





A Geno Technology, Inc. (USA) brand name

# AlbuminOUT™

(Cat. # 786-251, 786-252, 786-251T)



INTRODUCTION	3
ITEM(S) INCLUDED	3
STORAGE CONDITION	3
ADDITIONAL ITEMS REQUIRED	3
PROTOCOL	4
ELUTION OF COLUMN BOND ALBUMIN AND OTHER PROTEINS (OPTIONAL)	5
REMOVAL OF ALBUMIN FROM OTHER SPECIES:	5
RELATED PRODUCTS	5

#### INTRODUCTION

Samples that contain a large abundance of albumin, such as plasma and cerebrospinal fluid, tend to mask identification and discovery of other less abundant proteins in 2D gel electrophoresis. AlbuminOUT™ has been specifically developed for substantial removal of albumin from such samples. The albumin removal method is based on binding of albumin with Cibachron Blue dye. AlbuminOUT™ has been optimized for removal of human albumin from samples. AlbuminOUT™ is rapid spin column method, each column contains (0.2ml) dye bond resins with capacity > 2mg human albumin per column. AlbuminOUT™ will remove over >98% albumin from 5-50µl human plasma. Other nucleotide binding proteins may also be captured and removed by Cibachron Blue dye resins. AlbuminOUT™ may also be used for removal of albumin from other species. AlbuminOUT™ is suitable for processing 25 (Cat # 786-251) and 50 (Cat# 786-252) samples respectively.

## ITEM(S) INCLUDED

Description	Cat. # 786-251	Cat. # 786-252	Cat. # 786-251T
Albumin Binding Buffer	25ml	2x25ml	5ml
Albumin Elution Buffer	6ml	2x 6ml	2ml
AlbuminOUT <sup>™</sup> Spin Column	25	50	4
Collection Tubes (2ml)	50	100	10

## STORAGE CONDITION

The kit is shipped at ambient temperature. Upon arrival, store it at 4°C. When stored and used as recommended, the kit is good for use for 12 months.

## ADDITIONAL ITEMS REQUIRED

- Micro centrifuge
- 1.5ml collection tubes.

#### **PROTOCOL**

Please note that albumin content in samples may vary. Typically, each column will bind and remove >98% albumin (~2mg albumin) from 5-50 $\mu$ l human serum. Other factors, such as salt concentration, may influence the albumin retention capacity of the column, therefore, the total human serum load on the column must not exceed 50 $\mu$ l (undiluted human serum) - whenever possible, keep the human serum load under 50 $\mu$ l.

- Transfer 50μl Albumin Binding Buffer into a 1.5ml tube. Add 5-50μl sample containing albumin (serum or cerebrospinal fluid) into the tube. Mix the content and centrifuge 10,000xg for 5 minutes in a cold centrifuge. Collect the clear supernatant and store in an ice-bucket until used.
- Spin the AlbuminOUT<sup>™</sup> Spin Column for a brief 5 seconds. Break off and open the bottom plug of the column. Reposition the column in a 2ml Collection Tube provided with this kit.
- 3. Add 100µl Albumin Binding Buffer and spin the AlbuminOUT<sup>™</sup> Spin Column at 1000xg for 10-15 seconds. Repeat this step two more times. Empty the collection tube and replace the column in the collection tube.
- Load the diluted sample into the AlbuminOUT<sup>™</sup> Spin Column. Incubate the column for 1-2 minutes at room temperature.
- Collect any flow-through and re-apply to the AlbuminOUT<sup>™</sup> Spin Column. Incubate the column for 1-2 minutes at room temperature.
- 6. Centrifuge the AlbuminOUT<sup>™</sup> Spin Column at 1000xg for 5 seconds. Collect the albumin free flow-through for further processing.

**NOTE:** The flow through sample contains a salt concentration may not be suitable for 2D gel analysis. Dialyze the albumin free sample using 1-4Kd molecular weight cut off Tube-O-Dialyzer (Cat. # 786-141-143) against a salt free buffer. Alternatively, clean the sample with Perfect-FOCUS (Cat #786-124) before 2D gel analysis.

## Elution Of Column Bond Albumin And Other Proteins (Optional)

- 1. Wash the column 3 times with Albumin Binding Buffer, 200µl each wash.
- 2. Add 200µl Albumin Binding Buffer and spin the column at 1000xg for 5 seconds. Repeat this step twice. Empty the collection tube and replace the column in the collection tube.
- 3. Add 100-200µl Albumin Elution Buffer. Incubate the column for 5 minutes at room temperature. Centrifuge the spin column at 1000xg for 5 seconds.
- 4. Collect the flow-through. The flow-through predominantly contains albumin. Contamination with other proteins cannot be ruled-out. Other nucleotide binding proteins may also be captured by Cibachron Blue dye resins (AlbuminOUT Spin Column) and co-eluted with albumin.

**NOTE:**The eluted albumin contains a salt concentration may not be suitable for 2D gel analysis. Dialyze the eluent albumin using 1-4Kd molecular weight cut off Tube-O-Dialyzer (Cat.# 786-141-143) against a salt free buffer. Alternatively, clean the sample with Perfect-FOCUS (Cat#786-124) before 2D gel analysis.

#### REMOVAL OF ALBUMIN FROM OTHER SPECIES:

AlbuminOUT<sup>™</sup> may be used for removal of albumin from other sources such as pig, dog, sheep, rabbit, rat, and bovine. Binding efficiency of non-human source may vary between 1-2 mg albumin/column; therefore, it is important to perform a test before use. The protocol, as outlined above, is suitable for removal of albumin from any species or sample. In most cases, reducing the undiluted sample load by 30%-50% would be sufficient.

#### 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.

This page is intentionally left blank

This page is intentionally left blank



www.GBiosciences.com