
ProteinChip® 8-Well Bioprocessor User Guidelines

Catalog #C50-30008

Cleaning

- Soak the ProteinChip 8-well bioprocessor reservoir and gasket in a solution of 1% Triton X-100 in 1x PBS, pH 7.2 for at least 2 hours but not more than 24 hours
- Immediately before use, rinse the reservoir and gasket with deionized (DI) water to wash off the excess Triton X-100
- Air-dry the reservoir and gasket

Assembly

- Insert the ProteinChip array into the base with the logo extending beyond the base
- Mount the gasket on the ProteinChip bioprocessor reservoir, taking care to align the holes in the gasket with their corresponding pegs; the gasket fits in only one orientation
- Assemble the bioprocessor by lining up the holes in the base with the pegs on the reservoir
- Insert the assembly into the bracket to clamp the unit together snugly, but not too tightly; if the assembly is forced into the bracket, the gasket may stretch or be displaced, causing leakage and/or damage to the gasket

Leakage

- Test the ProteinChip bioprocessor for leakage by adding 550 µl of sample buffer to every other well
- Watch for leakage for 10 minutes. The meniscus should not fall during the 10 minutes. If there is leakage, there may be cross talk between adjacent wells and the bioprocessor should be dismantled and reassembled. If this fails to stop the leakage, the gasket may need to be replaced
- Drain the wells
- Repeat test for adjacent wells

Sample

- Load the sample into each well (50–150 µl/well)
- Seal the ProteinChip bioprocessor reservoir with cellophane tape or plate sealing tape to prevent sample evaporation

Incubation

- Secure the bioprocessor on the vortex plate; shake at 400–600 rpm

Note: Care should be taken to prevent the sample from splashing out of the wells. Too high a shake speed may cause sample to splash too high up the walls of each well.

Disassembly

- After incubation and washes are finished, drain each well with a pipet to prevent contamination of the wells
- To separate the reservoir and ProteinChip array assembly from the bracket, insert a solid cylindrical object, such as the base of a pen, into the hole at the back end of the bioprocessor and tap it gently on a hard surface. The reservoir and ProteinChip array assembly should loosen itself from the bracket
- While holding the ProteinChip array in place, carefully remove the reservoir and gasket
- Remove the ProteinChip array

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000