

High-Resolution Separation of Cysteine-Linked Antibody-Drug Conjugate Mimics Using Hydrophobic Interaction Chromatography

Julia Baek and Xiaodong Liu, Thermo Fisher Scientific, Sunnyvale, CA, USA

Key Words

Hydrophobic interaction chromatography, HIC, monoclonal antibody, mAb, antibody-drug conjugate, ADC, drug-to-antibody ratio, DAR, MAbPac HIC-Butyl

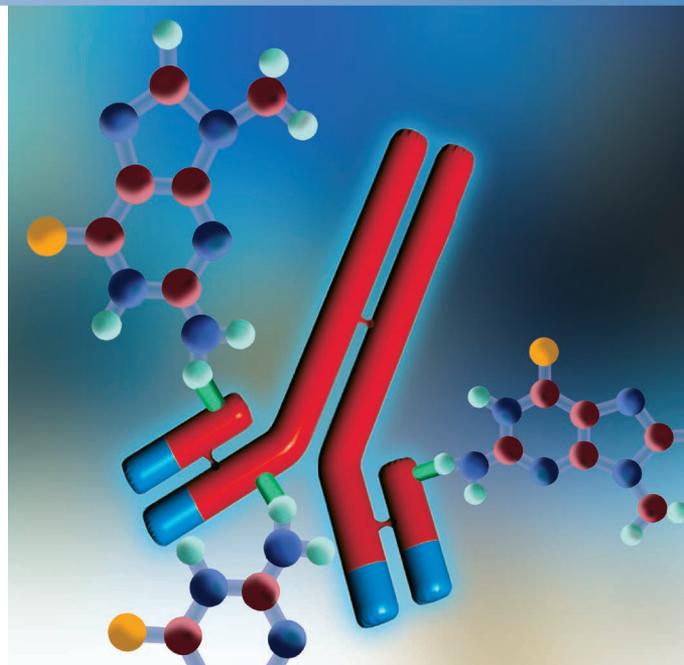
Goal

To demonstrate the analysis of antibody-drug conjugate (ADC) mimics using a Thermo Scientific™ MAbPac™ HIC-Butyl column.

Introduction

Monoclonal antibody drugs are the fastest growing class of biotherapeutics and have become a major part of the biopharmaceutical market. Monoclonal antibodies have demonstrated their effectiveness against autoimmune disorders, cardiovascular diseases, infectious diseases, and cancer.¹ Monoclonal antibodies selectively recognize antigens that are present on the tumor cells and exert their cytotoxic effect by receptor agonist activity, immune response, or vascular and stromal cell ablation. However, most of these mAbs are used in combination with chemotherapy, and many others have shown a lack of clinical efficacy. The success and the limitation of monoclonal antibodies have fueled the development of another form of antibody based therapeutics—antibody-drug conjugates (ADCs).² Antibody-drug conjugates utilize the exquisite selectivity of the antibody to achieve targeted delivery of cytotoxic drugs. ADCs have gained tremendous interest among pharmaceutical companies due to their significantly improved clinical efficacy over native monoclonal antibodies.³

The conjugation of drugs often results in an ADC molecule that is heterogeneous with respect to both the distribution and loading of cytotoxic drugs on the mAb. The number of drugs attached to the mAb has been shown to directly affect the safety and the efficacy of the drug. Unconjugated mAbs have significantly lower potency, and the ADCs with high drug load are subject to rapid renal clearance. Therefore, it is critical to fully characterize and monitor the heterogeneity of ADCs during development and production.⁴



Hydrophobic interaction chromatography (HIC) is a technique for separation of proteins, including monoclonal antibodies, by hydrophobicity. The HIC mobile phase usually consists of a salting-out agent, which at high concentration retains the protein by increasing hydrophobic interaction between the protein and the stationary phase. Bound proteins are eluted by decreasing the salt concentration.

HIC has been widely used as an orthogonal method to size-exclusion chromatography and ion-exchange chromatography for the characterization of mAb heterogeneity. Analysis of succinimides, antibody fragments, oxidated mAbs, and C-terminal lysine modifications were successfully carried out to monitor the stability and, in some cases, the potency of the drug.^{5,6} HIC is also suitable for the separation of ADCs since attachment of cytotoxin alters the hydrophobicity of the antibody. The least hydrophobic unconjugated antibody

elutes first and as the number of drugs attached increases the elution time increases. Therefore, HIC is considered to be the method of choice to characterize the distribution of ADC molecules with different drug-to-antibody ratios (DARs).⁷

The MAbPac HIC-Butyl column is a polymer-based HIC column designed for separation of mAbs and variants, including ADCs. The hydrophilic nature of polymer particles and the optimal density of butyl functional groups lead to excellent biocompatibility, low carryover, and high resolution. Here we describe the analysis of ADC samples that were conjugated via interchain disulfide bonds on the MAbPac HIC-Butyl column.

