



449PR

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

A Geno Technology, Inc. (USA) brand name

Tube-O-Array™ Dialyzer

(Cat. #786-145A)



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

INTRODUCTION

Tube-O-Array™ Dialyzer is specifically developed for dialysis-equilibration of samples prior to 2D-gel analysis or other applications. Optimize up to 12 samples at a time. It consists of an Array Dialyzer Tray for holding 12 Tube-O-DIALYZER assemblies for simple and convenient equilibration-dialysis of the samples. It is suitable for 20µl to 2.5ml samples (Patent Pending).

ITEM(S) SUPPLIED (Cat. # 786-145A)

Description	Size
Array Dialyzer Tray	1
Micro-Dialysis Cup	12
Floats (Medi)	2 x 6
Floats (Micro)	12
Stirring Balls	100

NOTE- Depending on your application individual Tube-O-DIALYZER™ are available separately.

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store it at room temperature.

ADDITIONAL ITEM(S) REQUIRED

- **Tube-O-DIALYZER™** with dialysis membrane cut off MWCO 1,000, 4,000, 8,000, 15,000, and 50,000. For 20µl to 250µl samples use Micro and for 200µl to 2.5ml samples use Medi size Tube-O-DIALYZER™ .
- **Centrifuge Tube Adaptor** [Cat # 786-145] Designed to assist centrifugation of Tube-O-DIALYZER™ in a bench top centrifuge. Centrifuge Tube Adaptor can be used for both Micro and Medi size Tube-O-DIALYZER™ .

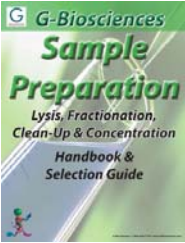
PROTOCOL

1. Select an appropriate MWCO of the Tube-O-DIALYZER™ . For reducing conductivity or de-salting, use 4kd Tube-O-DIALYZER™ . Follow the instructions supplied with the Tube-O-DIALYZER™ . Place your sample in the Tube-O-DIALYZER™ (for 20µl to 250µl samples use Micro and for 200µl to 2.5ml samples use Medi size Tube-O-DIALYZER™). Position the dialysis cap on the tube and tighten it firmly.
2. Secure a float and position the Tube-O-DIALYZER™ into the micro dialysis cups containing at least 2ml dialysis or equilibration buffer and 5-6 stirring balls.
3. Position the dialysis assembly into the Array DIALYZER™ tray. Place the Array DIALYZER™ tray on a shaker plate. Please note that the shaker must rotate at least 100-150 rotations per minute.

Dialysis Time: Dialysis time will depend on the nature of sample, MWCO of Tube-O-DIALYZER™, and the volume of the sample and the dialysis buffer. As a guide, the sample should be dialyzed 1-3 h (longer dialysis time may be necessary in some applications). For faster equilibration and better results, dialysis buffer should also be replaced at least once during dialysis.

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: 5/30/2014 CMH



www.GBiosciences.com