
Macro-Prep[®] t-Butyl and Methyl Hydrophobic Interaction Chromatography Media

Instruction Manual

Please read these instructions prior to using Macro-Prep hydrophobic interaction chromatography media. If you have any questions or comments regarding these instructions, please contact your local Bio-Rad Laboratories representative.

The Bio-Rad logo consists of the words "BIO-RAD" in a bold, white, sans-serif font. The text is centered within a black, rounded rectangular background.

Table of Contents

Section 1	Introduction	1
Section 2	Intended Use	1
Section 3	Media Screening.....	1
Section 4	Technical Specifications and Characteristics	2
Section 5	Preparation for Use	3
Section 6	Column Packing.....	3
Section 7	Column Packing Evaluation.....	5
Section 8	Operation and Maintenance	7
Section 9	Ordering Information	10

Section 1

Introduction

All Bio-Rad products are manufactured under an externally approved ISO 9001 quality system. Bio-Rad Laboratories guarantees product quality and performance of unopened product for 1 year from date of shipment. Regulatory support files are available for Macro-Prep hydrophobic interaction chromatography (HIC) media. If you need assistance validating the use of Macro-Prep media in a production process, contact your local Bio-Rad Process Chromatography representative.

Section 2

Intended Use

Macro-Prep HIC media are designed specifically for intermediate purification steps that remove host-cell contaminants from partially purified targets. Due to their rigidity and unique surface chemistry, these media are particularly suited for HIC operations requiring high throughput and high recovery of target.

Both Macro-Prep t-butyl and Macro-Prep methyl HIC media operate on a mechanism of interaction that is based on hydrophobicity and charge.

The t-butyl and methyl groups are mildly hydrophobic. Depending on the chosen pH of loading and elution buffers, the carboxyl groups can be exploited to ionically repel target molecules while the hydrophobic groups retain contaminants. This is an ideal strategy to minimize product losses commonly experienced with HIC media due to denaturation of proteins when exposed to hydrophobic surfaces.

Section 3

Media Screening

For ease of screening of Bio-Rad HIC media, both forms of Macro-Prep HIC media are provided as prepacked 1 ml and 5 ml cartridges. In addition, these media are part of the media sampler pack, which also includes ion exchangers and ceramic hydroxyapatite media. See Ordering Information in Section 9 at the end of this instruction manual.

With Macro-Prep HIC media it is important to test both salt and pH conditions. Since HIC is usually used after an affinity or ion exchange step, it may be necessary to adjust the salt concentration and/or the pH of the sample load to obtain optimal purification and recovery with Macro-Prep HIC media.

Section 4 Technical Specifications and Characteristics

Macro-Prep t-butyl and Macro-Prep methyl media are based on rigid, epoxy-activated methacrylate beads derivatized with t-butyl or methyl groups, respectively, and carboxyl groups (Figure 1). These products are shipped as a slurry of 50% settled media in 20% ethanol. Table 1 lists the properties of the two media. The rigidity of these media allows high flow rates without bed compression (Figure 2). Testing in small laboratory-scale columns (10 mm in diameter x 100 mm in height) showed no bed compression at a pressure of 7 bar and a flow rate of 40 ml/min (3,000 cm/hr). This property of the media provides for high throughput during loading, column washing, equilibration, and cleaning.

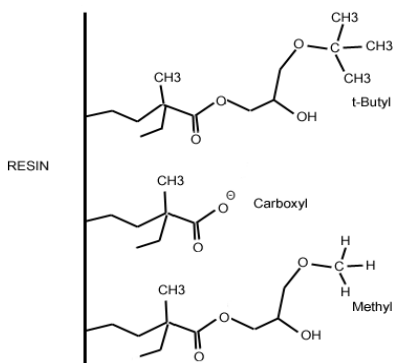


Fig. 1. Macro-Prep HIC: Macro-Prep t-butyl contains t-butyl and carboxyl groups; Macro-Prep methyl contains methyl and carboxyl groups.

