

Purification of Murine IgG₁ Using UNOsphere™ S and CHT™ Ceramic Hydroxyapatite Chromatography

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

Introduction

Immunoglobulins (IgGs) from a variety of sources have been purified using the conventional process composed of a protein A chromatography step along with ion exchange chromatography and hydrophobic interaction chromatography steps (Jiskoot et al. 1989, Godfrey et al. 1993, Shadle et al. 1995, Ford et al. 2001). Protein A chromatography is normally applied as the first step in the process, yielding antibodies with very high purity in a single step. However, the disadvantages of protein A chromatography are: 1) antibody isomers are not well separated from different species; 2) protein A leakage requires additional chromatographic steps for protein A removal; and 3) protein A resins are substantially more expensive than ion exchange and ceramic hydroxyapatite supports.

Here we present an alternative process for antibody purification consisting of UNOsphere S chromatography and CHT ceramic hydroxyapatite chromatography. This process is simpler than the process utilizing protein A and avoids the leakage of a contaminating ligand (Jiskoot et al. 1989).

UNOsphere S support is a strong cation exchanger made from acrylamido and vinylic copolymers. It exhibits high protein binding capacity and low column backpressure at high linear velocity. Thus, it is a suitable capture resin in the downstream purification process. CHT ceramic hydroxyapatite support, with Ca²⁺ ions and PO₄³⁻ ions in its spherical and ceramic structure, provides excellent resolution and selectivity for protein isolation.

Methods and Results

The BioLogic DuoFlow™ workstation was used to develop the processes. An analytical column (0.7 x 5 cm) was loaded with UNOsphere S support for small-scale process development.

A preparative column (2.2 x 20 cm) was used in process scale-up. All SDS-PAGE was performed using Ready Gel® precast gels, the Criterion™ cell, and the PowerPac™ 3000 power supply. Ceramic hydroxyapatite (Type I, 10 μm and 20 μm) media were used for this study.

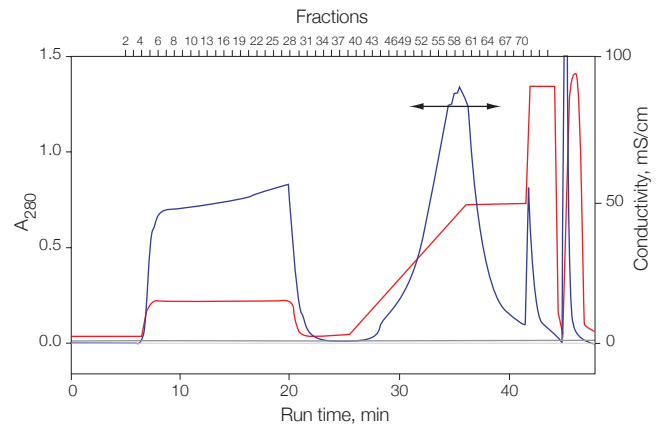


Fig. 1. Chromatogram of murine IgG₁ purification on the UNOsphere S column during small-scale process development. Diluted cell culture (20 ml) was loaded onto an UNOsphere S 0.7 x 5 cm column at a flow rate of 600 cm³/hr in 20 mM phosphate-citrate buffer, pH 4.0. The sample was eluted in 10 column volumes (CV) of a 0–0.5 M NaCl gradient, followed by 5 CV of 1 M NaCl in the same buffer. The column was then cleaned in 1 M NaOH. The double-headed arrow indicates the fractions (1 ml each) containing IgG₁ (see Figure 2). Blue trace, A₂₈₀; red trace, conductivity profile.

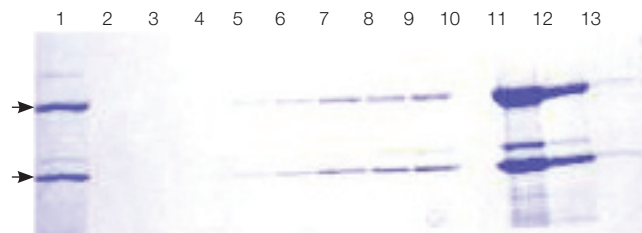


Fig. 2. SDS-PAGE analysis of murine IgG₁ purification by UNOsphere S chromatography during small-scale process development. Lane 1, sample load; lanes 2–9, flow-through fractions 4, 6, 8, 11, 12, 13, 21, 29; lane 10, wash fraction 40; lanes 11, 12, elution fractions 56, 61; lane 13, fraction after cleaning column with NaOH. Arrows indicate IgG₁ bands.

