



CHT

CHROMATOGRAPHY

CHT™ Ceramic Hydroxyapatite

Ca²⁺, (PO₄)³⁻

- High throughput
- High physical and chemical stability
- Unparalleled selectivity
- High reproducibility
- 3 particle sizes for any application

High-Resolution Protein Separations Using CHT Ceramic Hydroxyapatite

Summary

Traditional crystalline hydroxyapatite is widely known for its unique separation properties and unparalleled selectivity. The fragile nature of the crystalline material has, however, limited its usefulness, particularly for large-scale processes. CHT ceramic hydroxyapatite overcomes the physical and chemical limitations of traditional crystalline hydroxyapatite and provides the throughput, stability, and reproducibility required for industrial biopharmaceutical manufacturing.

CHT ceramic hydroxyapatite is a chemically pure form of hydroxyapatite that has been sintered at high temperatures to yield a physically and chemically robust support. It possesses high protein binding capacities and can be used reproducibly for hundreds of cycles in large columns run at high flow rates. It is available in two distinct material types,

Type I and Type II (see table), and three particle sizes, 20, 40, and 80 μm (see figure). Both types retain the separation properties of crystalline hydroxyapatite as well as possessing some unique properties of their own. Type I CHT has a higher protein binding capacity than Type II, and in particular, a higher capacity for acidic proteins. Type II often provides superior selectivity and resolution in the separation of many proteins, particularly various species, subspecies, and classes of antibodies, as well as nucleic acids.

The two types are often evaluated side by side to determine which material provides the maximum benefit in a given separation. Existing protocols that have been developed on crystalline hydroxyapatite can often be applied directly to the ceramic material with little or no modification.

Table. Properties of CHT ceramic hydroxyapatite.

Characteristic	Type I	Type II
Functional groups	Ca ²⁺ , (PO ₄) ³⁻	Ca ²⁺ , (PO ₄) ³⁻
Dynamic binding capacity	>25 mg lysozyme per gram CHT	>12.5 mg lysozyme per gram CHT
Typical IgG binding capacities at 500 cm/hr	10–50 mg/ml	5–25 mg/ml
Particle sizes available	20, 40, and 80 μm (nominal)	20, 40, and 80 μm (nominal)
Nominal pore diameter	600–800 Å	800–1,000 Å
Maximum operating pressure	60 bar (900 psi)	100 bar (1,500 psi)
Recommended operating flow rates	10–1,500 cm/hr	10–1,500 cm/hr
pH stability	5.5–14	5.5–14
Regeneration	in 0.4–1.0 M phosphate buffer	in 0.4–1.0 M phosphate buffer
Sanitization	in 1–2 N NaOH	in 1–2 N NaOH
Autoclavability	20 min at 121°C	20 min at 121°C
Chemical Compatibility		
1 N NaOH	>24 hr	>24 hr
8 M urea	>24 hr	>24 hr
8 M guanidine-HCl	>24 hr	>24 hr
Ethanol or methanol	>24 hr	>24 hr
100% acetonitrile	>24 hr	>24 hr

Note: A small amount of an appropriate counterion such as 0.1–10 mM NaPO₄⁻ should be added to all unbuffered solutions.



BIO-RAD

Mechanism of Action and Standard Chromatography

CHT ceramic hydroxyapatite interacts with biomolecules by multiple modes. Electrostatic interactions occur between positively charged calcium ions and negatively charged phosphate groups on the support and charged moieties on the biomolecule's surface. Much stronger coordination complexes can form between clusters of carboxyl groups (on proteins) and calcium sites on CHT. Repulsion effects and the geometric charge distribution on CHT provide unique selectivity. Typically, acidic, basic, and neutral proteins are bound to hydroxyapatite using a low ionic strength phosphate buffer. Elution is accomplished through the use of a phosphate gradient of increasing strength. Regeneration of the support with high-strength phosphate buffer is followed by sanitization with base. For more detailed information, refer to the instruction manual.

Storage and Shelf Life

CHT ceramic hydroxyapatite should be stored in 20% ethanol v/v containing a small amount of an appropriate counterion such as 0.1–10 mM NaPO_4^- . Storage in phosphate buffer containing a bacteriostatic agent such as sodium azide (NaN_3) is also possible. Ceramic hydroxyapatite can be stored indefinitely in dry form at room temperature when sealed in the original container.