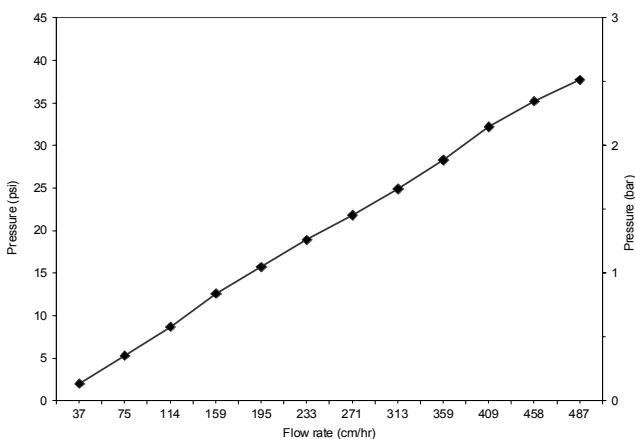


Fig. 2. Pressure vs. flow for Macro-Prep media. Media bed dimensions were 140 mm diameter by 250 mm length.

Table 1. Characteristics of Macro-Prep HIC Media

	Macro-Prep t-Butyl	Macro-Prep Methyl
Functional groups	COO- and t-butyl	COO- and methyl
Ionic capacity	Approximately 120 µeq/ml	Approximately 120 µeq/ml
Shipping counterion	Cl ⁻	Cl ⁻
Hydrophobicity	Mild	Mild
Median particle size (µm)	50	50
pH stability	1–10	1–10
Autoclavability (121°C, 30 min)	Yes	Yes
Shipping solution	20% ethanol	20% ethanol
Working flow-rate range	100–600 cm/hr	100–600 cm/hr
Sanitization	5 CV of 6 M guanidine-HCl, 100 cm/hr or 5 CV of 1% acetic acid in 0.12 M phosphoric acid, pH 1.5, 100 cm/hr	5 CV of 6 M guanidine-HCl, 100 cm/hr or 5 CV of 1% acetic acid in 0.12 M phosphoric acid, pH 1.5, 100 cm/hr
Storage at 4–40°C	20% ethanol or 1% acetic acid in 0.12 M phosphoric acid, pH 1.5	20% ethanol or 1% acetic acid in 0.12 M phosphoric acid, pH 1.5

Section 5 Preparation for Use

Macro-Prep HIC media are supplied fully hydrated in 20% ethanol as a 50% slurry. Small volumes of Macro-Prep HIC media are easily washed in a Büchner funnel with 4–5 column volumes (CV) of a low ionic strength buffer (<20 mM). For large process-scale volumes, pack the column in a low ionic strength buffer (<20 mM) or water and then rinse with at least 3 CV of the low ionic strength buffer or water before transferring to the high-salt buffer. Alternatively, decant the ethanol shipping solution and resuspend/decant 3–4 times in the low ionic strength buffer to be used for column packing.

Section 6 Column Packing

Columns can be packed in different ways depending on the media and type of column and equipment to be used. Macro-Prep HIC media can be packed using pressure, flow, or suction packing methods. To pack highly efficient columns, it is recommended to use a 50% slurry. Some general guidelines for packing small and large columns are given below. Make sure to read and follow the instructions for the column to be used. If unsure about a particular column, call your local Bio-Rad Laboratories representative. Throughout this instruction, flow rate is expressed linearly in centimeters per hour (cm/hr). The relationship between linear and volumetric flow rate as delivered by a pump is:

Linear flow rate (cm/hr) X Cross-sectional area of the column (cm²) = Volumetric flow rate

Example: 300 cm/hr in a 14 cm diameter column = 300 cm/hr X πr^2 or 300 cm/hr X (3.14 X 7²) cm² or 46,158 cm³/hr or 769.3 ml/min.

Packing Small Columns

The following slurry packing method was designed to pack 25 ml of Macro-Prep HIC media into 5–15 mm ID, low- to medium-pressure lab columns. All buffers should be degassed prior to packing. Use a packing reservoir that holds ~100 ml.

1. Prepare degassed buffer or water to be used as the packing solution.
2. Macro-Prep is shipped as a 50% slurry, so measure 50 ml of suspended slurry into a 100 ml graduated cylinder. Allow the media to settle. Decant the shipping solution from the settled media.
3. Add 25–50 ml of degassed packing buffer to the media.
4. Seal the cylinder and rotate it to suspend the media. Caution: Do not mix with a magnetic stir bar, to avoid formation of small fines. Larger volumes of slurry may be mixed with an overhead stirrer at low to moderate speed.
5. Fill the column with approximately 2 cm of packing buffer. Pour all of the media slurry into the column with the packing reservoir attached.
6. Insert the column flow adaptor and start flow at ~100 cm/hr. Gradually increase the flow rate from 100 cm/hr to the flow rate at which the maximum allowable pressure for the column is reached, or to the flow rate at which the maximum pressure for Macro-Prep is reached (7 bar or 102 psi). As the bed compresses with increasing flow rate, stop the flow and adjust the flow adaptor so it penetrates 1–2 mm into the media bed. Repeat the flow and adjustment cycle until the bed no longer compresses.
7. Attach the column to your chromatography system and purge with starting buffer at the maximum flow rate to be used during operation. You should not see any further bed compression. If you do, adjust the flow adaptor 1–2 mm into the media bed.
8. The column is now ready for use.