Small-Scale Process Development

UNOsphere S chromatography was used as the capture step for murine IgG₁ isolation. The binding capacity of UNOsphere S support for murine IgG₁ was optimized at pH values from 7.0 to 4.0. Decreasing the pH in the cell culture increased the binding capacity of UNOsphere S support for murine IgG₁ (data not shown). The murine IgG₁ was stable for at least 6 hr in buffer at pH 4.0 (data not shown), so the cell culture was adjusted to pH 4.0. The cell culture was then loaded onto the UNOsphere S column as described in Figure 1. The murine IgG₁ was eluted in the NaCl gradient. The fractions were analyzed by SDS-PAGE (Figure 2). It was found that Phenol Red and some contaminating proteins were removed in the flow-through pool by UNOsphere S chromatography.

For ceramic hydroxyapatite chromatography, the murine IgG₁ fractions were pooled and diluted 5-fold with 1 mM sodium phosphate buffer, pH 6.8. The murine IgG₁ elution pool from the UNOsphere S column could not be loaded directly on the CHT supports because the ionic strength of the pooled solution was too high to allow murine IgG₁ to bind. The diluted sample was loaded onto the CHT column at 300 cm/hr. Some contaminants were removed by a NaCl gradient. The murine IgG₁ was eluted by a phosphate gradient (Figure 3).

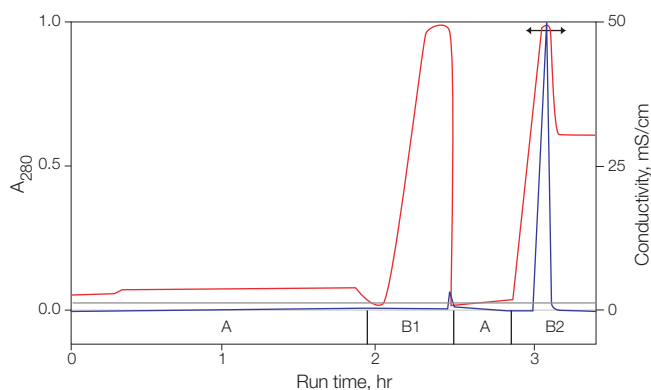


Fig. 3. Chromatogram of murine IgG₁ purification by CHT chromatography during small-scale process development. The murine IgG₁ pool (100 ml) from the UNOsphere S column (fractions indicated with arrow in Figure 1) was loaded onto a CHT-I 10 μ m, 0.7 x 5 cm column at a flow rate of 300 cm/hr in 1 mM sodium phosphate, pH 6.8 (buffer A). The sample was then eluted with a gradient of 0–0.5 M NaCl in buffer A over 10 CV followed by 5 CV of 0.5 M NaCl in buffer A (step B1), and 10 CV of buffer A. Finally, the column was eluted with 10 CV of a gradient of 1 mM–0.4 M sodium phosphate buffer, pH 6.8 (step B2). The arrow indicates the pooled fractions containing IgG₁. Blue trace, A₂₈₀; red trace, conductivity.

The fraction pool was analyzed by SDS-PAGE (Figure 4). We estimated that >95% purity of murine IgG₁ was achieved in this purification process.

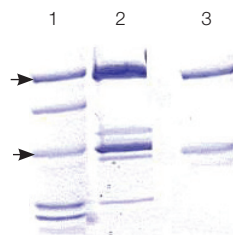


Fig. 4. SDS-PAGE analysis of the murine IgG₁ elution pool from the CHT ceramic hydroxyapatite column during small-scale process development. Lane 1, protein molecular weight standards (250, 150, 100, 75, 50, 37, 25, 15, and 10 kD); lane 2, UNOsphere S elution pool; lane 3, CHT elution pool. Arrows indicate IgG₁ bands.

Process Scale-Up

A 2.2 x 20 cm column was packed with 60 ml of UNOsphere S to a bed height of 16 cm and run as described (Figure 5). Some contaminants were removed, and the murine IgG₁ could be isolated in the process at this scale. The results indicated that the UNOsphere S purification process could be scaled up at least 30 times.

The murine IgG₁ pool from the UNOsphere S column was further purified by CHT chromatography as described in Figure 6. Murine IgG₁ was eluted in a phosphate gradient applied after some contaminants were removed by a NaCl gradient.

