I. Introduction

This protocol is provided for **Capturem Protein G 96-Well Plate** (Cat. No. 635726), a single-use disposable plate for simple, rapid purification of antibodies from animal sera, ascites fluid, cell culture media, and other sources. Each well can hold up to 1 ml of sample and requires a minimum elution volume of 200 µl. Depending upon the sample type, species, and antibody isotype the binding capacity varies. More information about bed volume and capacities can be found on <u>our website</u>.

II. Materials and Reagents

A. Components

• 1 Capturem Protein G 96-Well Plate (Cat. No. 635726)

B. Additional Materials Required

1. Purification Buffers

This kit is compatible with all standard buffers typically used for antibody purification, such as phosphate- and acetate-based buffers. We highly **recommend using a binding buffer with an optimal pH of 5,** preferably Protein G IgG binding buffer from Thermo Fisher Scientific, Cat. No. 21019. Loading the sample (e.g. serum matrix) without prior dilution is not recommended.

- Equilibration/Binding/Wash Buffer: To obtain optimal and reproducible performance, we recommend using Pierce Protein G IgG binding buffer. You can also use 20–100 mM sodium acetate containing 0.15–2 M NaCl, pH 5.0, or 100–500 mM sodium phosphate containing 0.15–2 M NaCl, pH 7.0.
- Wash Buffer 2: Dulbecco's PBS Buffer w/o Ca²⁺, Mg²⁺ (pH 7.5)
- Elution Buffer: 0.1 M glycine, pH 2.5–3.0
- Neutralization Buffer: 1 M Tris, pH 8.5

2. Collection Plates

- 96-well collection plates: Each purification will require four standard 96-well collection
 plates. These plates should be used throughout the protocol to collect flowthrough and
 wash samples that will be saved for SDS-PAGE analysis and/or colorimetric protein
 assays (e.g., Bradford assays). Alternatively, if only the eluates are to be analyzed, the
 flowthrough from the equilibration and wash steps can be discarded. In this case, only
 two collection plates are required.
- **96-well plate filtration device:** The Capturem Protein G 96-Well Plate is compatible with standard 96-well plate filtration devices, such as the NucleoVac 96 Vacuum Manifold (Cat. No. 740681).
- **96-well plate centrifuge:** The Capturem Protein G 96-well Plate is compatible with standard 96-well plate centrifuges (e.g. Eppendorf Centrifuge 5804 R with Deepwellplate rotor, A-2-DWP).
- Multi-channel pipette

III. Sample Preparation

- 1. Follow the standard protocols for preparing antibody samples. We recommend diluting the antibody sample with binding buffer in the range of 1:2–1:15 sample to binding buffer. A 1:2 1:4 dilution is recommended for hybridoma samples. We suggest a dilution of 1:15 for serum samples, which may differ depending on the species and may require further optimization to obtain high yields and purity. Loading the sample (e.g. serum matrix) without prior dilution is not recommended.
- 2. Samples with visible precipitates must be clarified by centrifugation or filtered through a 0.8-micron filter before loading onto the plate.

IV. Antibody Purification

Purification of antibodies from the samples prepared in Section III may be performed using vacuum filtration (Protocol A) or centrifugation in a standard 96-well plate centrifuge (Protocol B).

A. Protocol: Antibody Purification Using Vacuum Filtration

- 1. Assemble the Capturem Protein G 96-Well Plate in a 96-well plate filtration device together with a 96-well collection plate (supplied by the user—see Section II.B).
- 2. Add 800 μl of Equilibration/Binding/Wash Buffer to each well of the Capturem Protein G 96-Well Plate using a multi-channel pipette, to equilibrate the wells. We recommend Protein G IgG binding buffer from Thermo Fisher Scientific. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Remove the flowthrough and discard it along with the collection plate. Reassemble the Capturem Protein G 96-Well Plate in the 96-well plate filtration device together with a new collection plate.

