



## Centrifuge Tube Adaptor For Tube-O-Dialyzer™

### 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 up right position or inverted position for centrifugation.

### ITEM(S) SUPPLIED

Cat# 786-145

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

1. **NI™-Protein Assay (Cat # 786-005):** A protein assay that is free from interference of common laboratory agents, including reducing agents, detergents, dyes, EDTA, etc.
2. **Protein Standards:** We supply Bovine Gamma Globulin (2mg/ml) [Cat#786-007] and Bovine Serum Albumin (2mg/ml) [Cat#786-006] as protein standards.
3. **Pre-Diluted Protein Standard (Cat# 786-114):** BSA protein standard in an easy-to-use pre-diluted format, supplied as 6x5ml standards ranging from 0.1-1mg/ml.
4. **Cuvettes (Cat# 786-009):** For users convenience, we also supply semi-micro cuvettes (1.5ml) suitable for use in NI-Protein Assay. Call us for more information and price.
5. **Assay Tubes (Cat# 786-008):** For users convenience, we also supply assay tubes (2ml). We have extensively tested these assay tubes to work with NI Protein Assay, particularly in precipitation steps using UPPA.

**NOTE:** For other related products, visit our web site at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Rev 05.21.09/MM

