

# High-Resolution Separation of a Fusion Protein

Zhang Tingting,<sup>1</sup> Dai Zhenyu,<sup>1</sup> Xu Qun,<sup>1</sup> and Jeffrey Rohrer<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, Shanghai, People's Republic of China;

<sup>2</sup>Thermo Fisher Scientific, Sunnyvale, CA, USA

## Key Words

MABPac HIC-10 Column, Hydrophobic Interaction Chromatography (HIC), HPLC, Protein Drug, Isomer, Impurity

## Goal

To develop a high-resolution HIC method for the separation of a fusion protein from a truncated version and a structural isomer

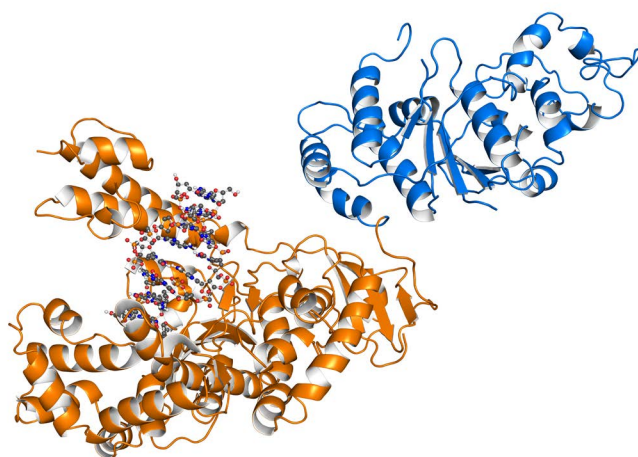
## Introduction

Fusion proteins are proteins created through the joining of two or more genes that originated from separate proteins. Fusion proteins created through genetic engineering impart properties from each of the parent proteins. Fusion protein therapeutics (e.g., Fc-fusion proteins) have proven successful in clinical treatments.<sup>1</sup>

Because fusion protein drugs are protein therapeutics, they require high-resolution methods for characterization and quality control techniques that are significantly different and more challenging than those used for small molecule drugs. HIC provides unique separation power based on protein surface hydrophobicity. Compared to other commercially available HIC columns, the Thermo Scientific™ ProPac™ HIC-10 column has demonstrated superior separation power for discriminating post-translational modifications of monoclonal antibodies, especially those due to methionine oxidation.<sup>2</sup> Here, a new HIC column, the Thermo Scientific™ MABPac™ HIC-10, provides impressive separation of a fusion protein from both a truncated version and a structural isomer—resolutions that cannot be achieved using existing commercially available HIC columns.

## Equipment

- Thermo Scientific™ Dionex™ UltiMate™ 3000 Dual Rapid Separation Liquid Chromatography (RSLC) system,\* including:
  - DGP-3600AB Biocompatible Dual Ternary Low-Pressure Proportioning Analytical Pump System (P/N 5037.0014) with SRD-3600 Integrated Solvent and Degasser Rack, 6 Channels (P/N 5035.9230)
  - WPS-3000TRS Wellplate Sampler, Thermostatted (P/N 5840.0020)
  - TCC-3000SD Thermostatted Column Compartment (P/N 5730.0010)
  - DAD-3000 Diode Array Detector, Without Flow Cell (P/N 5082.0010)



Or

- UltiMate 3000 Dual Bio LC system,\* including:
  - DGP-3600BM Biocompatible Dual-Gradient Micro Pump (P/N 5042.0066)
  - WPS-3000TBFC Thermostatted Biocompatible Pulled-Loop Well Plate Autosampler with Integrated Fraction Collection (P/N 5825.0020)
  - TCC-3000SD Thermostatted Column Compartment (P/N 5730.0010)
  - DAD-3000 Diode Array Detector, Without Flow Cell (P/N 5082.0010); with Analytical Flow Cell for DAD-3000 and MWD-3000 Series, SST, 13 µL Volume, 10 mm Path Length (P/N 6082.0100)
- Thermo Scientific™ Dionex™ Chromeleon™ Chromatography Data System software, version 7.2 or higher
- Thermo Scientific™ Orion™ 2-Star Benchtop pH Meter

\* This application was successfully evaluated on the two UltiMate systems listed.

