



457PR-02

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# UPPA-PROTEIN Concentrate™

For Concentrating Dilute Protein Solution

(Cat. # 786-120, 786-121)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

INTRODUCTION ..... 3

APPLICATIONS ..... 3

ITEM(S) SUPPLIED ..... 3

STORAGE CONDITIONS ..... 3

ADDITIONAL ITEMS REQUIRED ..... 3

PROTOCOLS ..... 4

PROCESSING LARGE SAMPLES ..... 5

RELATED PRODUCTS ..... 6

## INTRODUCTION

PROTEIN-Concentrate™ kit uses a proprietary reagent, Universal Protein Precipitation Agent (UPPA; Patent Pending). Protein solutions as dilute as 1ng/ml can be quantitatively concentrated into a small volume. Protein precipitation is not affected by the presence of detergents, chaotropics, or other common laboratory agents. After precipitation, the precipitate is washed to remove salts and other agents which produces protein samples of conductivity ~ 40-60µS - ideal for critical analysis. The protein precipitate is suspended in a small volume buffer. If the protocol is followed correctly, the recovery is generally 100%.

The PROTEIN-Concentrate™ kit is supplied in two sizes: the Micro Kit is for concentrating up to a total of 10 ml dilute protein solution, either single or multiple procedures, and the Medi Kit is for concentration up to a total of 30 ml dilute protein solution, either single or multiple procedures. Additional volumes of any reagent may be purchased separately.

## APPLICATIONS

Suitable for concentrating proteins for running gels, raising antibodies, protein purification, protein assays, and other applications. This kit is not suitable for those proteins which may lose some of their biological activities when precipitated, for such proteins use either Column-PROTEIN Concentration™ kit (Cat. #786-126) or OrgoSol PROTEIN Concentration™ kit (Cat. # 786-125).

## ITEM(S) SUPPLIED

Cat. #	786-120 (Micro)	786-121 (Medi)
UPPA™- I	30ml	100ml
UPPA™- II	30ml	100ml
OrgoSol Buffer™	50ml	2 x 50ml
UPC-Wash™	2.0ml	2 x 2.0ml
SEED™	300µl	2 x 300µl
Solubilization Buffer-I	2.0ml	2 x 2.0ml
Solubilization Buffer-II	0.5ml	2 x 0.5ml

## STORAGE CONDITIONS

The kit is shipped at ambient temperature. Store all the components at room temperature upon arrival.

**NOTE:** Chill OrgoSol Buffer at -20°C for ~1hr or more before use- see step 10 of the protocol.

## ADDITIONAL ITEMS REQUIRED

Centrifuge, Centrifuge Tubes, Microfuge, & Spin Columns

## PROTOCOLS

**NOTE:** Perform the entire procedure in the cold (ice bucket) unless specified otherwise. Concentration should be performed in a centrifuge tube. For small volumes, use microfuge tubes. Always position microfuge-tubes in the centrifuge at the same orientation, i.e. cap-hinge facing out-ward. This will allow the pellet to remain glued to the same side of the tube during repeated centrifugations and minimize the loss of protein pellets.

1. Mix 1 volume of protein solution with 3 volumes of UPPA-I (See example below). Vortex the mixture and incubate at 4°C for 10-15 minutes.
2. Add 3 volumes of UPPA-II in to the mixture of protein and UPPA-I (See example below). Vortex and centrifuge the tube.

**Example:** For 0.1ml protein solution, add 0.3ml UPPA-I, incubate and then add 0.3ml UPPA-II. Also, review Processing large Samples Protocol below.

3. Centrifuge the tube at 15,000xg for 5 minutes to form a tight pellet.
4. As soon as the centrifuge stops, remove the tube from the centrifuge.

**NOTE:** Pellets should not be allowed to diffuse after centrifugation is complete.
5. Carefully and without disturbing the pellet, use a pipette tip and remove the entire supernatant.
6. Carefully re-position the tube in the centrifuge as before, i.e. cap-hinge facing out-ward. Centrifuge the tube again for 30 seconds. Use a pipette tip and remove & discard the remaining supernatant.
7. Add 40µl of UPC-Wash on top of the pellet (for larger sample size, add Wash 3-4 x times the size of the pellet).
8. Carefully re-position the tube in the centrifuge as before, i.e. cap-hinge facing outward. Centrifuge the tube again for 5 minutes. Use a pipette tip, remove and discard the Wash
9. Add 25µl of pure water on top of the pellet (i.e., add water just enough to cover the pellet - a volume equal to the size of the pellet). Vortex the tube.

**NOTE:** Pellets do not dissolve in water.

10. For each 0.1-0.3ml protein solution, add 1ml OrgoSol Buffer (pre-chilled at -20°C), and 5µl SEED.

**NOTE:** In addition, OrgoSol Buffer must be at least 10 fold in excess of the water added in Step 8.
11. Vortex to suspend the pellet. It is important that the pellet is fully suspended in OrgoSol Buffer.

**NOTE:** Pellets do not dissolve in OrgoSol Buffer.
12. Incubate the tube at -20°C for 30 minutes. Periodically vortex the tube, 20-30 seconds vortex each burst.

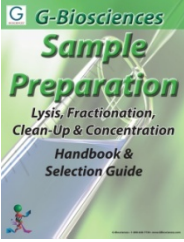
13. Centrifuge at 15,000xg for 5 minutes to form a tight pellet.
14. Remove and discard the supernatant. You will notice a white pellet in the tube. Air dry the pellet. On drying, the white pellet will turn translucent.  
***NOTE-** Do not over dry the pellets as these pellets may be difficult to dissolve.*
15. Suspend the pellet in an appropriate volume of Solubilization Buffer-I (5-50  $\mu$ l Solubilization Buffer-I). Vortex to suspend the pellet. Incubate for 2 minutes.
16. Add Solubilization Buffer-II. For each 5 $\mu$ l Solubilization Buffer-I used, add 1 $\mu$ l of Solubilization Buffer-II. Incubate for 5 minutes. After the pellet is dissolved, centrifuge and collect a clear protein solution. The protein solution at this stage contains 60mM Tris, pH 7-7.5. After dissolving the pellet, the protein solution may be mixed with SDS, Urea, Guanidine.HCl, SDS-PAGE gel loading buffer or other types of buffers and agents.
17. For buffer exchange, the protein suspension may be dialyzed or passed through a pre-equilibrated spin column.

### **PROCESSING LARGE SAMPLES**

Samples containing > 100 $\mu$ g protein produces large and tightly packed protein pellets which require a longer time to dissolve in Buffers. Grinding of the protein pellet with a pestle will accelerate solubilization of the pellet. We recommend use of microfuge tubes and tight fitting pestle for processing samples containing larger than 100 $\mu$ g protein. See related products for ordering information.

## 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](http://www.GBiosciences.com) or contact us.

Last saved: 5/19/2015 CMH

*This page is intentionally left blank*



[www.GBiosciences.com](http://www.GBiosciences.com)