chromatography

Protein A Removal From IgG on CHT[™] Ceramic Hydroxyapatite Support

Samuel G Franklin, PhD, Bio-Rad Laboratories, Inc., 2000 Alfred Nobel Drive, Hercules, CA 94547 USA

Introduction

Affinity chromatography with immobilized protein A from *Staphylococcus aureus* or from recombinant sources has been used for the purification of immunoglobulins since the 1970s (Gagnon 1996). Much of its popularity is due to the simplicity of the method itself, the high product purity and recovery, and protein A's selectivity to IgG class and species. Good clearance of DNA, virus, and endotoxin is often obtained, plus the immobilized protein A is relatively resistant to chemical regeneration and sanitization procedures.

Despite these advantages, conditioned media and cellular lysates frequently contain varying amounts of proteases, which can release fragments of protein A from immobilizing supports (Bloom et al. 1989). This is of great concern to the diagnostics and biopharmaceutical industries due to the wide range of immunomodulatory effects associated with protein A and its fragments (Gagnon 1996).

A number of chromatographic methods have been used to separate protein A from IgG. Cation exchange chromatography is attractive because conditions can be devised in which protein A does not bind to the support while IgG is retained. However, process development is often required and chaotropic agents may be needed to disrupt the protein A-IgG complex. Anion exchange chromatography has been used, but suffers from the disadvantage that the strength of the protein A-IgG association is enhanced by elevated pH. Since both IgG and protein A are retained, coelution of protein A fragments with IgG can occur (Bloom et al. 1989). Use of hydrophobic interaction chromatography for protein A-IgG separation has also been reported. Process development is again required since high salt concentrations increase the affinity of protein A for IgG. Coelution of protein A and IgG is common, reducing contaminant clearance and product yield.

Mariani et al. (1989) developed a simple method for the purification of IgG₁ in the presence of protein A on hydroxyapatite (Bio-Rad HPHT). They started with an ascites fluid ammonium sulfate fraction. In recent years, HPHT has been replaced by the more robust CHT ceramic hydroxyapatite, which was designed for process-scale separations. In this report, we show that CHT can resolve IgG₁ from an IgG₁-protein A complex in unfractionated conditioned media. We have also optimized conditions for the separation of protein A and an IgG₄ by cation exchange chromatography on UNOsphere[™] S support.

Results and Discussion

Initial studies in our laboratory indicated that recombinant protein A and IgGs from various sources exhibited virtually identical retention times on CHT under several elution conditions. We noted, however, that when small amounts of recombinant protein A were mixed with mouse ascites fluid containing IgG_1 , an additional peak of higher retention time than any other component in the system was generated. This finding inspired more detailed studies.

We obtained a conditioned medium containing about 0.24 mg/ml of mouse lgG_1 . When this medium or recombinant protein A was applied directly to a CHT column at pH 7.0 and eluted with a sodium phosphate gradient, peaks with essentially identical retention times were obtained (Figure 1, pink and green traces). However, upon mixing increasing amounts of recombinant protein A with the lgG_1 -containing medium, an additional peak was seen with a higher retention time than any other in the chromatogram. In concert, the lgG_1 peak appeared to diminish in intensity until, at the highest recombinant protein A concentration, most of the lgG_1 peak was depleted. This suggested that a stable complex between lgG_1 and protein A was being formed in these experiments and that this complex had a higher retention time than the other species present.



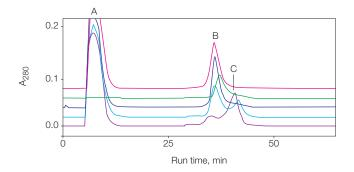


Fig. 1. CHT chromatography of conditioned medium containing murine IgG_1 and spiked with increasing amounts of recombinant protein A. Pink line, murine IgG_1 , 6.0 mg; green, recombinant protein A, 4.2 mg; dark blue, 0.48 mg $IgG_1 + 0.05$ mg protein A; light blue, 0.48 mg $IgG_1 + 0.1$ mg protein A; purple, 0.48 mg $IgG_1 + 0.2$ mg protein A. Proteins were mixed in a total volume of 6.0 ml buffer A (5 mM sodium phosphate buffer, pH 7.0). The column was a 7 x 50 mm Bio-Scale[™] column packed with CHT Type I, 20 µm (catalog #158-2000) equilibrated with buffer A. Elution was with a 20 column volume gradient to 200 mM sodium phosphate buffer, pH 7.0. at a flow rate of 1.5 ml/min. Detection was by absorption at 280 nm on a BioLogic[™] HR system. The fractions at the apex of peaks A, B, and C were analyzed by SDS-PAGE (Figure 2). Scale offset for comparison.

Mariani et al. (1989) did not attempt to demonstrate that the additional peak they had observed on HPHT upon mixing partially purified IgG_1 and protein A was actually a complex of the two proteins. To show this, we picked a ratio of IgG_1 to recombinant protein A of about 0.48 to 0.15 mg, which we had calculated would produce IgG_1 and IgG_1 -protein A complex peaks of approximately equal intensity on elution from the CHT column. The peaks were well resolved and of nearly equal height.

Fractions were collected and fractions from each of the three main peaks (A, B, and C) were examined with SDS-PAGE under reducing and nonreducing conditions. The gel was stained with an experimental fluorescent stain that is about 50 times more sensitive than Coomassie Blue (Figure 2). Lanes 2 and 3 of this gel show a commercially purified recombinant protein A standard, which appears to contain numerous trace impurities or degradation products. Lanes 4 and 5 contain the conditioned medium and show that IgG_1 is the main component. The unbound peak A (lanes 6 and 7)

appears to contain little or no protein and is probably indicator dye and other medium additives. Lanes 8 and 9 are from peak B and show this fraction to be the purified IgG_1 as expected. However, lanes 10 and 11 are from peak C, the putative IgG_1 -protein A complex, which show two bands in addition to the IgG_1 components. One of these bands comigrates with the main protein A band. The other band has a slightly higher mobility and is probably a proteolytic fragment of the recombinant protein A.

