



077PR

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

A Geno Technology, Inc. (USA) brand name

Centrifuge Tube Adaptor

For Tube-O-Dialyzer™

(Cat. #786-145)



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

This page is intentionally left blank

INTRODUCTION

Our Centrifuge Tube Adaptor for Tube-O-Dialyzer™ is designed to assist centrifugation of Tube-O-Dialyzer™ using a bench top centrifuge. The adaptor can be used for both Micro and Medi size Tube-O-Dialyzer™. The Tube-O-Dialyzer™ can be positioned into the adaptor either upright position or inverted position for centrifugation.

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

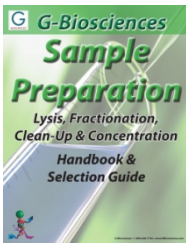
Description	Size
Centrifuge Tube Adaptor for Tube-O-Dialyzer™	Qty: 2

INSTRUCTION FOR USE

1. Use a bench top centrifuge fitted with a 50 ml bucket. Use two adaptors to balance the centrifuge bucket.
2. Use either the narrow or wide bore side of the Tube-O-Dialyzer™ Adaptor that is appropriate for the Tube-O-Dialyzer™ in use. Place the adaptors into the centrifuge buckets.
3. Position the Tube-O-Dialyzer™ in the adaptor, either in upright or inverted position depending on the nature of use.
4. Centrifuge at 500 – 1,000xg for 5-10 seconds.
5. Remove the Tube-O-Dialyzer™ from the adaptor at the end of centrifugation.

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: 8/20/2013 CMH



www.GBiosciences.com