NOTE: The required pressure may vary between 0.2–0.8 bar, depending on the sample. When performing vacuum filtration, use the following conversion factors to convert between different vacuum pressure units:

Pressure conversion factors:

1 mbar = $100 \text{ Pa} = 0.750 \text{ mm Hg} = 14.504 \text{ x } 10^{-3} \text{ psi} = 0.987 \text{ x } 10^{-3} \text{ atm}$

- 3. Load 200–1,000 µl of diluted antibody (from Section III) into each well of the equilibrated Capturem Protein G 96-Well Plate. We recommend Protein G IgG binding buffer from Thermo Fisher Scientific. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis. Reassemble the Capturem Protein G 96-Well Plate in the 96-well plate filtration device together with a new collection plate.
- 4. Add 300–800 µl of Equilibration/Binding/Wash Buffer to each well of the Capturem Protein G 96-Well Plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis.
- 5. Wash a second time by adding 400–800 μl of **Wash Buffer 2** (PBS Buffer w/o Ca²⁺, Mg²⁺, pH 7.5) to each well of the Capturem Protein G 96-Well Plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis.

NOTE: If significant background due to non-specific binding of other proteins is observed, please, include after step 5 an additional wash step by adding 400–800 µl Wash Buffer 2 (PBS w/o Ca, Mg). Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis.

- 6. Elute your antibody from the Capturem Protein G 96-Well Plate as follows:
 - a. Add 5–30 µl of Neutralization Buffer (1/10 the volume of Elution Buffer to be used) to each well of a new collection plate and reassemble the collection plate together with the Capturem plate in the 96-well plate filtration device for the first elution.
 - b. Add $50-300 \,\mu$ l of Elution Buffer to each well of the Capturem Protein G 96-Well Plate. Use a vacuum ($\sim 0.3-0.4$ bar) to run the buffer through the Capturem plate wells into the collection plate.
 - c. Perform a second elution by repeating Steps a and b above using a new collection plate. The wells of the two collection plates should contain your eluted antibodies, which are now ready for analysis.
- 7. Measure the amount of antibody in both eluates using UV absorbance at 280 nm or another colorimetric protein analysis method. One OD₂₈₀ unit typically corresponds to 0.73 mg/ml of IgG. The purity of the eluted antibodies can be analyzed by SDS-PAGE, size exclusion chromatography, etc.

NOTE: Most of the antibody is recovered in the two elution steps.

B. Protocol: Antibody Purification Using Centrifugation

- 1. Place the Capturem Protein G 96-Well Plate securely on top of a 96-well collection plate (supplied by the user—see Section II.B).
- 2. Add 800 μl of Equilibration/Binding/Wash Buffer, preferably Protein G IgG binding buffer from Thermo Fisher Scientific, to each well of the Capturem Protein G 96-Well Plate using a multi-channel pipette, to equilibrate the wells. Centrifuge at 2,000g for 2 min at room temperature. Remove the flowthrough and discard it along with the collection plate. Place the Capturem Protein G 96-Well Plate securely on top of a new 96-well collection plate.

NOTE: When performing centrifugation, use the following formula to convert rpm to centrifugal force units (RCF or g) if the centrifuge does not automatically provide this information:

Centrifugal force conversion formula:

RCF or $g = (1.12) \times (R) \times (rpm/1,000)^2$

Where R is the distance (in millimeters) from the center of the rotor to the end of the spin bucket when held horizontally and away from the center.

If a well does not drain entirely, centrifuge again at 2,000g for 2 min.

