

NUCLEIC ACID PURIFICATION

PCR Kleen™ Spin Columns

- Purify PCR products >200 bp from reaction mixture components in 4 min
- Provide purified DNA ready for downstream applications
- Accommodate sample loads of 25–100 µl
- Require no matrices or binding steps

Rapid PCR Product Purification

Amplification of DNA by the polymerase chain reaction (PCR) is a central technique in many molecular biology laboratories. Purification of the initial DNA template is a critical first step. PCR Kleen spin columns are a proven, superior choice for elimination of PCR impurities. These prepacked spin columns allow convenient purification of DNA and PCR products >200 bp directly from reaction mixtures. An easy, rapid 4 min protocol eliminates nucleotides (dNTPs, ddNTPs), enzymes (nucleases, polymerases), ethanol, salts, primers, and primer-dimers <32 bp in length. Use of the columns also eliminates the need to cut DNA fragments out of agarose gels.

Convenient Method

PCR Kleen spin columns eliminate the need for time-consuming matrix binding and wash buffer steps, gravity columns, dialysis membranes, electroelution devices, phenol extraction, and alcohol precipitation to isolate PCR products.

Rapid Procedure

Simply transfer 25–100 µl of the amplification mixture to a PCR Kleen spin column and centrifuge in a microcentrifuge for 3 min. Purified DNA fragments are eluted into the collection tube, eliminating contaminants for easy disposal.

Efficient Separation

The columns rapidly recover DNA >200 bp from reaction mixture components (enzymes, nucleotides, primers, and primer-dimers <32 bp in length) that can decrease downstream cloning efficiencies (Figure 1).

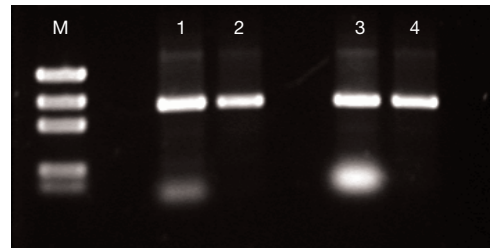


Fig. 1. Effective, rapid removal of primer-dimers from PCR reactions. Two PCR reaction mixtures were purified with PCR Kleen spin columns. Lane 1, a mixture of 2 µg of a 660 bp DNA fragment with 100 ng of a 25-mer primer-dimer; lane 3, a mixture of 2 µg of a 660 bp DNA fragment with 500 ng of a 25-mer primer; lane 2, primer-dimer purification; lane 4, primer purification; M, DNA markers. Reaction products and purified samples (10 µl each) were analyzed on a 1.5% agarose gel.

Effective Results

Purified DNA is immediately available for PCR, subcloning, restriction digestion, ligation, nucleotide sequencing, and other downstream applications.

BIO-RAD

Specifications

Method	Size exclusion chromatography in spin column
Application	Removal of nucleic acids <32 bases and other reaction mixture components from PCR products >200 bp
Capacity	25–100 µl
Equilibration buffer	10 mM Tris, 1 mM EDTA, pH 7.4
Sample preparation time	<4 min
Yield	50–80% recovery

Ordering Information

Catalog #	Description
732-6300	PCR Kleen Spin Columns, 25

For more information, request bulletin 2950, or go to www.bio-rad.com/nasampleprep/

Practice of the polymerase chain reaction (PCR) may require a license.

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