

PCR Purification Kit

DNA Purification by silica-gel membrane adsorption

DNA Cleanup

Cat.-No.	Amount
PP-201S	50 preparations
PP-201L	250 preparations

For *in vitro* use only
 Quality guaranteed for 12 months
 Store at room temperature

Kit contents

Binding Buffer

Washing Buffer (before use, add 96-99% Ethanol as indicated on the bottle)

Elution Buffer

Spin Columns

2 ml Collection Tubes

To be provided by you

96-99% Ethanol

Isopropanol (for high yield sample preparation)

1.5 ml microtubes

Description

PCR Purification Kit is designed for the work-up of PCR reactions (removal of primer dimers, primers, nucleotides, proteins, salt, agarose, ethidium bromide, and other impurities). The preparation is based on a silica-membrane technology for binding DNA in high-salt and elution in low-salt buffer. The kit provides a simple and efficient way to purify linear or circular DNA and is optimized for working with DNA amounts of up to 50 µg. It requires no organic extractions or precipitation and guarantees high and stable recovery rates.

Preparation procedure

The DNA purification follows a simple binding, washing, and eluting procedure. Before start, add 96-99% Ethanol as indicated on the bottle to the Washing Buffer. The additional use of Isopropanol is recommended for fragments smaller than 200 bp or larger than 5 kbp.

1a Standard Sample Preparation

For DNA fragment sizes in the range of 200 bp to 5 kbp:

- Add 5 volumes of *Binding Buffer* to 1 volume of DNA sample and mix well. For example, if the volume of your DNA sample is 50 µl, add 250 µl Binding Buffer.

1b High Yield Sample Preparation

For DNA fragment sizes smaller than 200 bp or larger than 5 kbp:

- Add 3 volumes *Binding Buffer* and 2 volumes of Isopropanol to the PCR sample. For example, if the volume of your DNA sample is 50 µl, add 150 µl Binding Buffer and 100 µl *Isopropanol*.

2. Column Loading

- Place a *Spin Column* into a 2 ml collection tube
- Apply the sample mixture from step 1 into the *Spin Column*.
- Centrifuge at 10,000g for 30 sec in a micro-centrifuge.
- Discard the flow-through.

4. Column Washing

- Add 750 µl of *Washing Buffer* to the *Spin Column*.
- Centrifuge at 10,000 g for 2 min.
- Discard the flow-through.

5. Elution

- Place the *Spin Column* into a clean 1.5 ml microtube (not provided in the kit)
- Add 30-50 µl *Elution Buffer* or dd-water to the center of the column membrane
- Incubate at room temperature for 1 min.
- Centrifuge at 10,000 g for 1 min to elute DNA.