

# nucleic acid purification

## INSTAGENE™ MATRIX AND KITS

- **QUALITY-CONTROLLED CHELATING MATRIX FOR REPRODUCIBLE PCR RESULTS**
- **PCR-READY DNA FROM TISSUES AND CELLS IN 30-45 MINUTES**
- **INEXPENSIVE — NO SPIN COLUMNS OR MEMBRANES REQUIRED**
- **CONVENIENT KITS FOR CELLS AND TISSUE, WHOLE BLOOD AND DRY BLOOD**

## PCR\* Ready DNA in the Palm of Your Hand

### InstaGene Matrix

- Specially formulated, quality-controlled Chelex® resin for superior PCR results
- Reproducible from lot to lot
- Offered in ready-to-use, smaller particle slurry that won't clog pipette tips
- Fast and affordable; takes only 45 minutes

### InstaGene Whole Blood Kit

- Formulated and quality-controlled for optimum results with whole blood
- Special buffer speeds up cell lysis step
- Works with commonly used anticoagulants such as heparin, citrate, EDTA
- Fast and affordable; takes only 30 minutes

### InstaGene Genomic DNA Kit

- Yields consistent 50 Kb or larger DNA from whole blood
- Ideal for long PCR and Southern blotting
- Only 30 minutes of hands-on time
- Works with commonly used anticoagulants such as heparin, citrate, EDTA

### InstaGene Dry Blood Kit

- Optimum results with dried blood on filter paper
- Works well with 1/8" or 1/4" paper punches
- Fast and affordable; takes only 45 minutes



**BIO-RAD**

## InstaGene Matrix

Treatment of samples with Chelex 100 is the standard procedure for PCR sample preparation from bacteria, whole blood, mammalian cells and animal tissue. The InstaGene matrix makes Chelex easier to handle and is QC'd to ensure reproducibility from lot to lot. It's a ready-to-use slurry made with a small particle size for simple pipetting, and includes a stirring bar for easy resuspension.

