



186PR-01

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

A Geno Technology, Inc. (USA) brand name

GET™ CLEAN DNA

For Removal of Restriction Enzymes & PCR Clean Up

(Cat. # 786-356, 786-357)



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

INTRODUCTION

The GET™ CLEAN DNA kit uses spin columns 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. The GET™ Spin Columns have enhanced binding properties and therefore the protocols take as little as 5-10 minutes and yield high DNA recovery. Ideal applications include PCR clean up and restriction enzyme removal from plasmid DNA prior to *in-vitro* transcription.

ITEM(S) SUPPLIED

Description	Cat. # 786-356 50 Preps	Cat. # 786-357 100 Preps
Binding Buffer	30ml	2 x 30ml
DNA Wash	20ml	20ml
GET™ Spin Columns (Mini)	50	2 x 50
TE buffer	2 x 1.5ml	10ml

STORAGE CONDITIONS

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

PREPARATION BEFORE USE

1. Wash preparation: Add 80ml absolute ethanol to the DNA Wash. DNA WASH should be cooled on ice before use (alternatively, store wash at -20°C).
2. Preheat TE buffer to 55-60°C.

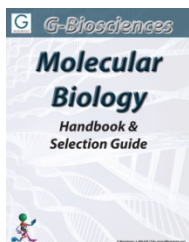
PROTOCOL

1. Add 5 volumes of Binding Buffer to 1 volume of the DNA sample to be cleaned and mix. It is not necessary to remove mineral oil or kerosene from PCR reactions. For example, add 500µl Binding Buffer to 100µl of PCR sample.
2. To bind the DNA, apply the sample to the spin column and centrifuge for 30-60 seconds at 14,000-16,000g.
3. Discard the flow-through. Place the column back into the same collection tube.
4. Add 750µl of DNA wash to the column and centrifuge for 30-60 seconds at 14,000-16,000g.
5. Discard flow-through and place column back in the collection tube. Centrifuge column at 14,000-16,000g for an additional one minute. This ensures that all DNA wash is removed.
6. Place column in a clean collection tube and elute the DNA by adding 50µl prewarmed TE buffer or water to the center of the membrane in the column and incubate at room temperature for 1 minute. Centrifuge for 1 minute at 14,000-16,000g.

NOTE: Alternatively, the sample can be eluted twice using half the recommended volume each time (2 x 25µl). For DNA fragments or plasmids >5kb increase TE buffer incubation to 5-10 minutes at 55-65°C.

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: 8/6/2012 CMH



www.GBiosciences.com