



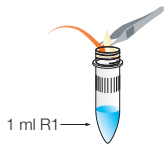
Aquadien DNA Extraction Kit, 3578121

Short Protocol for Dirty/Clogging Samples

Quick Guide



- Pipet 1 ml of R1 solution into a microtube
- Place a polycarbonate membrane filter on a sterilized filtration apparatus mounted on an air pump or vacuum flask
- Filter 100 ml–1 L of water



- Carefully fold the membrane in half 3 times to obtain a cone
- Using tweezers, place the membrane in the microtube containing 1 ml of R1 solution

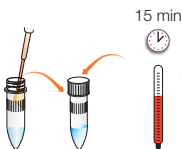


1,300 rpm

95°C

15 min

- Incubate at 95°C for 15 min at 1,300 rpm in a heating thermoshaker



15 min

4 ± 2°C



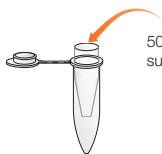
15 min

12,000 x g

4°C

- Carefully take out the membrane, pressing it to the walls of the tube to recover all the solution
- Add 100 µl of cool Aquadien W2 Wash Solution (catalog #3578119) and vortex for 5 sec
- Incubate at 4 ± 2°C for 15 min
- Centrifuge at 12,000 x g at 4°C for 15 min

The DNA is contained in the supernatant.



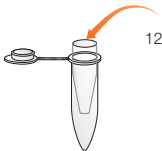
500 µl supernatant



6,000 x g

10 min

- Place a purification column in a collector vial
- Transfer 500 µl of the supernatant to the purification column
- Centrifuge at 6,000 x g for 10 min



125 µl R2

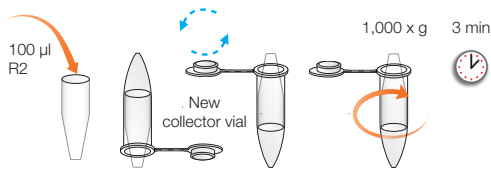


6,000 x g

5 min

- Add 125 µl of R2 solution
- Centrifuge at 6,000 x g for 5 min

continues



- Add 100 µl of R2 solution to the purification column; throw away the collector vial
- Cover the purification column with a clean collector vial and turn the whole unit upside down
- Centrifuge at 1,000 x g for 3 min
- Throw away the purification column



100 µl of purified DNA solution is obtained.

- Use 5 µl of the extracted DNA solution for real-time PCR analysis

For detailed instructions, review the kit user guide.

Visit bio-rad.com/legionella for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. AQUADIEN is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