## Genomic DNA Kit

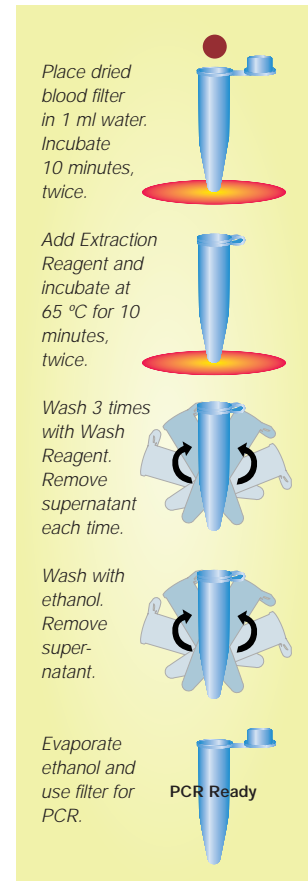
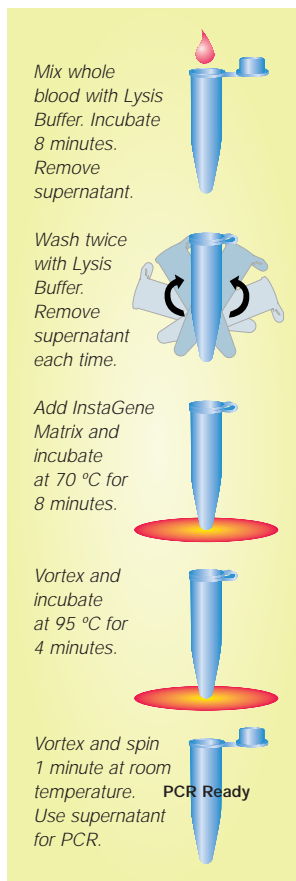
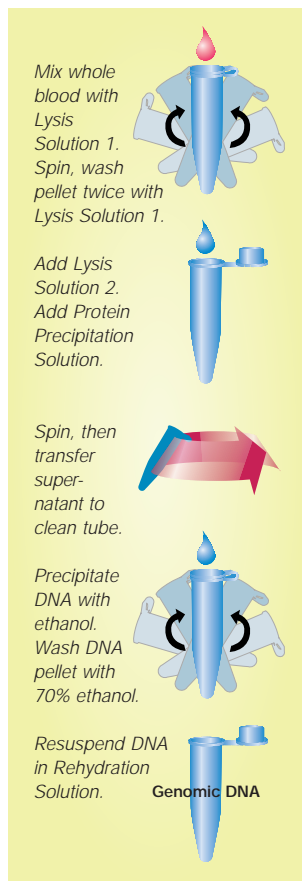
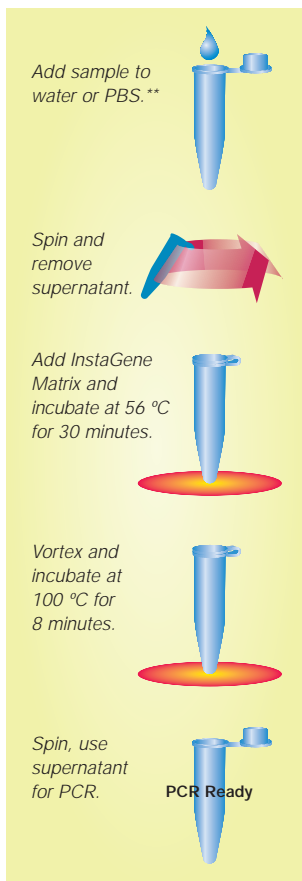
The molecular diagnostics laboratory is often interested in preparing DNA from blood that can be used for different diagnostic procedures, including short and long PCR and Southern blotting. These applications require a method that consistently yields pure DNA over 50 Kb in size. The InstaGene genomic DNA kit meets this criteria quickly and efficiently, requiring only 30 minutes of hands-on time.

## Whole Blood Kit

In order to increase the efficiency of the cell lysis step, Bio-Rad has designed a lysis buffer to pair up with the ready-to-use InstaGene matrix. The InstaGene whole blood kit offers an extremely efficient and fast method for preparing PCR-ready DNA templates from whole blood. The protocol for whole blood is presented below to illustrate the speed and simplicity of this method.

## Dry Blood Kit

While the whole blood kit contains the InstaGene matrix, the InstaGene dry blood kit does not. Instead, an extraction and a washing reagent are supplied. The DNA remains bound to the filter paper throughout the procedure and is amplified in situ. The resulting protocol is fast and simple, as illustrated below.



\* The Polymerase Chain Reaction (PCR) process is covered by patents owned by Hoffman-LaRoche.

\*\* See manual for specific protocols.

## Ordering Information

Catalog #	Description
732-6030	InstaGene Matrix, 20 ml (enough for 100 samples)
732-6028	InstaGene Genomic DNA Kit, one bottle each of Lysis Solution 1, Lysis Solution 2, Protein Precipitation Solution, Rehydration Solution (enough for 100 samples when using 300 µl of blood)
732-6211	InstaGene Whole Blood Kit, includes one bottle of InstaGene Matrix and one bottle of InstaGene Lysis Buffer (enough for 100 samples)
732-6212	InstaGene Dry Blood Kit, includes one bottle of Extraction Reagent and one bottle of Wash Reagent (enough for 125 samples)

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

Website [www.bio-rad.com](http://www.bio-rad.com) U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771 China 86-10-62051850 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France (01) 43 90 46 90 Germany 089 318 84-0 Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 Korea 82-2-3473-4460 The Netherlands 31 318-540666 New Zealand 64-9-4152280 Singapore 65-2729877 Spain (91) 661 70 85 Sweden 46 (0)8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800-181134