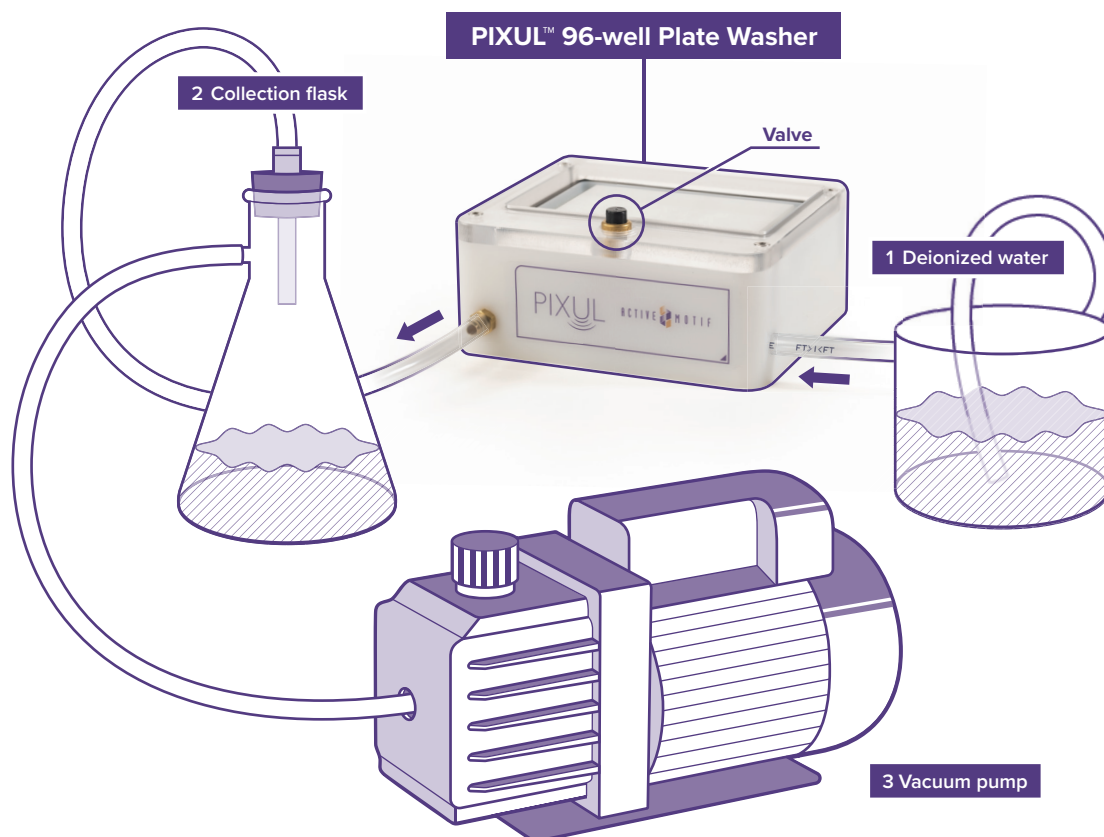
Cat. No. 53148

Additional Materials Required

- Vacuum pump
- Collection flask
- Deionized water
- Beaker

Instructions

1. Connect tubing to front port on PIXUL 96-well Plate Washer. This outgoing tubing will go to the **collection flask**.
2. Set up **beaker with deionized water** at the incoming tubing end, and a **collection flask** connected to a **vacuum pump** at the outgoing tubing end as shown in the illustration.
3. Fill PIXUL 96-well Plate Washer with **deionized water**.
4. Take the sonicated PIXUL 96-well Plate with Plate Sealer from a PIXUL sonication run and place it on top of the PIXUL 96-well Plate Washer.
5. Turn on the **vacuum pump**.
6. Press and hold the **valve button** on the PIXUL 96-well Plate Washer for approximately 10 seconds. Press gently on the PIXUL 96-well Plate to ensure a good seal. Water will flow under the PIXUL 96-well Plate to rinse any excess Coupling Fluid away from the sonication run.
7. When **valve button** is released, simply pull the PIXUL 96-well Plate off the PIXUL 96-well Plate Washer.



Technical Services

If you need assistance at any time, please call or send an e-mail to Active Motif Technical Service at one of the locations listed below.

Active Motif North America

Toll free: 877.222.9543
Direct: 760.431.1263
Fax: 760.431.1351
E-mail: tech_service@activemotif.com

Active Motif Europe

UK Free Phone: 0800/169 31 47
France Free Phone: 0800/90 99 79
Germany Free Phone: 0800/181 99 10
Direct: +32 (0)2 653 0001
Fax: +32 (0)2 653 0050
E-mail: eurotech@activemotif.com

Active Motif Japan

Direct: +81 (0)3 5225 3638
Fax: +81 (0)3 5261 8733
E-mail: japantech@activemotif.com

Active Motif China

Direct: (86)-21-20926090
Cell Phone: 18521362870
E-mail: techchina@activemotif.com

Visit Active Motif online at activemotif.com