Note: For optimal long-term use, do not exceed 75% of the maximum pressure attained during the column packing.

Packing Large Columns

In large columns, Macro-Prep should be packed using a 50% slurry at constant pressure up to a maximum of 7 bar (102 psi). Pack with a constant flow rate that is approximately 25% greater than the anticipated maximum to be used during operation. Given the variety of industrial column hardware and packing skids, we recommend use of standard operating procedures for packing provided by the column vendor.

Bio-Rad EasyPack™ and GelTec™ columns are supplied with a movable piston. They are packed by pumping the packing solution through the media bed at constant pressure or flow.

1. Packing begins with preparation of the tubing and the column. Clean and rinse the column well with packing buffer before packing. The rinse is also used to displace air from the column and its exit tubing. Make sure there is no air trapped in the bottom flow cell. After removal of all the air from the bottom flow cell and exit tubing, leave about 2 cm of liquid in the bottom of the column. Close the outlet valve(s).
2. In a suitable container, mix the packing buffer with the media to form a 50% slurry.
3. Pour or pump the slurry into the column.
4. Insert the movable piston and lower it to the surface of the liquid, making sure no air is trapped under the piston.
5. Pressurize the inflatable seal to 2 bar (29 psi), so when lowering the piston into the slurry, the buffer exits the top of the column through the inlet line. Be sure the inlet line is completely filled with packing buffer to avoid air being pumped back into the column.
6. Connect a pump and a pressure gauge, open all inlet and outlet valves, and start packing at constant flow rate or pressure. Keep the flow rate or pressure constant throughout the packing. Check the pressure at the column inlet. Do not exceed the pressure limit of column or media.
7. Lower the piston as the bed compresses.
8. When the media bed no longer compresses, mark the bed height on the column tube, close the outlet valve, and stop the pump. The bed will start to expand in the column.
9. Lower the piston to within 1 cm of the surface of the media bed. Fully inflate the seal (4 bar), start the pump, open the valves, and continue packing.
10. Repeat steps 8 and 9 until there is a maximum of 1 cm between media bed surface and piston when the media bed is stable.
11. Close the bottom valve, stop the pump, disconnect the column inlet, and lower the piston to the gel bed surface. The column is ready for testing of column packing efficiency.

Section 7

Column Packing Evaluation

To check the packing efficiency and to monitor the column's status over time, packing efficiency and asymmetry should be tested immediately after packing and at regular intervals to ensure reproducible performance. The values to be determined for the packed column are the height equivalent theoretical plate (HETP), number of theoretical plates (N), and the asymmetry factor (A_s). The method used involves applying a test sample of a low molecular weight substance

that has no interaction with the media, e.g., 2 M NaCl. When using salt as the test substance, use a concentration of 2 M NaCl in water with 0.5 M NaCl in water as the elution buffer. Concentrated buffer solutions, e.g., 10x buffer concentrate, is also used. HETP varies depending on the test conditions. Since it is used as a reference value, it is important to keep conditions and equipment constant so results can be compared over time. Changes in solute, solvent, elution buffer, sample volume, flow rate, flow path, temperature, etc. influence results. For optimal results, ensure that the sample volume does not exceed 2.5% of the column volume, and maintain the flow rate between 75 and 150 cm/hr. If an acceptance limit is defined in relation to column performance, the column HETP and/or calculated plate number, N , can be used as acceptance criteria for column use.

