



## Proteomic Protein Control

### INTRODUCTION

Proteomic protein controls are specifically prepared as controls for standardizing 2D electrophoresis methods and protocols. Protein preparations are substantially free from non-protein agents such as nucleic acids, detergents, salts, lipids, natural products, common laboratory agents and has low conductivity. Proteomic protein controls are supplied as dry protein pellets (2 x 2mg/vial each) in ready-to-use form. Simply re-hydrate and dissolve protein pellets in an appropriate buffer of your choice and its ready to use. The *Proteomic Protein Controls* are supplied individually and as a set, which contains one vial (2.0 mg protein) each of Animal cells, *E. coli*, Yeast, and Plant proteins.

### ITEM(S) SUPPLIED

Cat #	Description	Size
PSC-01	<i>Proteomic Protein Control-Yeast</i>	2 x 2mg Protein & 2 Pestles
PSC-02	<i>Proteomic Protein Control- Animal Cells</i>	2 x 2mg Protein & 2 Pestles
PSC-03	<i>Proteomic Protein Control- E. coli</i>	2 x 2mg Protein & 2 Pestles
PSC-04	<i>Proteomic Protein Control- Plant</i>	2 x 2mg Protein & 2 Pestles
PSC-05	<i>Proteomic Protein Control Set of Yeast, Animal cell, E. coli and Plant proteins</i>	4 x 2mg Protein (1 from each species) & 4 Pestles

### STORAGE CONDITIONS

It is shipped at ambient temperature. Store the supplied (un-hydrated) protein vials at -20°C. After hydration, store at -70°C. (**Note:** Avoid frequent freezing and thawing). When stored and handled properly, stable for 12 months.

### Protein Hydration & Solubilization

Centrifuge the supplied tube for 30 seconds to collect the contents in the bottom of the tube.

To solubilize the supplied protein, first add 100µl solubilization buffer of your choice containing detergent(s) and or chaotropic agent(s) into the provided protein pellet tube. The solubilization buffer must also contain a reducing agent such as 2-10mM TCEP (Cat. #786-230) or DTT up to 100mM for achieving highest solubility. Use the provided pestle to dis-aggregate and mix the protein pellet. (**Note:** *effective grinding & mixing of pellet shortens the protein solubilization time*)

Rinse the pestle with the solubilization buffer into the tube and adjust the final volume with the same buffer to give 2mg/ml protein concentration or as needed. Incubate the tube at RT and periodically vortex it until the protein solution is clear.

### RELATED PRODUCTS

**1. Focus Protease Arrest (Cat # 786-108F):** A protease cocktail specifically developed for sample preparation for 2D-studies and provides 95-98% inhibition of protease activity.

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

Rev 05.29.09/MM

