

Version: 02

Update: 09/30/2022

## Protein A Resin FF Prepacked Column

### Cat. No. L00680

|     |                              |   |
|-----|------------------------------|---|
| I   | Product Description .....    | 1 |
| II  | Buffer Preparation .....     | 2 |
| III | Purification Procedure ..... | 2 |
| IV  | Storage .....                | 2 |
| V   | Troubleshooting .....        | 3 |
| VI  | Ordering Information.....    | 3 |

### I Product Description

GenScript Protein A Resin FF is an affinity chromatography media designed for easy, one-step purification of classes, subclasses and fragments of immunoglobulins from biological fluids and from cell culture media. Protein A Resin FF can also be used for immunoprecipitation of proteins, protein complexes or antigens. The recombinant protein A ligand is coupled to 4% highly cross-linked agarose. The coupling is optimized to have a high binding capacity for immunoglobulins. The static binding capacity of Protein A Resin FF is greater than 40 mg human IgG/ ml settled resin. The dynamic binding capacity will vary depending on several factors such as target antibody, flow rate, etc. Table 1 lists the characteristics of Protein A Resin FF Prepacked Column.

Protein A, a bacterial cell wall protein isolated from *Staphylococcus aureus*, binds to mammalian IgGs mainly at their Fc region. Native protein A has five IgG binding domains and many other domains with unknown functions. Recombinant protein A contains five high affinity IgG binding domains with other non-essential domains removed to reduce nonspecific binding. Since only the Fc region is involved in binding to recombinant protein A, the Fab region is available for binding antigens.

**Table 1. Characteristics of Protein A Resin FF Prepacked Column**

|                                        |                                                                        |
|----------------------------------------|------------------------------------------------------------------------|
| Resin Volume                           | 1 ml x 2(L00680-12);5ml x 1(L00680-51);5ml x 5(L00680-55)              |
| Ligand                                 | Recombinant <i>Streptococcal</i> protein A expressed in <i>E. coli</i> |
| Number of IgG binding sites per ligand | 5                                                                      |
| M.W. of ligand                         | Approximately 34 kDa                                                   |
| PI of ligand                           | 5.17                                                                   |
| Degree of substitution                 | Approximately 5 mg protein A / ml settled resin                        |
| Static binding capacity                | > 40 mg human IgG/ ml settled resin                                    |
| Matrix component                       | Agarose, 4% highly cross-linked                                        |
| Average particle size                  | 90 µm (45-165 µm)                                                      |
| Storage solution                       | 20% ethanol                                                            |
| Storage conditions                     | 2-8 °C                                                                 |

### II Buffer Preparation

Water and chemicals used for buffer preparation should be of the highest purity. It is recommended to filter the buffers by passing them through a 0.45 µm filter before use.

Binding/Wash Buffer: 0.15 M NaCl, 20 mM Na<sub>2</sub>HPO<sub>4</sub>, pH 7.0

Elution Buffer: 0.1 M glycine, pH 3.0

Neutralization Buffer: 1 M Tris-HCl, pH 8.5

### III Purification Procedure

This procedure is optimized for a column of 1 ml bed volume. The volumes of the reagents can be scaled up or down according to the size of the column.

- **Sample Preparation**

To insure that proper ionic strength and pH are maintained for optimal binding, it is necessary to dilute serum samples, ascites fluid or cell culture supernatant at least 1:1 with Binding/Wash Buffer. Alternatively, the sample may be dialyzed overnight against Binding/Wash Buffer.

- **Column Purification**

1. Add the sample into the column and drain the flow-through with a flow speed of about 1 ml/min. Collect the flow-through to measure the binding efficiency to the resin, i.e. by SDS-PAGE.
2. Wash the column with 30 ml Binding/Wash Buffer and drain the buffer with a flow speed of about 2 ml/min, or until the absorbance of the effluent at 280 nm is stable.
3. Elute the antibody with 10-15 ml Elution Buffer and drain the eluate with a flow speed of about 1 ml/min. Collect the eluate containing the target immunoglobulin and immediately neutralize to pH 7.4 with Neutralization Buffer (1/10 volume of total eluate).

- **Regeneration of Column**

Regenerate the column by washing the resin with 10 ml Elution Buffer followed by equilibration with 5 ml Binding/Wash buffer. Columns can be regenerated up to 10 times without significant loss of binding capacity.

### IV Storage

Store regenerated Protein A Resin FF Prepacked Column in Binding/Wash Buffer containing 20% ethanol at 2°C to 8°C. **Do not freeze.**

### V Trouble Shooting

| <b>Problem</b>                                                                                     | <b>Possible Cause</b>                                                             | <b>Solution</b>                                                                                                            |
|----------------------------------------------------------------------------------------------------|-----------------------------------------------------------------------------------|----------------------------------------------------------------------------------------------------------------------------|
| The flow rate of the column is very low (<0.5 ml/minute).                                          | Tiny air bubbles from buffer or particles from sample are blocking the gel pores. | De-gas buffers and samples. Do not allow the column to dry.                                                                |
| A considerable amount of sample has been loaded, but no specific antibody of interest is detected. | The concentration of the antibody of interest is very low.                        | Purify the antibody using the specific antigen coupled to a resin (i.e., High-Affinity Iodoacetyl Resin, Cat. No. L00403). |
| The antibody is degraded.                                                                          | The antibody is sensitive to low-pH elution buffer                                | Neutralize the eluted fractions with Neutralization Buffer immediately after elution.                                      |
| No antibody is detected in any elution fraction.                                                   | The IgG subclass does not bind to protein A.                                      | Try other affinity chromatography media to purify the antibody, such as Protein G Resin or Protein L Resin.                |

## VI Ordering Information

| <b>Product Name</b>                 | <b>Cat. No.</b> |
|-------------------------------------|-----------------|
| Monofinity A Resin                  | L00433          |
| Protein A Resin FF                  | L00464          |
| Protein G Resin FF                  | L00664          |
| Protein L Resin                     | L00239          |
| Protein G Resin FF Prepacked Column | L00681          |
| Protein A MagBeads MX               | L00672          |
| Protein G MagBeads MX               | L00673          |

**For research and manufacturing use. Direct human use, including taking orally and injection are forbidden.**

生产商：南京金斯瑞生物科技有限公司 江苏省南京市江宁区科学园雍熙路 28 号

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China