- 3. Load 200–1,000 µl of diluted antibody (from Section III) into each well of the equilibrated Capturem Protein G 96-Well Plate. Centrifuge at 2,000g for 2 min at room temperature. Save the collection plate containing the flowthrough for antibody analysis. Place the Capturem Protein G 96-Well Plate securely on top of a new 96-well collection plate.
- 4. Add 300–800 μl of Equilibration/Binding/Wash Buffer to each well of the Capturem Protein G 96-Well Plate. Centrifuge at 2,000g for 2 min at room temperature. Save the collection plate containing the flowthrough for antibody analysis.
- 5. Wash a second time by adding 400–800 μl of **Wash Buffer 2** (PBS Buffer w/o Ca²⁺, Mg²⁺, pH 7.5) to each well of the Capturem Protein G 96-Well Plate. Centrifuge at 2,000g for 2 min at room temperature. Save the collection plate containing the flowthrough for antibody analysis.

NOTE: If significant background due to non-specific binding of other proteins is observed, please, include after step 5 an additional wash step by adding 400–800 µl Wash Buffer 2 (PBS w/o Ca, Mg). Centrifuge at 2,000g for 2 min. at room temperature to drain the wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis.

- 6. Elute your antibody from the Capturem Protein G 96-Well Plate as follows:
 - a. Add 10– $30~\mu l$ of Neutralization Buffer (1/10 the volume of Elution Buffer to be used) to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
 - b. Add 100–300 µl of Elution Buffer to each well of the Capturem Protein G 96-Well Plate.
 Centrifuge at 2,000g for 2 min at room temperature.
 - c. Perform a second elution by repeating Steps a and b above using a new collection plate. The wells of the two collection plates should contain your eluted antibodies, which are now ready for analysis.

7. Measure the amount of antibody in both eluates using UV absorbance at 280 nm or another colorimetric protein analysis method. One OD₂₈₀ unit typically corresponds to 0.73 mg/ml of IgG. The purity of the eluted antibodies can be analyzed by SDS-PAGE, size exclusion chromatography, etc.

NOTE: Most of the antibody is recovered in the two elution steps.

Appendix A. Troubleshooting Guide

Table 1. Troubleshooting Guide

Problem	Possible Explanation	Solution
Background bands/poor purity	Non-specific binding of proteins to membrane	Add an additional wash step with Wash Buffer 2 (PBS) or TBS
Minimal recovery	The sample contains more antibody than the Protein G 96-Well Plate has capacity for.	Reduce the amount of sample added to each well. If you need to purify more antibody, consider using Capturem Protein G 24 well or Capturem Protein G Maxi kit which has a higher capacity (1-2mg).
Low yield of purified antibody	Poor binding affinity of the IgG subtype used. Different subtype classes and species bind differently to Protein G. Please, refer to <i>J. Chromatogr A.</i> 2007 ,1160, 44-55.	Verify that Protein G based purification is suitable for the IgG subtype class and species used. Alternatively, consider using Protein A based purification instead.
Low yield of purified antibody	pH of the binding buffer was altered by the hybridoma, serum or cell culture medium used.	Make sure the pH of the buffer is in optimal range (pH 5). We recommend Protein G IgG buffer from ThermoFisher Scientific as binding buffer.
Low yield of purified antibody	Suboptimal dilution	Make sure to use the optimal dilution of your sample. For hybridoma medium 1:2 to 1:4 were found to work well, but it is sample dependent.
Antibody does not elute	Elution conditions too mild	 Follow the instructions use recommended elution buffer (0.1 M glycine, pH 2.5–3.0) Make sure the pH of the elution buffer is 2.5-3.0 if you are using different buffer than the one recommended.
Spin column does not fully drain	Clogging due to particles or very viscous sample	 Pre-clarify the solution by centrifugation or using a 0.8-micron filter Repeat spin at 2,000g

Contact Us		
Customer Service/Ordering	Technical Support	
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)	
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)	
web: takarabio.com	web: takarabio.com	
e-mail: ordersUS@takarabio.com	e-mail: techUS@takarabio.com	

Notice to Purchaser

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Takara Bio USA, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <u>takarabio.com</u>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

© 2017 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at <u>takarabio.com</u>.

This document has been reviewed and approved by the Quality Department.