



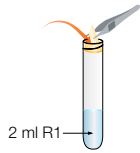
Aquadien DNA Extraction Kit, 3578121

Standard Protocol

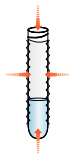
Quick Guide



- 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



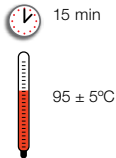
2 ml R1



20 sec

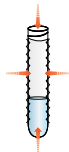


- Carefully fold the membrane in half 3 times to obtain a cone
- Using tweezers, place the membrane in the cryotube containing 2 ml of R1 solution
- Vortex for 20 sec



15 min

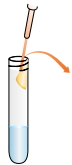
95 ± 5°C



20 sec



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

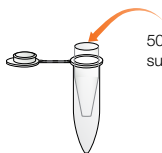


20 min



- Carefully take out the membrane, pressing it to the walls of the tube to recover all the solution
- Allow the cryotube to sit for 20 min at room temperature. The resin of the R1 reagent forms a pellet at the bottom of the cryotube

The DNA is contained in the 1.6 ml of supernatant



500 µl supernatant



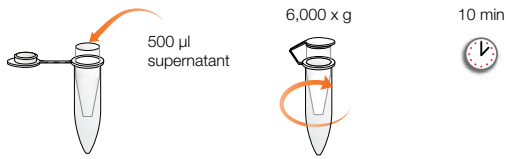
6,000 x g

10 min

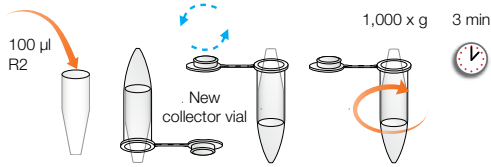


- Place a purification column in a collector vial
- Transfer 500 µl of the supernatant to the purification column
- Centrifuge at 6,000 x g for 10 min
- Empty the collector vial
- Remove column, empty collector vial, then replace column

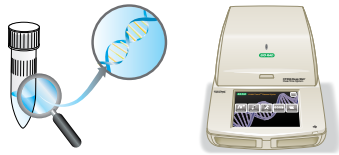
continues



- Transfer another 500 µl of the supernatant to the purification column
- Centrifuge at 6,000 x g for 10 min



- 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

**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