**Thermo**  
SCIENTIFIC

Conditions	
Columns:	ProPac HIC-10 Analytical, 5 $\mu$ m, 4.6 $\times$ 100 mm (P/N 063655) MABPac HIC-10 Analytical, 5 $\mu$ m, 4.6 $\times$ 100 mm (P/N 088480)
Mobile Phases:	A. 2.0 M ammonium sulfate, 0.1 M sodium phosphate, pH 7.0/isopropanol (93:7 v/v) B. 0.1 M sodium phosphate, pH 7.0/isopropanol (93:7 v/v)
Gradient:	Routine, 0–30 min, 5–100% B Shallow, 0–5 min, 30% B; 5–30 min, 30–60% B
Flow Rate:	1.0 mL/min
Temperature:	30 $^{\circ}$ C
Detection:	UV, absorbance at 214 nm

## Results and Discussion

The fusion protein sample was donated by a customer and had three fractions: the fusion protein, the fusion protein minus several amino acid residues, and the fusion protein with a different tertiary structure. As shown in Figure 1, these three fractions were not resolved using a popular commercially available HIC column.

Figure 2 shows separations of the fusion protein sample using the ProPac HIC-10 and MABPac HIC-10 columns under the routine gradient conditions. Three peaks were observed when using the ProPac HIC-10 column; their separation was further improved using the MABPac HIC-10 column.

Experiments showed that a lower starting salt concentration and shallower gradient improved the separation. As shown in Figure 3 and Table 1, the three components were better resolved using the shallower gradient.

For complicated fusion protein drug samples, it is critical to obtain pure fractions for further characterization. These samples often contain structural forms of the protein. High resolution is also needed for quality control of therapeutic proteins. The MABPac HIC-10 column provides another option for the separation of challenging fusion protein samples.

Table 1. Peak resolution (Rs).

Column	Peak 1	Peak 2	Peak 3
ProPac HIC-10	—	NA*	NA*
MABPac HIC-10 (Routine Gradient)	—	1.60	1.53
MABPac HIC-10 (Shallow Gradient)	—	1.66	1.61

\*NA = not applicable

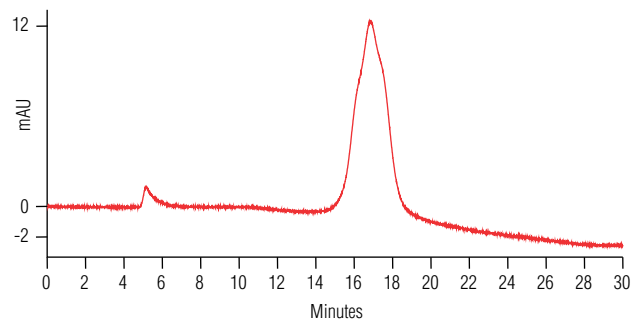


Figure 1. The fusion protein sample separated using a commercially available HIC column (chromatogram provided by the customer without conditions).

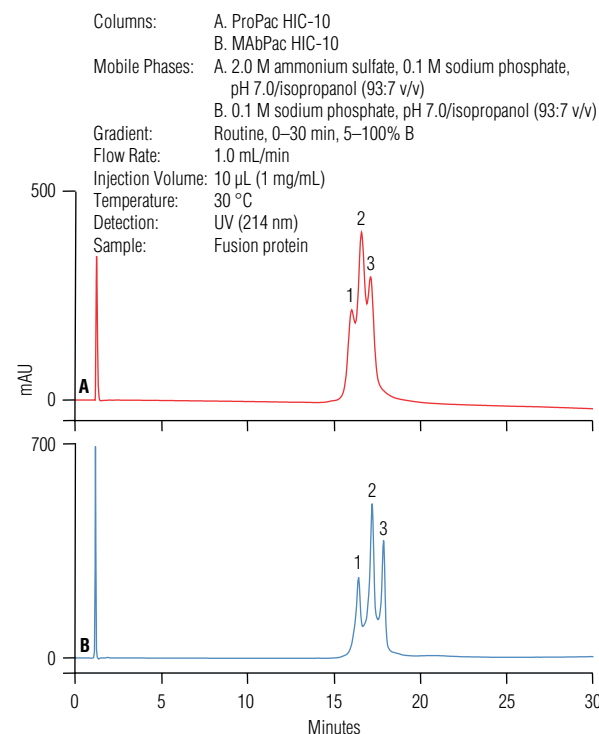


Figure 2. The fusion protein sample separated using (A) a ProPac HIC-10 column and (B) a MABPac HIC-10 column with a routine gradient.

