

# Removal of Aggregate from an IgG<sub>4</sub> Product Using CHT™ Ceramic Hydroxyapatite Resin

Tech  
Note

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

## Process Separations

Bulletin 2940

### Introduction

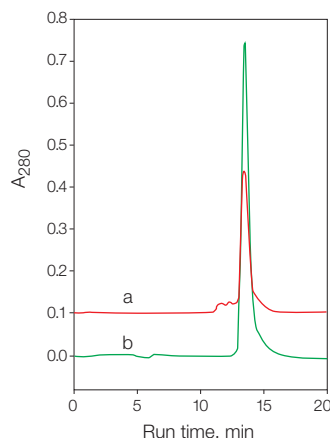
The presence of even trace amounts of aggregate in immunoglobulin biopharmaceuticals is of significant concern during all phases of drug development, from process design to quality control. The formation of aggregates can adversely impact process economics by causing decreased product yield, peak broadening (requiring the addition of polishing steps), and loss of activity (Gagnon 1996). Aggregate formation can also adversely affect product safety because it can cause complement activation or even anaphylaxis upon injection (Ritchie and Navolotskaia 1996, WHO Report 1995).

Most chromatographic techniques would be expected to be able to remove aggregated or multimeric species of immunoglobulin from the monomer. However, techniques such as ion exchange and hydrophobic interaction chromatography may also induce the formation of additional aggregate or multimer due to increased protein concentration or changes in salt concentration, pH, or both required for elution. Size exclusion chromatography does not suffer from these disadvantages, but does result in significant dilution of the product. If large quantities of product are required, large size exclusion columns, concentration steps, and support equipment will add considerably to production costs. In some cases, additional product loss due to denaturation may occur during dilution.

CHT Ceramic Hydroxyapatite exhibits electrostatic, repulsive, and coordinate covalent bond formation in interacting with protein species. We have explored its potential in removing aggregate from a purified IgG<sub>4</sub> biopharmaceutical known to contain aggregate, and report here on a simple and effective process to do so.

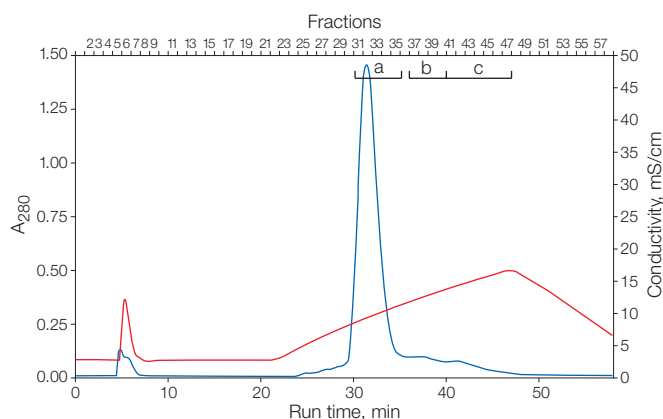
### Results

Size exclusion chromatography analysis of the IgG<sub>4</sub> sample (derived from chromatogram a, Figure 1) indicated that it contained 86.6% monomer of 141.9 kD and 10.1% aggregate, which consisted of nearly equal amounts of 485.3 kD and >1.1 MD species. Possible degradation products with apparent molecular weights below 3,000 were also present (about 1.0% of total protein).



**Fig. 1. Size exclusion chromatography analysis of IgG<sub>4</sub> samples.**

Chromatogram a (—), starting material; chromatogram b (—), main peak (Figure 2, peak a) from the CHT Column (scales are offset for comparison). Column, Bio-Sil® SEC 250-5, 7.8 x 300 mm (catalog #125-0062); buffer, 0.1 M sodium phosphate, pH 6.8, containing 0.15 M NaCl; flow rate, 1.0 ml/min. Molecular weights were calculated by calibration with standards (IgG, ovalbumin, myoglobin, vitamin B<sub>12</sub>, catalog #151-1901)



**Fig. 2. CHT chromatography of IgG<sub>4</sub> sample containing aggregate.**

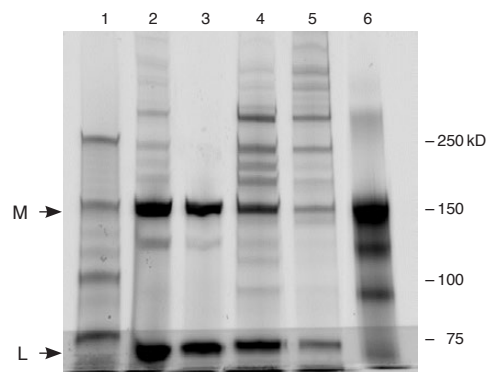
Chromatography was performed on a BioLogic DuoFlow™ 10 System (catalog #760-0037). Sample (12 mg) was loaded onto a 7 x 52 mm Bio-Scale™ MT Column (catalog #751-0081) containing CHT, 20 μm (catalog #158-2000) in 20 mM sodium phosphate, pH 7.0; elution was from 20 to 160 mM sodium phosphate in 20 column volumes. The flow rate was 1.5 ml/min, and the fraction size was 1.5 ml. Peaks were pooled as indicated in the figure: a, fractions 31–35; b, fractions 37–40; c, fractions 41–47. A<sub>280</sub> (—); conductivity (—).