Experimental

Chemicals and Reagents

- Deionized (DI) water, 18.2 M Ω -cm resistivity
- 2-Propanol (Fisher Scientific P/N A461-4)
- Sodium phosphate monobasic monohydrate (NaH₂PO₄•H₂O, $\geq 98.0\%$)
- Ammonium sulfate [(NH₄)₂SO₄, $\geq 99.0\%$]

Sample Handling Equipment

Polypropylene, 0.3 mL vials (P/N 055428)

Sample Preparation

Cysteine-conjugated ADC mimic samples were provided by a customer. 10 mg/mL ADC mimic samples were diluted in half with 2 M (NH₄)₂SO₄, 100 mM NaH₂PO₄, pH 7.0 solution. 25 mg/mL control mAb was diluted five-fold with 2 M (NH₄)₂SO₄, 100 mM NaH₂PO₄, pH 7.0 solution.

LC Separation

The LC separation conditions were as follows:

Instrumentation	Thermo Scientific™ Dionex™ UltiMate™ 3000 BioRSLC system equipped with:		
	SR-3000 Solvent Rack (without degasser) (P/N 5035.9200)		
	LPG-3400RS Biocompatible Quaternary Rapid Separation Pump (P/N 5040.0036)		
	WPS-3000TBRS Biocompatible Rapid Separation Thermostatted Autosampler (P/N 5841.0020)		
	TCC-3000RS Rapid Separation Thermostatted Column Compartment (P/N 5730.0000)		
	VWD-3400RS Rapid Separation Variable Wavelength Detector) equipped with a micro flow cell (P/N 5074.0010)		
Column	MAbPac HIC-Butyl, 4.6 × 100 mm (P/N 088558)		
Mobile phase A	1.5 M ammonium sulfate, 50 mM sodium phosphate, pH 7.0 / 2-propanol (95:5 v/v) Dissolve 6.55 g of NaH ₂ PO ₄ •H ₂ O and 188.30 g of (NH ₄) ₂ SO ₄ in 750 mL DI water, adjust the pH to 7.0 using 50% sodium hydroxide (NaOH) solution. Bring the volume to 950 mL with DI water. Then bring the volume to 1000 mL with 2-propanol.		
Mobile phase B	50 mM sodium phosphate, pH 7.0 / 2-propanol (80:20 v/v) Dissolve 5.52 g of NaH ₂ PO ₄ •H ₂ O in 750 mL DI water, adjust the pH to 7.0 using 50% NaOH solution. Bring the volume to 800 mL with DI water. Then bring the volume to 1000 mL with 2-propanol.		
Gradient	Time (min)	%A	%B
	-5.0	100	0
	0.0	100	0
	1.0	100	0
	15.0	0	100
	20.0	0	100
Flow rate	1.0 mL/min		
Run time	20 min		
Injection volume	5 μ L		
Temperature	25 °C		
UV detector wavelength	280 nm		

Data Processing

Thermo Scientific™ Dionex™ Chromeleon™ 6.8 Chromatography Data System

Results and Discussion

The ADC mimics used in this work were conjugates between a drug mimic and mAb via the sulfhydryl group of interchain cysteine residues (Figure 1a). This procedure results in a mixture of drug-loaded antibody species with 0 to 8 drugs (Figure 1b). The stoichiometry of the drug mimic was varied in the conjugation reaction. Samples are labeled low, moderate, and high based on the average number of the drug mimic attached to the mAb. The typical separation of cysteine-conjugated ADC mimic on a MAbPac HIC-Butyl column is demonstrated in Figure 2. HIC peak identities were assigned by examining the UV spectra of each peak using a diode array detector using a literature method.⁷

Concentration of organic solvent in mobile phases and temperature are critical factors for the resolution and peak shape when analyzing protein samples in HIC. The best peak shape and separation was achieved using 5% 2-propanol in mobile phase A and 20% 2-propanol in mobile phase B at 25 °C. Without organic solvent, DAR species were not well resolved, while addition of more 2-propanol in mobile phase A or mobile phase B resulted in a very broad peak for the 8-drug form, probably due to denaturing of the heavy and light chain. All the ADC species with 0 to 8 drugs were observed with three additional peaks for the 6-drug species. This indicates there are at least four hydrophobic variants of the 6-drug species.

