

ab156894

PCR Clean-up Kit

Instructions for Use

For efficient clean-up of PCR products from various sources.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

Table of Contents	1
1. Background	2
2. Assay Summary	3
3. Materials Supplied	3
4. Storage and Stability	4
5. Materials Needed Not Supplied	4
6. Reagent Preparation	4
7. Assay Procedure	4

1. Background

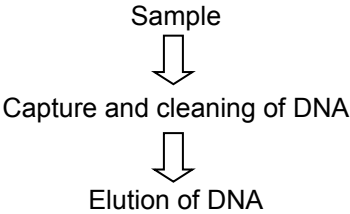
The PCR Clean-up Kit uses a unique procedure and composition to efficiently clean-up PCR products from various sources including conventional or real time thermal cycling PCR and isothermal gene amplification (PCR). The kit has the following features:

- The fastest procedure available (< 2 minutes).
- Ultra-purity and recovery of small PCR products with efficient removal of contaminants (salts, proteins, enzymes, nucleotides, and primers).
- Use of preparatory binding solution and non chaotropic or other irritant reagents.

The PCR Clean-up Kit allows cleaning of PCR product sizes from 200 bp to 10 kb; DNA quantity from 0.1 ng to 2 µg, optimal at between 1 ng and 1 µg. Recovery rate of DNA is greater than 70%. Cleaned PCR product is ready for direct use in the following applications: ligation/transformation, sequencing, restriction digestion, microarray analysis, labeling, microinjection, and *in vitro* transcription.

ab156894 simply applies our proprietary binding buffer to PCR products. After transferring the sample to the specially designed F-Spin Column and washing with the washing buffer, the PCR products are easily recovered in 8-20 µl elution buffer. Purified PCR products are ready for down-stream application.

2. Assay Summary



3. Materials Supplied

Item	50 tests	100 tests	Storage (Before Preparation)
PC1 (Binding Buffer)	16 ml	2 x 16 ml	RT
PC2 (Wash Buffer)	2.5 ml	2 x 2.5 ml	RT
PC3 (DNA Elution Solution)	1 ml	2 ml	RT
F-Spin Column	50	100	RT
F-Collection Tube	50	100	RT

4. Storage and Stability

PCR Clean-up Kit can be stored at room temperature (15-22°C).

5. Materials Needed Not Supplied

- Pipette
- 100% ethanol
- Microcentrifuge capable of 12000 rpm

6. Reagent Preparation

Prepare PC2 (Final Wash Buffer):

Add 20 ml of 100% Ethanol to each bottle of PC2 (Wash Buffer)

7. Assay Procedure

Note: Always cap spin columns before placing them in the microcentrifuge.

- a) Add 2 volumes of PC1 (Binding Buffer) to each volume of the PCR product sample, and mix well. (For example, add 200 μ l of PC1 (Binding Buffer) to 100 μ l of PCR product sample. If required volume is less than 100 μ l, add 100 μ l of PC1 (Binding Buffer).)
- b) Transfer mixture to the column. Centrifuge at 12,000 rpm for 15 seconds. Discard the flowthrough. Replace the column to the collection tube

Note: *maximum volume of the column is 600 μ l*

- c) Add 200 μ l of the Modified PC2 (Final Wash Buffer) to the column and centrifuge at 12,000 rpm for 15 seconds. Discard the flowthrough and replace the column to the collection tube.
- d) Add 200 μ l of the Modified PC2 (Final Wash Buffer) to the column and centrifuge at 12,000 rpm for 30 seconds.
- e) Place the column in a new 1.5 ml vial. Add 8-18 μ l of PC3 (DNA Elution Solution) directly to the column filter, centrifuge at 12,000 rpm for 20 seconds to elute cleaned DNA.
- f) DNA is now ready for use or storage at -20°C .

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)