

# InstaGene Matrix: Prepare DNA Templates for PCR Without Phenol or Chloroform Extractions and Without Deproteinization Steps

Prepare DNA from blood, cultured cells, or bacteria:

- Incubate the 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 Laboratories, Inc, catalog #2239350 and 2239351).

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

Visit [bio-rad.com/InstaGene](https://www.bio-rad.com/InstaGene) for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc.  
All trademarks used herein are the property of their respective owner.  
© 2026 Bio-Rad Laboratories, Inc.



**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** 55 11 3065 7550  
**Canada** 1 800 361 1808 **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 **Greece** 30 210 7774396 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23  
**India** 91 124 4029300 **Israel** 000 800 00 24 67 23 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23  
**Mexico** 52 55 5488 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** 7 495 721 14 04 **Singapore** 65 6415-3170 **South Africa** 27 21 531 7504 **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** 662 651 8311 **United Arab Emirates** 971 4 818 7300  
**United Kingdom** 00 800 00 24 67 23