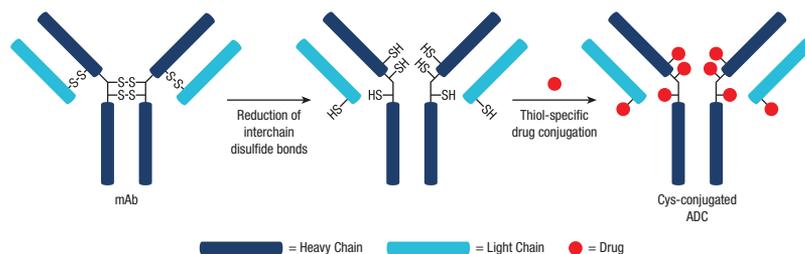


Figure 1a. Schematic representation of conjugation of drug mimic via interchain cysteine residues.

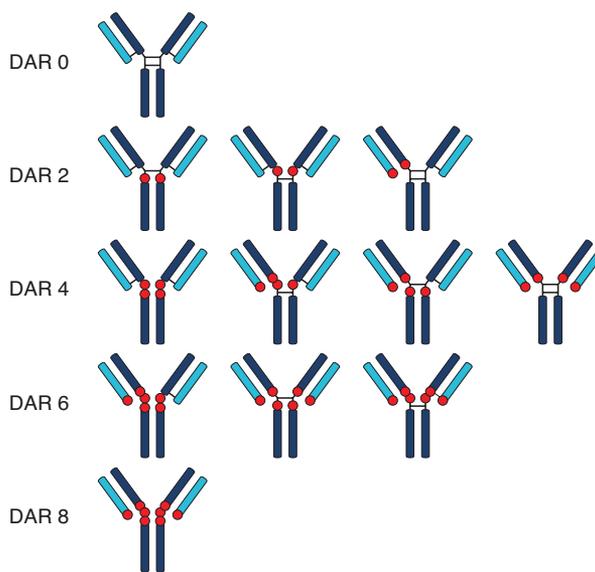


Figure 1b. Heterogeneity of cysteine-conjugated ADCs.

Column: **MAbPac HIC-Butyl**, 5 μ m
 Format: 4.6 \times 100 mm
 Mobile Phase A: 1.5 M ammonium sulfate, 50 mM sodium phosphate, pH 7.0 / 2-propanol (95:5 v/v)
 Mobile Phase B: 50 mM sodium phosphate, pH 7.0 / 2-propanol (80:20 v/v)
 Gradient:

Time (min)	%A	%B
-5.0	100	0
0.0	100	0
1.0	100	0
15.0	0	100
20.0	0	100

 Flow Rate: 1.0 mL/min
 Inj. Volume: 5 μ L
 Temp.: 25 °C
 Detection: UV (280 nm)
 Sample: Cys-conjugated ADC mimic (5 mg/mL)

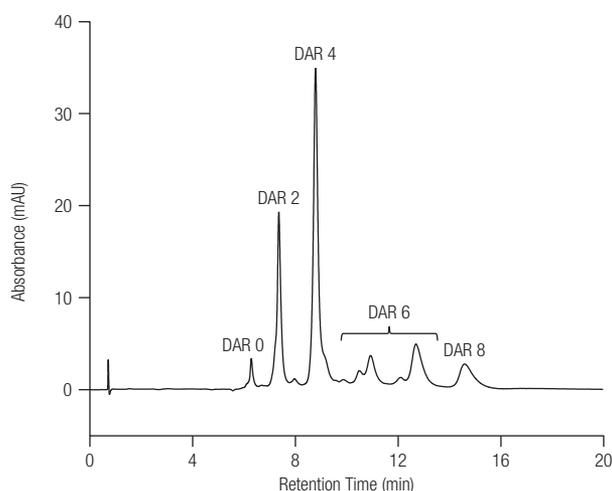


Figure 2. Separation of cysteine-conjugated ADC mimic.

Figure 3 compares the analysis of cysteine-conjugated ADC mimics with different drug loads assuming the UV absorption of the drug mimic is minimal at 280 nm. The mimic with high drug load shows no unconjugated mAb and a higher intensity of the 8-drug form while the low drug load sample contains a significant amount of unconjugated mAb, 2-drug form, and 4-drug form.