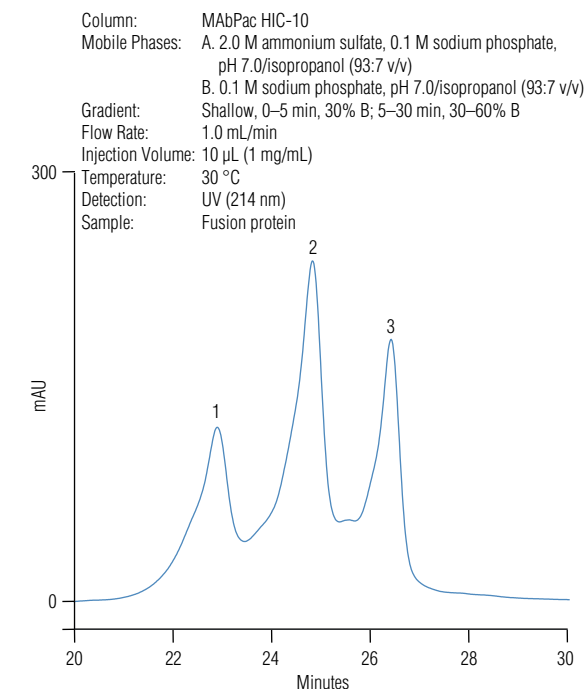


Figure 3. The fusion protein sample separated using a MABPac HIC-10 column with a shallow gradient.

## References

1. Beck, A.; Reichert, J.M. Therapeutic Fc-Fusion Proteins and Peptides as Successful Alternatives to Antibodies. *MAbs*, 2011, 3 (5), 415–416.
2. Valliere-Douglass, J.; Wallace, A.; Balland, A. Separation of Populations of Antibody Variants by Fine Tuning of Hydrophobic-Interaction Chromatography Operating Conditions. *J. Chromatogr., A* 2008, 1214, 81–89.

[www.thermofisher.com/chromatography](http://www.thermofisher.com/chromatography)

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

<b>Africa</b> +43 1 333 50 34 0	<b>Denmark</b> +45 70 23 62 60	<b>Japan</b> +81 6 6885 1213	<b>Russia/CIS</b> +43 1 333 50 34 0
<b>Australia</b> +61 3 9757 4300	<b>Europe-Other</b> +43 1 333 50 34 0	<b>Korea</b> +82 2 3420 8600	<b>Singapore</b> +65 6289 1190
<b>Austria</b> +43 810 282 206	<b>Finland</b> +358 9 3291 0200	<b>Latin America</b> +1 561 688 8700	<b>Sweden</b> +46 8 556 468 00
<b>Belgium</b> +32 53 73 42 41	<b>France</b> +33 1 60 92 48 00	<b>Middle East</b> +43 1 333 50 34 0	<b>Switzerland</b> +41 61 716 77 00
<b>Brazil</b> +55 11 3731 5140	<b>Germany</b> +49 6103 408 1014	<b>Netherlands</b> +31 76 579 55 55	<b>Taiwan</b> +886 2 8751 6655
<b>Canada</b> +1 800 530 8447	<b>India</b> +91 22 6742 9494	<b>New Zealand</b> +64 9 980 6700	<b>UK/Ireland</b> +44 1442 233555
<b>China</b> 800 810 5118 (free call domestic) 400 650 5118	<b>Italy</b> +39 02 950 591	<b>Norway</b> +46 8 556 468 00	<b>USA</b> +1 800 532 4752

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand