



## *PinkRESIN*<sup>™</sup> Resin for DNA Cleaning

### INTRODUCTION

The *pinkRESIN*<sup>™</sup> removes excess salts, enzymes, unincorporated nucleotides, and primer-dimers from DNA preparations. DNA fragments larger than approximately 100 base pairs are isolated. Because *pinkRESIN*<sup>™</sup> has enhanced DNA 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.

### ITEM(S) SUPPLIED                      Cat # 786-89

<i>pinkRESIN</i> <sup>™</sup>	1ml x 1
-------------------------------	---------

### ITEMS NEEDED BUT NOT SUPPLIED

- DNA Binding Buffer
- Wash Buffer
- TE Buffer
- Wide bore pipet tips

### STORAGE CONDITIONS

It is shipped at ambient temperature. Store it at ambient temperature. This product is stable for 1 year.

### PROTOCOL

**1. Bind DNA to *pinkRESIN*<sup>™</sup>:** Add equal volume of a DNA Binding Buffer to each sample (plasmid DNA preparation, PCR reaction, etc.). Suspend *pinkRESIN*<sup>™</sup> by vigorous vortexing and add 10µl of the resin per 200µl of the sample using a wide bore pipet tip (if the sample contains more than 2µg DNA, add 5µl more *pinkRESIN*<sup>™</sup> for each µg DNA). Incubate on ice for 1-3 minutes. Centrifuge 10-30 seconds at high speed in a microcentrifuge. Remove the supernatant by decanting or pipetting (save supernatant if the sample is critical). Use the small remaining volume (20-30µl) in the tube to resuspend the *pinkRESIN*<sup>™</sup> pellet. Spin column users continue to step 2a.

**2. Wash:** Add 500µl ice cold DNA Wash and fully resuspend the *pinkRESIN*<sup>™</sup>. Centrifuge 10-30 seconds at high speed. Decant or pipet off the wash. Add 500µl DNA Wash and spin 10-30 seconds. Remove the second wash and let the pellet air dry for 1-2 minutes or until all traces of the wash are gone.

**3. Recover DNA:** Add pre-warmed (55-60°C) TE buffer or DNase-free water (*not supplied*) to the microcentrifuge tube and resuspend the *pinkRESIN*<sup>™</sup> pellet. Incubate at room temperature 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*<sup>™</sup>. Centrifuge 10-30 seconds to pellet the resin and transfer the sample containing cleaned DNA to a clean tube.

### Spin Column Protocol

**2a.** Add 300µl DNA Wash to the tube and fully resuspend the *pinkRESIN*<sup>™</sup>. Place a micro spin column on a capless 1.5 or 2 ml centrifuge tube. Pipet or pour the slurry into the spin column and centrifuge for 10-30 seconds at high speed. Add 300µl more DNA Wash to the tube and resuspend any remaining resin. Add this to the spin column and spin 10-30 seconds (there is no need to remove the first wash from the lower tube). Pour out the first two washes and place the micro spin column back in the capless tube. Add 600µl DNA Wash to the spin column and centrifuge. Drain the lower tube. A final spin is recommended to remove any wash from the side of the spin column.

**3a.** When all traces of DNA Wash are gone, place the spin column on a clean tube. Add pre-warmed (55-60°C) TE buffer or molecular grade water (see step 3 for recommended volumes) to the spin column and swirl to resuspend the resin. Place the spin column in a clean tube and centrifuge for 10-30 seconds to elute the DNA from *pinkRESIN*<sup>™</sup>.



### **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 (step 3) 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**

**I.NUCLEIC dotMETRIC™ (Cat # 786-61)**: Allows DNA, RNA and oligonucleotide concentrations to be measured using as little as 1µl of sample. Measurements take 2 minutes and are perfect for monitoring pinkCLEANUP™ yields or any other use requiring a minimal waste of sample.

**Note:** For other related products, visit our web site at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.