Column:	MABPac HIC-Butyl , 5 μ m		
Dimension:	4.6 \times 100 mm		
Mobile Phase A:	1.5 M ammonium sulfate, 50 mM sodium phosphate, pH 7.0 / 2-propanol (95:5 v/v)		
Mobile Phase B:	50 mM sodium phosphate, pH 7.0 / 2-propanol (80:20 v/v)		
Gradient:	Time (min)	%A	%B
	-5.0	100	0
	0.0	100	0
	1.0	100	0
	15.0	0	100
	20.0	0	100
Flow Rate:	1.0 mL/min		
Inj. Volume:	5 μ L		
Temp.:	25 $^{\circ}$ C		
Detection:	UV (280 nm)		
Sample:	a. Unconjugated mAb (5 mg/mL)		
	b. Cys-conjugated ADC mimic (low load; 5 mg/mL)		
	c. Cys-conjugated ADC mimic (moderate load; 5 mg/mL)		
	d. Cys-conjugated ADC mimic (high load; 5 mg/mL)		

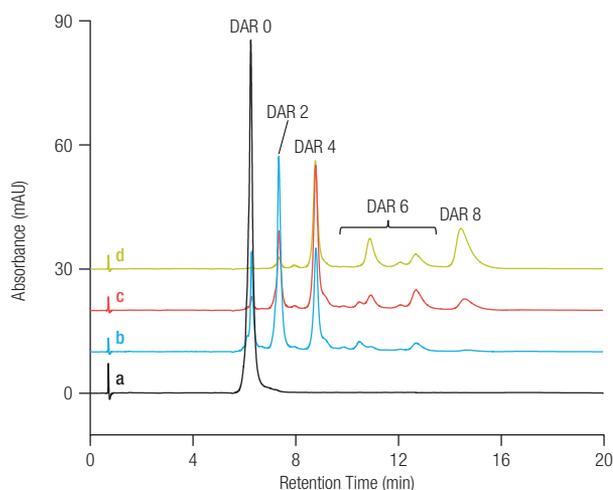


Figure 3. Comparison of cysteine-conjugated ADC mimics with different drug loads

Conclusion

- The MABPac HIC-Butyl column provides high resolution for unconjugated mAbs and cysteine-linked ADCs with different drug-to-antibody ratios (DARs).
- The drug load distribution can be monitored using the MABPac HIC-Butyl column.

References

- 1 Leavy, O. Therapeutic antibodies: past, present and future, *Nature Review Immunology*, **2010**, *10*, 297.
- 2 Sievers, E.L. *et al.* Antibody-Drug Conjugates in Cancer Therapy, *Annu. Rev. Med.*, **2013**, *64*, 15-29.
- 3 Thayer, A.M. Building antibody-drug conjugates, *C&EN*, **2014**, *92*, 13-21.
- 4 Hamblett, K.J. *et al.* Effects of Drug Loading on the Antitumor Activity of a Monoclonal Antibody Drug Conjugate, *Clin. Cancer Res.*, **2004**, *10*, 7063-70.
- 5 Valliere-Douglass, J. *et al.* Separation and characterization of an IgG2 antibody containing a cyclic imide in CDR1 of light chain by hydrophobic interaction chromatography and mass spectrometry, *Anal. Chem.*, **2008**, *80*, 3168-3174.
- 6 Valliere-Douglass, J. *et al.* Separation of populations of antibody variants by fine tuning of hydrophobic-interaction chromatography operating conditions, *J. Chromatogr. A*, **2008**, *1214*, 81-89.
- 7 Ouyang, J. Drug-to-antibody ratio (DAR) and drug load distribution by hydrophobic interaction chromatography and reversed phase high-performance liquid chromatography, *Methods in Molecular Biology*, **2013**, *1045*, 275-283.

For Research Use Only

thermoscientific.com/biolc

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other 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.



USA and Canada +1 800 332 3331
Australia 1300 735 292 (free call domestic)
China 800 810 5118 (free call domestic)
 400 650 5118
France +33 (0)1 60 92 48 34
Germany +49 (0) 2423 9431 20 or 21
India +91 22 6742 9494
 +91 27 1766 2352

Japan 0120 753 670 (free call domestic)
 0120 753 671 (fax)
Korea +82 2 3420 8600
United Kingdom +44 (0) 1928 534 110
New Zealand 0800 933 966 (free call domestic)
Singapore +65 6289 1190
All Other Enquiries +44 (0) 1928 534 050

Technical Support

For advice and support, please visit our website:
www.thermoscientific.com/chromexpert

Thermo
 SCIENTIFIC

A Thermo Fisher Scientific Brand