If a UV absorbing substance is used as the test sample, use a UV absorbance monitor set at 280 nm. If 2 M NaCl is the probe, use a conductivity monitor; the elution buffer should be 500 mM NaCl. Apply the sample as close to the column inlet as possible to avoid dilution in the tubing and piping leading up to the column inlet. Calculate HETP and the number of theoretical plates (N) as follows:

$$\text{HETP} = L/N$$

$$N = 5.54(V_e/W_{1/2h})^2$$

L = Bed height (cm)

N = Number of theoretical plates

V_e = Peak elution volume or time

$W_{1/2h}$ = Peak width at peak's half height in volume or time

V_e and $W_{1/2h}$ should always be the same units

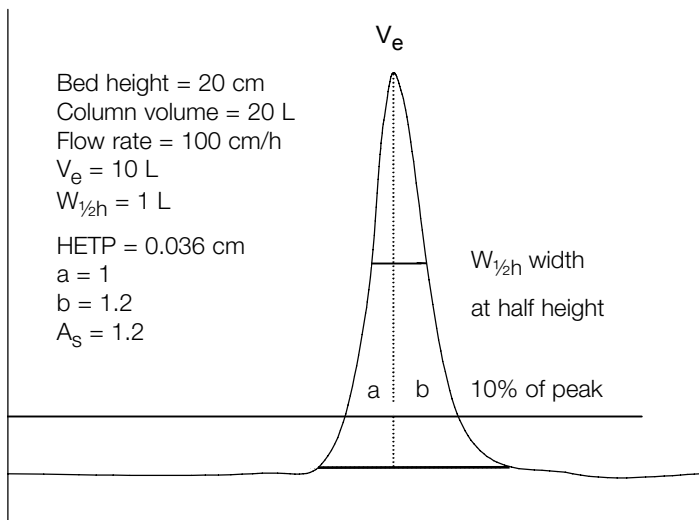


Fig. 3. A simulated chromatography profile from which HETP and A_s values are calculated.

Peaks should be symmetrical. The best achievable asymmetry factor is 1. Values of 0.8 to 1.8 are excellent for large columns >30 cm in diameter. If functional performance is acceptable, A_S factors more distant from 1 put less demand on the column packing procedure in production. A change in the shape of the peak is usually the first indication of deteriorating performance. Peak asymmetry factor calculation:

$$A_S = b/a$$

a = Front section of peak width at 10% of peak height bisected by line denoting V_e .

b = Latter section of peak width at 10% of peak height bisected by line denoting V_e .

Section 8 Operation and Maintenance

Macro-Prep media were designed to remove trace contaminants while operating at high throughput. Macro-Prep media should be run at the highest linear velocities and loading capacities while staying within the pressure limits of the column and chromatography system. For flow-through mode, a linear flow rate of 300 cm/hr in a 20 cm bed is a reasonable starting point. If the product is being retained on the Macro-Prep, flow rates of 100–200 cm/hr are recommended. Optimization can be achieved by changing the pH, the ionic strength, or additives in the elution buffer.

All buffers commonly used for cation exchange chromatography can be used with Macro-Prep HIC media (see Table 2).

Table 2. Common Buffers for Macro-Prep HIC.

Buffer	Useful pH Range
Citric acid	2.6–3.6
Lactic acid	3.6–4.3
Formic acid	3.8–4.3
Acetic acid	4.8–5.2
MES	5.5–6.7
PIPES	6.1–7.5
MOPS	6.5–7.9
Phosphate	6.7–7.6
Bicine	8.2–8.7
HEPES	6.8–8.2
TES	6.8–8.2
Tricine	7.4–8.8

Column Equilibration

Higher salt concentrations are used to enhance hydrophobic interactions. Normally, a column is equilibrated in a high-salt buffer, and the sample is loaded onto the column at the same pH and salt concentration. At small scale, the following salts are commonly used: ammonium sulfate (<pH 8), 1 M solutions of sodium sulfate, sodium chloride, potassium chloride, or sodium citrate. At larger scale, sodium citrate is a very practical salt. This salt can function as both a buffer and a salt during HIC. Lower molar quantities are required to reach specific ionic concentrations and the somewhat hydrophobic nature of the citrate ion often

minimizes product loss due to overly strong product interaction with the media. If the target molecule is in the flow-through, sodium citrate is often used at concentrations of 500–600 mM for loading.

Sample Preparation

Adjust salt concentration and pH as necessary for desired binding of target or contaminants. This is best done by mixing the correct amount of liquid concentrate of salt into the sample load and then adjusting the pH.