**BIO-RAD**

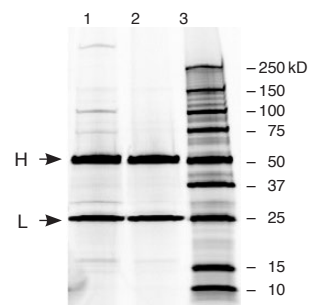
The IgG<sub>4</sub> sample was applied to a CHT Column and eluted with a linear phosphate gradient. A single major peak, flanked by two or three broader minor peaks, was obtained (Figure 2).

The major peak from the CHT Column (peak a) was analyzed on a size exclusion column and was found to consist of only monomeric IgG<sub>4</sub> (Figure 1, chromatogram b). Spectrophotometric analysis at 280 nm indicated that it represented about 72.4% of the protein applied to the CHT Column. The major CHT column pools were also analyzed by nonreducing 5% SDS-PAGE (Figure 3).

A major band comigrating with the 150 kD standard was presumed to be IgG<sub>4</sub> monomer. Some nine or ten components with molecular weights greater than that of the IgG<sub>4</sub> monomer were apparent in the CHT column load (Figure 3, lane 2) and were presumed to be aggregates. The major CHT column peak (lane 3) appeared to be free of these aggregates and



**Fig. 3.** Nonreducing SDS-PAGE analysis of CHT column pools. Lane 1, MW standards; lane 2, IgG<sub>4</sub> starting material; lane 3, main CHT column peak (peak a); lane 4, CHT peak b; lane 5, CHT peak c; lane 6, human IgG (Sigma). The gel was a 5% Ready Gel® Precast Gel (catalog #161-1213), and the stain was an experimental stain about 50-fold more sensitive than Coomassie Blue. Positions of the IgG<sub>4</sub> monomer (M) and light chain (L) are indicated by arrows.



**Fig. 4.** Reducing SDS-PAGE analysis of CHT column pools. Lane 1, IgG<sub>4</sub> starting material; lane 2, main CHT column pool (peak a); lane 3, MW standards. The gel was a 4–20% Ready Gel Precast Gel (catalog #161-1159); the stain was the same as in Figure 3. Positions of the IgG<sub>4</sub> heavy (H) and light (L) chains are indicated by arrows.

had a trace of a single lower molecular weight species, which may be the F(ab')<sub>2</sub> fragment. Lanes 4 and 5, which contained the later eluting minor peaks (b and c, respectively) from the CHT Column, contained an array of aggregates and some lower molecular weight species, which may be IgG<sub>4</sub> degradation products. A major band, presumed to be the IgG<sub>4</sub> light chain, ran slightly ahead of the 75 kD standard under these conditions. Some of the components observed in the IgG<sub>4</sub> fractions were also found in a commercially available human IgG preparation (lane 6).

A reducing 4–20% SDS gel was also run (Figure 4), which shows that some aggregated species were present under these conditions as well and may have arisen from covalent crosslinking. These components were also removed by CHT chromatography. The heavy and light chains migrated as expected in this gel.

## Discussion

The results reported here indicate that CHT column chromatography is capable of removing essentially all of the aggregation and degradation products found in a human IgG<sub>4</sub> biopharmaceutical in a single step.

The aggregates found in this preparation did not appear to be simple multimers of the 141.9 kD IgG<sub>4</sub> monomer. The absence of simple multimers may have been due to the formation of various complexes with the monomer, F(ab')<sub>2</sub> fragment, and other degradation products. Both disulfide and nonreducible crosslinks (possibly enzyme-mediated) appeared to exist in this preparation.

The CHT Column performed well at a relatively high sample load (6 mg/ml) and with a relatively short bed height (52 mm). No reaggregation was evident in these experiments. The procedure is simple and should be easily scaled up. This suggests CHT chromatography may be useful for immunoglobulin polishing in numerous circumstances.

## References

- Gagnon P (1996). Hydroxyapatite Chromatography. In Purification Tools for Monoclonal Antibodies (Arizona: Validated Biosystems, Inc.), p. 95.
- Ritchie RF and Navolotskaia O (1996). Serum Proteins in Clinical Medicine, Vol 1 (Maine: Foundation for Blood Research).
- Veterinary Public Health Unit (1995). Report of a WHO Consultation on Intradermal Application of Human Rabies Vaccines (Geneva, Switzerland: World Health Organization).

Information in this tech note was current as of the date it was written (2002), not necessarily the date this version (Rev B, 2015) was printed.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000