

Version: 02
 Update: 09/30/2022

Protein G Resin FF Prepacked Column

Cat No: L00681

I	Product Description	1
II	Operation	2
III	Purification Procedure	2
IV	Troubleshooting.....	2
V	Ordering Information	3

I Product Description

GenScript Protein G Resin FF is an affinity chromatography medium designed for easy, one-step purification of classes, subclasses and fragments of immunoglobulins from biological fluids and from cell culture media. The recombinant protein G ligand is coupled to 4% highly cross-linked agarose. The static binding capacity of Protein G Resin FF is greater than 20 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 G Resin FF Prepacked Column .

Protein G, a bacterial cell wall protein isolated from group G Streptococci, binds to mammalian IgGs mainly through Fc regions. Native protein G has 3 IgG binding domains and also sites for albumin and cell-surface binding. The latter have been eliminated from recombinant protein G to reduce nonspecific binding. Although protein G has very similar tertiary structures to protein A, their amino acid compositions differ significantly, resulting in different binding characteristics. Protein G can be used for purification of mammalian monoclonal and polyclonal IgGs that do not bind well to protein A. Protein G has greater affinity than protein A for most mammalian IgGs, especially for certain subclasses including human IgG3, mouse IgG1 and rat IgG2a. Unlike protein A, protein G does not bind to human IgM, IgD and IgA.

Table 1. Characteristics of Protein G Resin FF Prepacked Column

Resin Volume	1ml x 2(L00681-12);5ml x 1(L00681-51); 5ml x 5(L00681-55);
Ligand	Recombinant Streptococcal protein G lacking the albumin-binding sites expressed in <i>E. coli</i>
Number of IgG binding sites per ligand	3
M.W. of ligand	Approximately 22 kDa
PI of ligand	4.69
Degree of substitution	Approximately 2 mg protein G/ml settled resin
Static binding capacity	≥ 20 mg human IgG/ml settled resin
Matrix spherical	4% highly cross-linked agarose
Average particle size	90 μm (45-165 μm)
Recommended flow rate	50-300cm/h
Storage solution	20% ethanol; stored at 2-8 °C

II Operation

Buffer Preparation

Water and chemicals used for buffer preparation should be of high purity. We recommend filtering the buffers by passing them through a 0.45 µm filter before use.

Binding/Wash Buffer: 20 mM Na₂HPO₄, 0.15 M NaCl, pH 8.0

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) Apply the sample onto the column and drain the flow-through with a flow speed of about 1 ml/min. Collect the flow-through for measuring 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 immunoglobulins with 10-15 ml Elution Buffer and drain the eluate with a flow speed of about 1 ml/min. Collect the eluate 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.

Storage

Store regenerated Protein G Resin FF in 20% ethanol at 2°C to 8°C. **Do not freeze.**

IV Troubleshooting

Problem	Possible Cause	Solution
The flow rate of the column is very low (<0.5 ml/minute).	Tiny air bubbles from buffer or particles from sample block 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 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 G.	Try other affinity chromatography media to purify the antibody, such as Protein A Resin or Protein L Resin.

V Ordering Information

Product Name	Cat. No.
Monofinity A Resin	L00433
Protein A Resin FF	L00464
Protein G Resin FF	L00664
Protein L Resin	L00239
Protein A Resin FF Prepacked Column	L00680
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