

InstaGene™ Matrix: Prepare DNA Templates for PCR* With No Phenol/Chloroform Extractions and No Deproteinization Steps

Prepare DNA from blood, cultured cells, or bacteria:

- Just incubate sample with InstaGene matrix at 56°C for 15–30 min, then boil for 8 min.
- InstaGene matrix absorbs cell lysis products, producing an improved substrate for PCR amplification.

DNA Preparation From Whole Blood

Use whole blood samples that are frozen, refrigerated, or collected fresh.

1. Add 3–6 µl of whole blood to 1 ml of autoclaved double distilled water in a 1.5 ml microfuge tube. Mix by inverting the tube several times.
2. Incubate the tube at room temperature for 15–30 min.
3. Spin at 10,000–12,000 rpm for 2–3 min.
4. Carefully remove all but 20–30 µl of the supernatant. Do not disturb the pellet.
5. Add 200 µl of InstaGene matrix to the pellet and incubate at 56°C for 15–30 min.

Note: InstaGene matrix should be mixed at moderate speed on a magnetic stirrer to maintain the matrix in suspension. The pipet tip used should have a large bore, such as on a 1,000 µl pipet tip (Bio-Rad's TBR tips, catalog #223-9350 and 223-9351).

6. Vortex at high speed for 10 sec. Place the tube in a 100°C heating block or boiling water bath for 8 min.
7. Vortex at high speed for 10 sec. Spin at 10,000–12,000 rpm for 2–3 min.
8. Use 20 µl of the resulting supernatant per 50 µl PCR reaction. Store the remainder at -20°C. Repeat step 7 when reusing the InstaGene DNA preparation.

Note: It is important to store the prepared sample at -20°C when not in use.

DNA Preparation From Cultured Mammalian Cells

1. Pellet 200 µl of cells suspension from media in a microfuge tube. Spin at 10,000–12,000 rpm for 1 min and remove the supernatant.
2. Resuspend cells in 1 ml of 1x PBS and centrifuge for 1 min at 10,000–12,000 rpm.
3. Resuspend cells in autoclaved water at 20–30 cells/µl.
4. Add 20 µl of this cell suspension to 200 µl of InstaGene matrix. Incubate at 56°C for 15–30 min.

Note: InstaGene matrix should be mixed at moderate speed on a magnetic stirrer to maintain the matrix in suspension. The pipet tip used should have a large bore, such as on a 1,000 µl pipet tip.

5. Follow the procedure for DNA preparation from whole blood, steps 6–8.

DNA Preparation From Bacteria

The protocol described below is for the preparation of genomic DNA or episomal DNA from bacteria.

1. Pick an isolated bacterial colony and resuspend it in 1 ml of autoclaved water in a microfuge tube.
2. Centrifuge for 1 min at 10,000–12,000 rpm. Remove the supernatant.
3. Follow the DNA preparation procedure for whole blood, steps 5–8.

Ordering Information

Catalog #	Description
732-6030	InstaGene Matrix, 20 ml, 100 DNA preps

* The polymerase chain reaction (PCR) process is covered by patents owned by Hoffman-LaRoche. Practice of the PCR process requires a license.



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191 Canada (905) 712-2771 China (86-21) 63052255 Czech Republic (420) 2-4141 0532 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124) 6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30 Poland (48) 22-8126 672 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-2729877 South Africa 00 27 11 4428508 Spain 590 5200 Sweden 46 (0)8-55 51 27 00 Switzerland 061 717-9555 United Kingdom 0800-181134