

CHROMATOGRAPHY

Macro-Prep[®] HIC Support

- Reverse-phase selectivity with maintenance of biological activity
- Efficient capture of proteins from high-salt conditions

Increase Throughput and Efficiency

Summary

Hydrophobic interaction chromatography (HIC) separates proteins on the basis of relative hydrophobicity. HIC is a natural second step after either ion exchange or salt precipitation since the sample is applied in high-salt buffer. At high ionic strengths, hydrophobic sites of the protein interact with the alkyl groups of the support. Retention, selectivity, and biological activity are somewhat dependent on pH, type of salt used, and its concentration.

The Macro-Prep methyl HIC support is ideal for purification of proteins with strongly hydrophobic regions. The Macro-Prep t-butyl HIC support is ideal for purification of proteins with few or weakly hydrophobic regions. The two different ligands provide alternative selectivities for easier optimization of separation (Figure 1 and Table 1). The properties of the supports are summarized in Table 2. These methacrylate copolymer beads provide high resolution at very high flow rates. They can be sanitized quickly and efficiently in 0.15% peracetic acid (Figure 2), and are compatible with many common solutions useful in HIC (Table 3). Changes in pH or ionic strength of the buffer do not cause shrinking or swelling of the support.

Table 1. Comparison of retention times using Macro-Prep methyl and t-butyl ligands.

Peak	Protein	Retention Times, min	
		Methyl	t-Butyl
1	Cytochrome c	7.30	10.41
2	Ovalbumin	32.25	40.22
3	α -Amylase	62.81	71.50
4	Ferritin	81.49	90.90

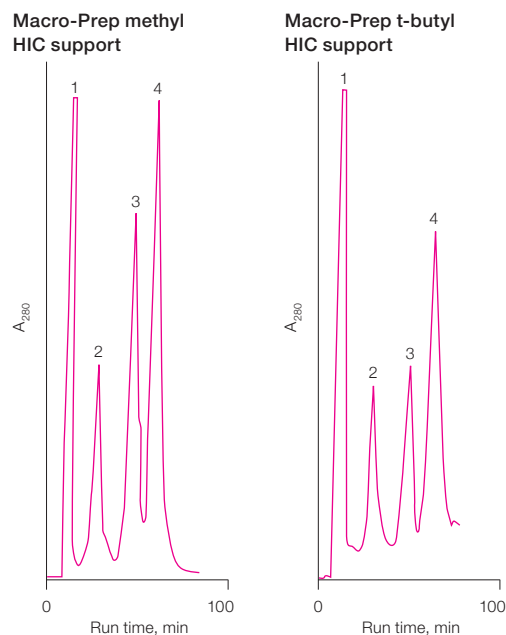


Fig. 1. Sample separation of proteins on Macro-Prep methyl and t-butyl supports. A 50 μ l sample (0.6 mg total protein) of cytochrome c (peak 1), ovalbumin (peak 2), α -amylase (peak 3), and ferritin (peak 4) was run on a 1 x 10 cm (4.5 ml) column at a linear flow rate of 38 cm/hr. The sample was eluted with 0.1 M Na_2PO_4 , pH 7.0, 1.85 M $(\text{NH}_4)_2\text{SO}_4$ for 10 min, followed by a gradient from 1.85 M to 0 M $(\text{NH}_4)_2\text{SO}_4$ over 90 min. The retention times for each protein are shown in Table 1 for comparison.



Table 2. Properties of Macro-Prep HIC supports.

	Methyl	t-Butyl
Type of support	HIC	HIC
Functional group	-CH ₃	-C(CH ₃) ₃
Binding capacity*	>25 mg/ml	>15 mg/ml
Nominal particle size	50 µm	50 µm
Max. linear flow rate	3,000 cm/h	3,000 cm/h
Autoclavability	121°C	121°C
pH stability	1–10	1–10
Regeneration	70% ethanol	70% ethanol
Sanitization	0.15% peracetic acid	0.15% peracetic acid

* Determined with human serum albumin (HSA).