Sample Load and Adsorption

The sample load is determined empirically by loading and evaluating breakthrough of target or contaminants. Since HIC is an adsorption technique, the sample volume is not a critical factor. Large volumes of very dilute feed such as cell culture supernatant and clarified lysates can be loaded onto the media without prior concentration. If the salt concentration of the sample load is lower than what is used to equilibrate the column before loading, the protein of interest can show low binding capacities, or can begin to elute from the column with unwanted contaminants.

Wash through

After loading of the column, follow with 2 to 3 CV of the high ionic strength buffer. This will wash out any unbound materials.

Elution

Elute target molecules with either a step or linear gradient. Usually, the salt concentration at which the desired product binds is predetermined at small scale. With this knowledge, the pH and salt concentration used in wash-through are adjusted to eliminate the maximum amount of contamination before starting elution of the target.

Regeneration

After each run, the packed bed should be washed with 3–5 CV of a low ionic strength buffer such as 20 mM sodium citrate or 20 mM HEPES. This is followed with 3–5 CV of water. If contamination remains, wash with 2–3 CV of 0.12 M phosphoric acid in 1% acetic acid, pH 1.5. For very difficult cleaning, follow the acid wash with a combination of nonionic detergents and then 30–70% ethanol.

Cleaning-in-Place (CIP)

If a column no longer yields reproducible results, the media may require a thorough CIP to remove strongly bound contaminants. Acceptable CIP agents include 25% acetic acid, 8 M urea, 1% Triton X-100, 70% ethanol, 30% isopropyl alcohol, 1 N HCl, and 6 M guanidine-HCl. Cleaning with NaOH is not recommended because the t-butyl and methyl groups are cleaved from the media under strongly basic conditions. If it is absolutely necessary to use NaOH, use 0.1 N for 1 hr, and then wash immediately with a low pH, high-salt buffer such as 600 mM sodium citrate, pH 5.

Sanitization and Long Term Storage

Sanitization is the reduction of bioburden, i.e. microorganisms and spores in the column. Before long-term storage, Macro-Prep HIC media should be cleaned and sanitized.

Procedure:

Cleaning: After elution, clean the column with 8 CV of 1% acetic acid in 0.12 M phosphoric acid, pH 1.5, <100 cm/hr. If not satisfactorily cleaned, continue to wash with 5 CV of 6 M guanidine-HCl, <100 cm/hr. If still not clean, wash the column with 5 CV of 20% ethanol, <100 cm/hr.

To sanitize and store: Wash the column with 5 CV of 1% acetic acid in 0.12 M phosphoric acid, pH 1.5. Store the column at 4–40°C.

Section 9

Ordering Information

Macro-Prep HIC media are available in a variety of package sizes, as prepacked Econo-Pac® cartridges for easy testing and small-scale purifications, and in the media sampler pack.

Macro-Prep Methyl Products

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
156-0082	Macro-Prep Methyl HIC Support, 5 L
156-0083	Macro-Prep Methyl HIC Support, 10 L
732-0051	Macro-Prep Methyl HIC Support, 1 x 5 ml Econo-Pac cartridge
732-0053	Macro-Prep Methyl HIC Support, 5 x 1 ml Econo-Pac cartridges
732-0055	Macro-Prep Methyl HIC Support, 5 x 5 ml Econo-Pac cartridges
158-0100	Media Sampler Pack, contains one 25 ml bottle each of UNOsphere™ Q and S ion exchangers, Macro-Prep methyl and t-butyl media, and Macro-Prep High Q, High S, DEAE, and CM ion exchangers, one 10 g bottle each of CHT™ ceramic hydroxyapatite types I and II

Macro-Prep t-Butyl Products

Catalog #	Description
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
156-0092	Macro-Prep t-Butyl HIC Support, 5 L
156-0093	Macro-Prep t-Butyl HIC Support, 10 L
732-0056	Macro-Prep t-Butyl HIC Support, 1 x 5 ml Econo-Pac cartridge
732-0057	Macro-Prep t-Butyl HIC Support, 5 x 5 ml Econo-Pac cartridges
732-0058	Macro-Prep t-Butyl HIC Support, 5 x 1 ml Econo-Pac cartridges
158-0100	Media Sampler Pack, as described above

Ask your representative for pricing details, larger volumes, and special packaging.

Triton is a trademark of Union Carbide Chemicals & Plastics Technology Corporation.

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Dr., Hercules, CA 94547 USA

510-741-1000

LIT486 Rev C