



263PR

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

A Geno Technology, Inc. (USA) brand name

Genomic Tube-O-DIALYZER™

For Purification of Genomic DNA

(Cat. # 786-142-45MC)



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

INTRODUCTION

Genomic Tube-O-DIALYZER™ allows dialysis of small genomic-DNA samples without taking the sample out of the tube and eliminates loss. A simple design converts a sample tube into a dialyzer– Genomic Tube-O-DIALYZER™ in which the tube cap is adapted with a dialysis membrane. Invert the Genomic-Tube-O-DIALYZER™ in a dialysis tank and dialyze your samples. Salts and other molecules rapidly exchange across the dialysis membrane mounted on the cap. When dialysis is complete, spin the tube for a brief 5 seconds and collect 100% of your sample. Sample collection by spinning allows clean and 100% recovery. There is no left over sample as in a dialysis bags.

After dialysis, replace the dialyzing cap with a Tube-O-Dialyzer™ Cap (Medi) supplied with each kit for storage of the sample. Genomic Tube-O-DIALYZER™ is supplied with 0.45µm size membrane, which allows rapid removal of impurities, RNA, small fragment DNA etc. and retains only genomic DNA.

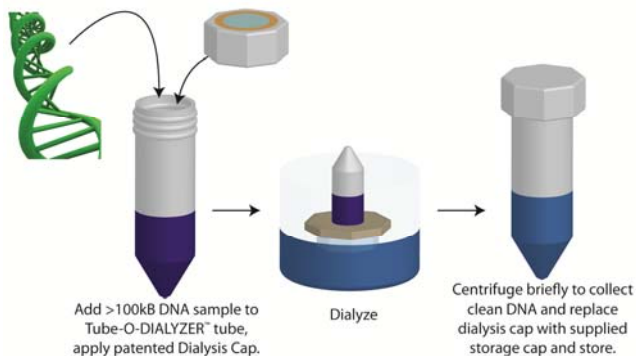
ITEM(S) SUPPLIED (CAT. # 786-142-45MC)

Description	Size
Genomic Tube-O-DIALYZER™ (0.45µm)	25
Tube-O-Dialyzer™ Caps (Medi)	25
Floats (Medi)	6
Forceps	1

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store at room temperature and is stable for 1 year.

PROTOCOL

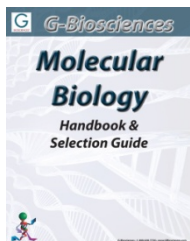


1. Immediately prior to use, remove the cap of the Genomic Tube-O-DIALYZER™. Be carefully not to disturb the membrane.

2. Place your sample in the Genomic Tube-O-DIALYZER™. Position the dialysis cap on the tube and tighten firmly.
3. Invert the Genomic Tube-O-DIALYZER™. Make sure the entire sample rests on the dialysis membrane. If sample is viscous, spin the Genomic Tube-O-DIALYZER™ in an inverted position (i.e. the dialysis membrane facing downward) for 5-6 seconds.
 - a. Place the Genomic Tube-O-DIALYZER™ in a 50ml centrifuge tube or Centrifuge Tube Adaptor (Cat. # 786-145) with the cap facing downward. Spin at 500-1000xg for 5-6 seconds.
 - b. Check the tube to make sure the entire sample has moved down on to the dialysis membrane of the Genomic Tube-O-DIALYZER™.
 - c. Hold the Genomic Tube-O-DIALYZER™ in an inverted position when removing from the centrifuge.
WARNING: *Spinning longer than necessary may rupture the membrane.*
4. Secure a float on the Tube-O-DIALYZER™ and float the assembly in an appropriate dialysis buffer.
5. Check that the dialysis membrane contacts the dialysis buffer. If there are large air-bubbles trapped underneath the dialysis membrane surface, tilt the tube or squirt buffer to remove the air bubbles. Gently, stir the dialysis buffer.
6. *Dialysis Time:* Dialysis time will depend on the nature of the sample, volume of sample as well as concentration of the sample solution. As a guide, the sample should be dialyzed 2-12 h. For efficient and complete dialysis we recommend inverting or gently tapping the Genomic Tube-O-DIALYZER™ 1-2 times during dialysis to mix the sample and replacing the dialysis buffer at least once.
7. After dialysis, remove the Genomic Tube-O-DIALYZER™ from the float and immediately spin the Genomic Tube-O-DIALYZER™ in an upright position for 5-6 seconds at 500-1000xg.
8. Discard the dialysis cap. Replace with a normal cap for storage. The genomic DNA sample is ready for use.

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.



www.GBiosciences.com