

TuMinute™ PCR Clean-Up Kit

Base Catalog # P-1005

PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The TuMinute™ 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.

KIT CONTENTS

Components	50 samples P-1005-1	100 samples P-1005-2
PC1 (Binding Buffer)	16 ml	2 x 16 ml
PC2 (Wash Buffer)	2.5 ml	2 x 2.5 ml
PC3 (DNA Elution Solution)	1 ml	2 ml
F-Spin Column	50	100
F-Collection Tube	50	100

SHIPPING & STORAGE

The *TuMinute*™ PCR Clean-Up Kit can be stored at room temperature (15-22°C) for up to 1 year.

GENERAL PRODUCT INFORMATION

Quality Control: EpigenTek guarantees the performance of all products in the manner described in our product instructions.

Product Updates: EpigenTek reserves the right to change or modify any product to enhance its performance and design. The information in this User Guide is subject to change at any time without notice. Be sure to use the latest User Guide for this kit which can be accessed online at www.epigentek.com/datasheet.

Usage Limitation: The *TuMinute*™ PCR Clean-Up Kit is for research use only and is not intended for diagnostic or therapeutic application.

Intellectual Property: *TuMinute*™ is a trademark of EpigenTek, Inc.

A BRIEF OVERVIEW

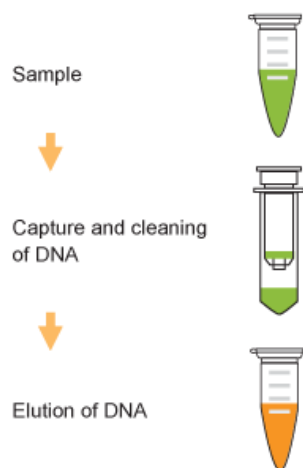
The *TuMinute*™ PCR Clean-Up Kit uses a unique procedure and composition to efficiently clean-up PCR products from various sources including conventional or real time thermal cycling PCR and isothermal gene amplification (PCR). The kit has the following features:

- 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.

PRINCIPLE & PROCEDURE

The *TuMinute*™ PCR Clean-Up Kit simply applies our proprietary binding buffer to PCR products. After transferring the sample to the specially designed Fast-Spin Column and washing with the

washing buffer, the PCR products are easily recovered in 8-20 μ l elution buffer. Purified PCR products are ready for down-stream application.



Schematic Procedure for Using the *TuMinute*™ PCR Clean-Up Kit

PROTOCOL

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

Before starting, add 20 ml of 100% ethanol to PC2 to make the final wash buffer.

1. Add 2 volumes of **PC1** to each volume of the PCR product sample, and mix well. (For example, add 200 μ l of **PC1** to 100 μ l of PCR product sample. If required volume is less than 100 μ l, add 100 μ l of **PC1**.)
2. 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.*)
3. 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. Add 200 μ l of the modified **PC2** (*final wash buffer*) to the column and centrifuge at 12,000 rpm for 30 seconds.
4. Place the column in a new 1.5 ml vial. Add 8-18 μ l of **PC3** directly to the column filter, centrifuge at 12,000 rpm for 20 seconds to elute cleaned DNA.

DNA is now ready for use or storage at -20°C .

RELATED PRODUCTS

P-1006 DNA Concentrator Kit