



028PR-02

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

A Geno Technology, Inc. (USA) brand name

Concentrator Powder

For Concentrating Solutions by Dialysis

(Cat. #786-144)



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

INTRODUCTION

Concentrator Powder is a high molecular weight polymer which will not migrate across the dialysis membrane. Simply transfer your dilute protein solution to a dialysis bag or dialysis device, such as our patented Tube-O-DIALYZER™ and then cover the membrane with Concentrator Powder. Concentrator Powder rapidly absorbs water from the sample and reduces the sample volume.

ITEM(S) SUPPLIED (Cat. # 786-144)

Description	Size
Concentrator Powder	150g

STORAGE CONDITIONS

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

ADDITIONAL ITEM(S) REQUIRED

- Dialysis Tubing or Device (*Tube-O-DIALYZER™*, Cat. # 786-610 to 786-624)

NOTE: Ensure that all dialysis membranes used have a MWCO <30kDa.

PROTOCOL

For Dialysis Tubing

1. Lay the dialysis tubing or bag containing the sample to be concentrated in a suitable tray.
2. Generously coat the dialysis membrane with the Concentrator Powder and incubate at 4°C to room temperature.
3. Visual inspect the dialysis bag during the incubation and peel off the Concentrator Powder once a suitable volume has been achieved.
4. The dialysis bag can be quickly rinsed in a suitable buffer to remove excess powder.

NOTE: The concentration can be accelerated by removing the dialysis bag from the Concentrator Powder every 30-60 minutes and then removing the wet concentrator from the dialysis membrane. A brief rinse with water or a suitable buffer may be performed to remove the Concentrator, afterward, remove all free solution from the membrane before returning the dialysis bag to the Concentrator Powder.

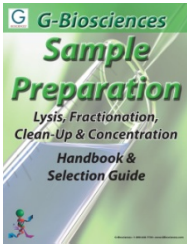
For Tube-O-DIALYZER™

1. Transfer 5-8gm Concentrator Powder into a suitable container.
2. Remove any excess solution from the outside of the TUBE-O-DIALYZER™.
3. Invert the Tube-O-DIALYZER™ into the Concentrator Powder, ensuring the cap of the Tube-O-DIALYZER™ is completely covered.
4. Incubate until desired volume and concentration is reached.
5. Quickly rinse in a suitable buffer to remove excess powder and then transfer your sample to a clean tube

NOTE: The concentration can be accelerated by removing the Tube-O-DIALYZER™ from the Concentrator Powder every 30-60 minutes and then removing the wet concentrator from the dialysis membrane. Rinse with water or a suitable buffer to remove the concentrator, afterward, remove all free solution from the membrane before inverting the Tube-O-DIALYZER™ back into the concentrator.

RELATED PRODUCTS

Download our Sample Preparation Handbook.



<http://info.gbiosciences.com/complete-protein-sample-preparation-handbook>

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

Last saved: 6/17/2013 CMH



www.GBiosciences.com