Table 3. Compatibility of Macro-Prep HIC supports.

	Methyl	t-Butyl
1% SDS	Yes	Yes
8 M guanidine-HCl	Yes	Yes
1 M HCl	Yes	Yes
100% ethanol	Yes	Yes

Recommended Procedure

The Macro-Prep HIC supports are easy to use. Rinse out the ethanol solution with 2–3 bed volumes of double distilled water and equilibrate in starting buffer. Samples are typically loaded on the column in the pH range of 4–9 and in a salt concentration of 0.5 to 2 M ammonium sulfate or NaCl. Protein binding increases with both ligand hydrophobicity and salt concentration. In theory, less salt is required to salt out a protein when the pH is close to its isoelectric point. Temperature influences the hydrophobic interaction; protein binding is increased at elevated temperatures (40°C) and reduced at low temperatures (4°C). The salt concentration is gradually decreased to elute bound protein. Elution is followed by stripping and regeneration steps of choice. For more detailed information, refer to the instruction manual.

Technical Assistance

The Bio-Rad Life Sciences Group and its design, development, and manufacture of chemicals and analytical instruments, have been assessed and registered by National Quality Assurance Limited against the provisions of BS EN ISO:9001:2000. For additional information and technical assistance, contact your Bio-Rad representative.

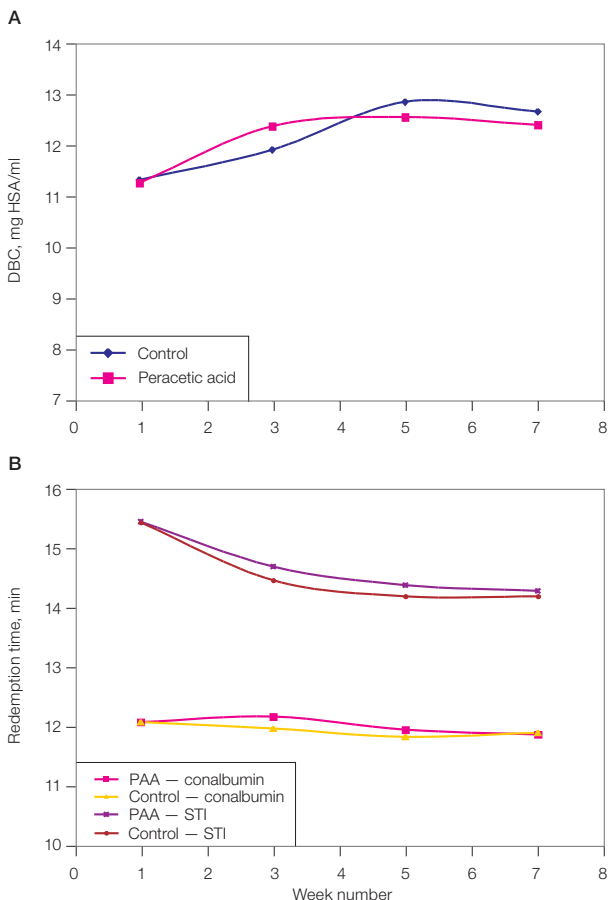


Fig. 2. Macro-Prep HIC supports can be sanitized extensively with 0.15% peracetic acid. A, Dynamic binding capacity; B, retention time.

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

Ordering Information

Catalog #	Description
158-0080	Macro-Prep Methyl HIC Support, 25 ml
156-0080	Macro-Prep Methyl HIC Support, 100 ml
156-0081	Macro-Prep Methyl HIC Support, 500 ml
158-0090	Macro-Prep t-Butyl HIC Support, 25 ml
156-0090	Macro-Prep t-Butyl HIC Support, 100 ml
156-0091	Macro-Prep t-Butyl HIC Support, 500 ml

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 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