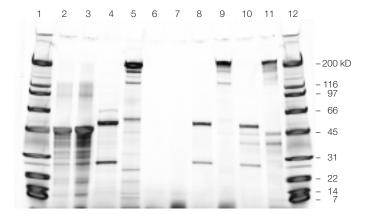


Fig. 2. SDS-PAGE analysis of CHT column fractions. Lanes 1 and 12, broad range molecular weight standards (catalog # 161-0317); lanes 2 and 3, protein A standard, reduced and unreduced; lanes 4 and 5, IgG₁ in conditioned medium, reduced and unreduced; lanes 6 and 7, peak A, reduced and unreduced; lanes 8 and 9, peak B, reduced and unreduced; lanes 10 and 11, peak C, reduced and unreduced. The gel was a 4–20% Ready Gel® precast gel.

In biopharmaceutical production processes, additional polishing steps are often included to ensure a pure product. We evaluated cation exchange chromatography on UNOsphere S as a potential additional polishing step.

Protein A has an isoelectric point of 5.1, and an examination of its calculated titration curve revealed that below pH 5.1, the negative charge on protein A increases rapidly, while above this pH, the change is much more gradual. Thus, protein A and a human IgG_4 both bind to UNOsphere S below pH 5.1. At pH 6.0, which does not favor interaction between protein A and IgG_4 , the two proteins could be separated (Figure 3). The proteins were run individually and mixed at various ratios, but retention times and peak shape did not vary significantly. This suggests that little or no protein A- IgG_4 complex was formed under these conditions (unpublished data).

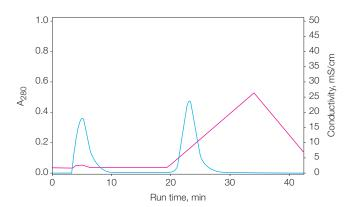


Fig. 3. Separation of recombinant protein A and human IgG_4 on UNOsphere S support. The protein A and IgG_4 (30 and 4.8 mg, respectively) were mixed in a total volume of 6.0 ml of buffer containing 50 mM MES, 1.0 mM NaH₂PO₄, pH 6.0. Proteins were applied to a 7 x 50 mm Bio-Scale column containing UNOsphere S (catalog # 156-0111). Elution was by a 20 column volume gradient from 0 to 0.5 M NaCl. Flow rate was 2.0 ml/min. Light blue line, A₂₈₀; pink line, conductivity. A BioLogic HR chromatography system was used.

Recovery of IgG from the protein A spiking experiment on CHT (see table) was about 64%. Almost 30% of the applied IgG was recovered in the protein A-IgG complex. About 85% of the IgG and 98% of the protein A applied to the UNOsphere S column was recovered.

We have shown that both supports offer high recovery and are effective in removing protein A from IgG from conditioned medium. These and other studies (Bio-Rad bulletins 2774 and 2780) also show high purity of the IgG recovered from a complex mixture using CHT and UNOsphere S. CHT and UNOsphere S chromatography are effective for protein A removal and as polishing steps in biopharmaceutical purification processes.

Table. Protein recovery from CHT and UNOsphere S columns.

	Protein Recovery (%)*		
Support	lgG	IgG-Protein A Complex	Protein A
CHT**	63.8	28.9 [‡]	ND
UNOsphere S [†]	84.8	ND	98.6

* Recovery calculated by A280 of all pooled fractions.

 ** The CHT column load was conditioned culture medium containing 0.48 mg $\rm IgG_1$ and 0.1 mg recombinant protein A. Conditions were as in Figure 1.

 † The UNOsphere S column load was 5.0 mg purified human IgG_4 and 30 mg recombinant protein A. Conditions were as in Figure 3.

 ‡ Expressed as $\% lgG_1$ recovered. This value was reduced by 10% to account for the protein A contribution to absorbance (the extinction coefficient of protein A is 10% that of lgG).

References

Bloom JW et al., Detection and reduction of protein A contamination in immobilized protein A purified monoclonal antibody preparations, J Immunol Methods 117, 83–89 (1989)

Gagnon P, Purification Tools for Monoclonal Antibodies, Chapter 9, Validated Biosystems, Inc., Tucson, AZ (1996)

Lai H and Franklin SG, Purification of transgenic antibody from corn seed using UNOsphere S and CHT ceramic hydroxyapatite supports, Bio-Rad bulletin 2774

Lai H and Franklin SG, Purification of murine IgG using UNOsphere S and CHT ceramic hydroxyapatite chromatography, Bio-Rad bulletin 2780

Mariani M et al., Purification of monoclonal antibodies for in vivo use: a twostep HPLC shows equivalent performance to immunoaffinity chromatography, Biochromatography 4, 149–155 (1989)

Coomassie is a trademark of Imperial Chemical Industries PLC.

Information in this tech note is current as of the date of writing (2002) and not necessarily the date of this revision or printing.



Bio-Rad Laboratories, Inc.

Life Science Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191 Canada (905) 712-2771 China (86-21) 63052255 Czech Republic (420) 2-4141 0532 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124) 6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30 Poland (48) 22-8126 672 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-2729877 South Africa 00 27 11 4428508 Spain 590 5200 Sweden 46 (0)8-55 51 27 00 Switzerland 061 717-9555 United Kingdom 0800-181134