Technical Assistance

All of the CHT ceramic hydroxyapatite supports have manufacturing processes registered with the United States Food and Drug Administration (FDA) by submission of a Type II Drug Master File (DMF). Regulatory support files are available upon request to companies entering into clinical trials. Bio-Rad Laboratories is an ISO 9001 registered corporation. For additional information and technical assistance, contact your local Bio-Rad office. In the USA and Canada, call 1-800-4BIORAD.

Visit us on the Web at www.bio-rad.com for more information on Bio-Rad's complete line of process chromatography supports and other products for life science research and production.

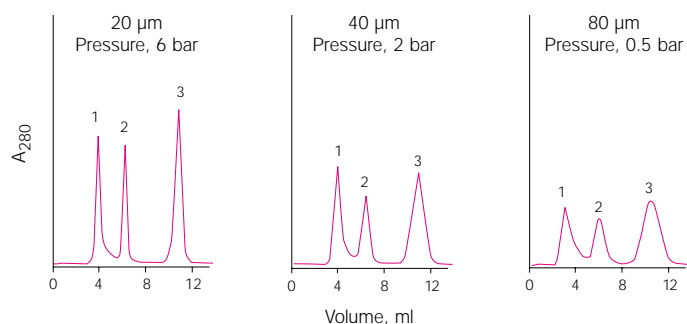


Figure. Effect of particle size on separation of proteins. A 10 µl sample of 10 mg/ml BSA (peak 1), 1.3 mg/ml lysozyme (peak 2), and 5 mg/ml cytochrome c (peak 3) was run on each 4 x 100 mm column packed with the indicated particle size at a flow rate of 478 cm/hr. The elution buffer was a linear gradient of 1–400 mM sodium phosphate, pH 6.8, over 15 min.

Ordering Information

Catalog #	Description
158-2000	CHT Ceramic Hydroxyapatite, Type I, 20 µm, 10 g
157-0020	CHT Ceramic Hydroxyapatite, Type I, 20 µm, 100 g
157-0021	CHT Ceramic Hydroxyapatite, Type I, 20 µm, 1 kg
157-0025	CHT Ceramic Hydroxyapatite, Type I, 20 µm, 5 kg
158-4000	CHT Ceramic Hydroxyapatite, Type I, 40 µm, 10 g
157-0040	CHT Ceramic Hydroxyapatite, Type I, 40 µm, 100 g
157-0041	CHT Ceramic Hydroxyapatite, Type I, 40 µm, 1 kg
157-0045	CHT Ceramic Hydroxyapatite, Type I, 40 µm, 5 kg
158-8000	CHT Ceramic Hydroxyapatite, Type I, 80 µm, 10 g
157-0080	CHT Ceramic Hydroxyapatite, Type I, 80 µm, 100 g
157-0081	CHT Ceramic Hydroxyapatite, Type I, 80 µm, 1 kg
157-0085	CHT Ceramic Hydroxyapatite, Type I, 80 µm, 5 kg
158-2200	CHT Ceramic Hydroxyapatite, Type II, 20 µm, 10 g
157-2000	CHT Ceramic Hydroxyapatite, Type II, 20 µm, 100 g
157-2100	CHT Ceramic Hydroxyapatite, Type II, 20 µm, 1 kg
157-2500	CHT Ceramic Hydroxyapatite, Type II, 20 µm, 5 kg
158-4200	CHT Ceramic Hydroxyapatite, Type II, 40 µm, 10 g
157-4000	CHT Ceramic Hydroxyapatite, Type II, 40 µm, 100 g
157-4100	CHT Ceramic Hydroxyapatite, Type II, 40 µm, 1 kg
157-4500	CHT Ceramic Hydroxyapatite, Type II, 40 µm, 5 kg
158-8200	CHT Ceramic Hydroxyapatite, Type II, 80 µm, 10 g
157-8000	CHT Ceramic Hydroxyapatite, Type II, 80 µm, 100 g
157-8100	CHT Ceramic Hydroxyapatite, Type II, 80 µm, 1 kg
157-8500	CHT Ceramic Hydroxyapatite, Type II, 80 µm, 5 kg

Larger quantities available on request.

BIO-RAD

**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 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 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