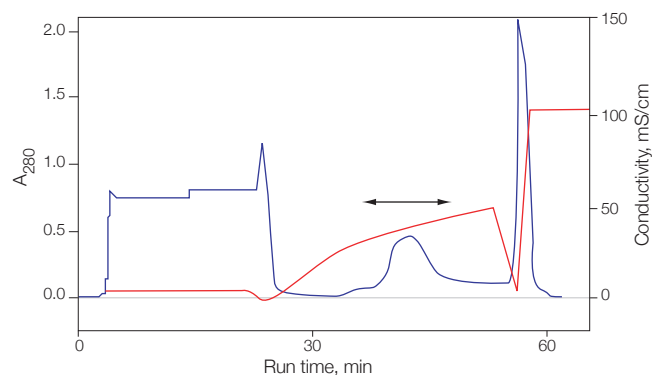


Fig. 5. Chromatogram of murine IgG₁ purification by UNOsphere S chromatography during scale-up. The column was equilibrated to pH 4.0 with 1 mM sodium acetate, pH 4.0 (buffer A). The pH of the cell culture sample was adjusted to 4.0 with concentrated acetic acid, then loaded onto the column at 300 cm/hr, the same linear velocity used in the small-scale purification process. Once 600 ml of pH-adjusted sample had been loaded, the column was washed with buffer A. The murine IgG₁ was eluted with 20 CV of a 0–0.5 M NaCl gradient in buffer A. The column was then cleaned with 10 CV of 1 M NaCl in buffer A followed by 5 CV of 0.5 M NaOH. The arrow indicates the pooled fractions containing IgG₁. Blue trace, A₂₈₀; red trace, conductivity.

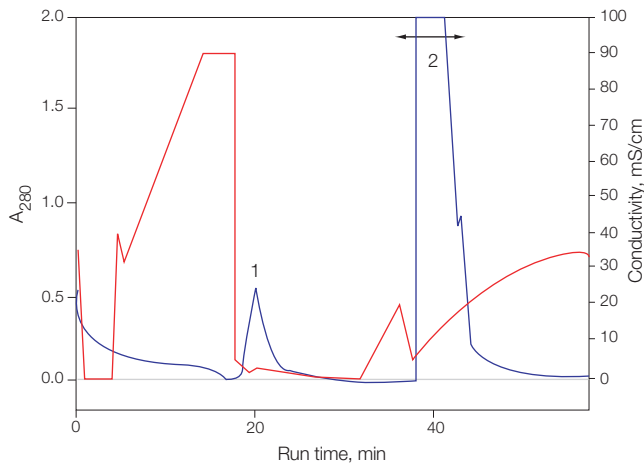


Fig. 6. Chromatogram of murine IgG₁ purification by CHT chromatography during scale-up. Murine IgG₁ fractions from the UNOsphere S column (indicated by arrow in Figure 5) were pooled, diluted 5 times to 1.85 L with 1 mM sodium phosphate buffer, pH 6.8 (buffer A), and loaded onto a 20 ml CHT Type I column (20 μ m, 1.1 x 10 cm) at 300 cm/hr. The column was washed with buffer A, then eluted with a 0–1 M NaCl gradient in buffer A for 10 CV followed by 1 CV of 1 M NaCl. The column was then washed again with buffer A for 5 CV. A 1 mM–0.4 M sodium phosphate gradient (pH 6.8) was run for 20 CV to elute the murine IgG₁. The arrow indicates the pooled fractions containing IgG₁. Blue trace, A₂₈₀; red trace, conductivity.

SDS-PAGE analysis showed that the scaled-up CHT purification process yields murine IgG₁ of similar purity to that in small-scale purification (Figure 7). The murine IgG₁ was isolated by phosphate gradient; however, some minor bands were observed in lane 3 when the gel was overloaded with 40 μ g of a murine IgG₁ elution pool. These bands suggest that there are still some degradation products or contaminants in the purified murine IgG₁. Average recovery in this scale-up purification process was about 80%. Data from flow cytometry analyses suggest that the murine IgG₁ solution from our purification process yields very similar immunospecificity to processes using protein A chromatography (data not shown).

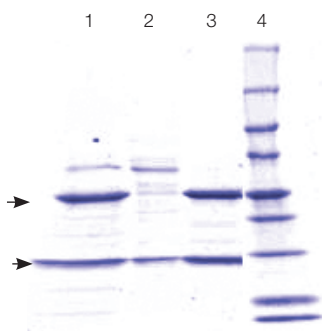


Fig. 7. SDS-PAGE analyses of CHT column fractions during scale-up. Lane 1, load; lane 2, peak 1 elution pool; lane 3, peak 2 elution pool; lane 4, molecular weight standards. Arrows indicate IgG₁ bands.

Acknowledgement

Special thanks to Paul Yazaki, PhD, Beckman Research Institute of the City of Hope National Medical Center, Duarte, CA, for his numerous contributions to this project.

References

- Ford CH et al., Affinity purification of novel bispecific antibodies recognising carcinoembryonic antigen and doxorubicin, *J Chromatogr B Biomed Sci and Appl* 754, 427–435 (2001)
- Godfrey MAJ et al., Assessment of the suitability of commercially available SpA affinity solid phases for the purification of murine monoclonal antibodies at process scale, *J Immunol Methods* 160, 97–105 (1993)
- Jiskoot W et al., Two-step purification of a murine monoclonal antibody intended for therapeutic application in man. Optimisation of purification conditions and scaling up, *J Immunol Methods* 124, 143–156 (1989)
- Shadle PJ et al., Antibody purification, US patent 5,429,746 (1995)

Information in this tech note was current as of the date of writing (2002) and not necessarily the date this version (rev A, 2006) was published.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000
