



371PR-02

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

PinkCLEANUP™

DNA Cleaning Kit

(Cat. # 786-87)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION 3

ITEM(S) SUPPLIED 3

STORAGE CONDITIONS 3

PREPARATION BEFORE USE 3

PROTOCOL 4

SPIN COLUMN PROTOCOL 5

APPLICATION NOTES 6

RELATED PRODUCTS 6

INTRODUCTION

The pinkCLEANUP™ kit uses G-Biosciences's revolutionary pinkRESIN™ to remove excess salts, enzymes, unincorporated nucleotides, and primer-dimers from DNA preparations. DNA fragments larger than approximately 100 base pairs are isolated. No toxic phenol/chloroform extractions or alcohol precipitations are needed. Because pinkRESIN™ has enhanced binding properties and the ability to resuspend easily, protocols take as little as 5-10 minutes. Ideal applications include PCR clean up and restriction enzyme removal from plasmid DNA prior to in-vitro transcription. Instructions are included for users wishing to use spin columns (Catalog # 786-88).

The kit is designed for 200 x 200µl samples containing <10ug DNA/ml.

ITEM(S) SUPPLIED (Cat # 786-87)

Description	Size
GET™ Plasmid Binding Buffer	2 x 30ml
DNA Wash	50ml*
pinkRESIN™	2 x 1ml
TE buffer	1 x 10ml
Wide bore pipette tips (Small)	200

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Store refrigerated at room temperature upon arrival, and is stable for 1 year.

PREPARATION BEFORE USE

- Wash preparation: Add 200ml absolute ethanol to DNA Wash. DNA WASH should be cooled on ice before use (alternatively, store wash at -20°C).
- Preheat TE buffer to 55-60°C

PROTOCOL

1. Add equal volume of Binding Buffer to the sample.
2. Suspend pinkRESIN™ by vigorous vortexing and add 10µl of the resin per 200µl of the sample using a wide bore pipette tip.
NOTE: *If the sample contains more than 2µg DNA, add an additional 5µl pinkRESIN™ for each 1µg DNA.*
3. Incubate on ice for 1-3 minutes and then centrifuge for 30 seconds at high speed in a microcentrifuge.
4. Remove the supernatant by decanting or pipetting. Use the small remaining volume (20-30µl) in the tube to resuspend the pinkRESIN™ pellet. Spin column users continue to Spin Column Protocol.
5. Add 500µl ice cold DNA Wash and fully resuspend the pinkRESIN™. Centrifuge for 30 seconds at high speed. Decant or pipette out the wash. Repeat the wash step one more time and then allow the pellet air dry for 1-2 minutes or until all traces of the wash are gone.
6. Add pre-warmed (55-60°C) TE buffer or DNase-free water to the microcentrifuge tube and resuspend the pinkRESIN™ pellet.
7. Incubate at room temperature for 1-2 minutes. The volume can be equal to or less than the original sample volume but should be at least twice the original volume of pinkRESIN™. Centrifuge 10-30 seconds to pellet the resin and transfer the sample containing cleaned DNA to a clean tube.

SPIN COLUMN PROTOCOL

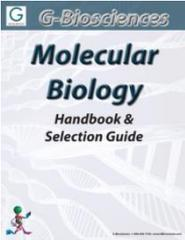
1. Add 300µl DNA Wash to tube and fully resuspend the pinkRESIN™.
2. Insert a micro spin column into a 1.5 or 2 ml centrifuge tube.
3. Add the slurry into the spin column and centrifuge for 10-30 seconds at high speed.
4. Add 300µl DNA Wash to the tube and resuspend any remaining resin. Add this to the spin column and spin 10-30 seconds.
5. Discard the first two washes and add 600µl DNA Wash to the spin column and centrifuge. Drain the lower tube. A final spin is recommended to remove any residual wash.
6. Place the spin column in a clean collection tube. Add pre-warmed (55-60°C) TE buffer or molecular grade water to the spin column and swirl to resuspend the resin. The volume can be equal to or less than the original sample volume but should be at least twice the original volume of pinkRESIN™.
7. Centrifuge for 10-30 seconds to elute the DNA from the pinkRESIN™.

APPLICATION NOTES

1. For DNA fragments or plasmids larger than 5kb increase TE buffer incubation to 5-10 minutes at 55-65°C.
2. Use two installments of TE Buffer to elute DNA from pinkRESIN. Each time, use approximately half the final volume.
3. For maximal yield combine the supernatant collected in step 1 with the remaining pinkRESIN™, suspend the pellet and continue with the protocol. Combine the final TE eluant with the original.

RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 8/8/2014 AB

This page is intentionally left blank



www.GBiosciences.com