



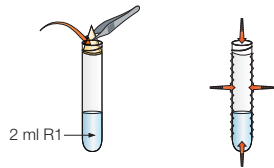
Aquadien DNA Extraction Kit, 3578121

Standard Protocol for Dirty/Clogging Samples

Quick Guide



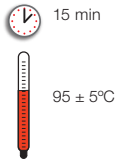
- Pipet 2 ml of R1 solution into a cryotube
- Place a polycarbonate membrane filter on a sterilized filtration apparatus mounted on an air pump or vacuum flask
- Filter 100 ml–1 L of water



20 sec

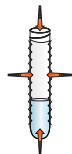


- Carefully fold the membrane in half 3 times to obtain a cone
- Using tweezers, place the membrane in the cryotube containing 2 ml of R1 solution
- Vortex for 20 sec



15 min

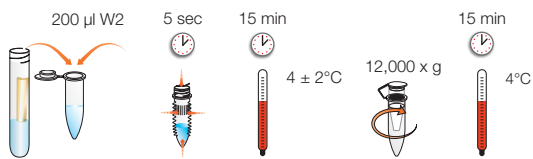
95 ± 5°C



20 sec



- Incubate at 95 ± 5°C for 15 min in a water bath
- Vortex for 20 sec



200 µl W2

5 sec

15 min

4 ± 2°C

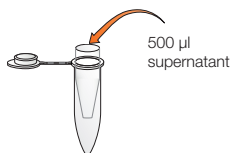
12,000 x g

15 min

4°C

- Transfer the sample (including the resin) to a 2 ml tube
- Add 200 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

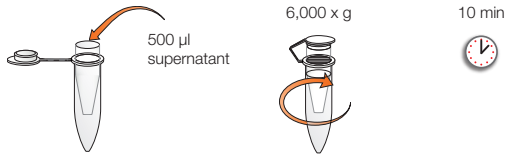


10 min

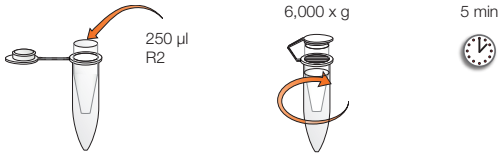


- Place a purification column in a collector vial
- Transfer 500 µl of the supernatant to the purification column
- Centrifuge at 6,000 x g for 10 min
- Remove column, empty collector vial, then replace column

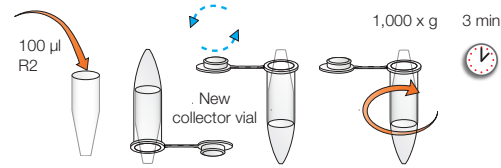
continues



- Transfer another 500 µl of the supernatant to the purification column
- Centrifuge at 6,000 x g for 10 min
- Remove column, empty collector vial, then replace column



- Add 250 µl of R2 solution
- Centrifuge at 6,000 x g for 5 min



- 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

