



## Introduction

Monoclonal antibodies (mAbs) represent a major category of therapeutic proteins. A number of bioanalytical tools are required to monitor the antibody heterogeneity. Liquid chromatography/mass spectrometry (LC/MS) is routine technology nowadays applied for characterization of these biomolecules. Often the preparation and storage reagents of mAbs contains nonvolatile salts, detergents, and solubilizing agents. The presence of these reagents causes adverse effects in mass spectrometry, suppressing ionization which in turn limits LC/MS application. Hence it is very important to remove these salts and detergents prior to MS analysis. In the present study, a new desalting cartridge is evaluated for 1D and 2D-LC/MS based protein applications. The reverse phase LC method for desalting was developed using a polymeric based desalting material. For effective desalting performance, (a) salt containing samples (1D), and (b) two multidimensional separation experiments (IEX → RP and affinity chromatography → RP) were tested. The desalting approach described in this work demonstrates that coupling of these desalting columns to mass spectrometry improves the protein spectra from salt and detergent contaminated solutions.

## Experimental

Monoclonal antibodies mAb1, and mAb2 were purchased from a local pharmacy and stored according to the manufacturers' instructions. All solvents used were LC grade.

| First-dimensional pump (IEX)        |  |
|-------------------------------------|--|
| Solvent A                           | Water  |
| Solvent B                           | NaCl (850.0 mM)  |
| Solvent C                           | NaH <sub>2</sub> PO <sub>4</sub> (41.0 mM)   |
| Solvent D                           | Na <sub>2</sub> HPO <sub>4</sub> (55.0 mM)   |
| Flow rate                           | 0.75 mL/min  |
| Gradient                            | 0 minutes: 30.3 %A, 0.0 %B, 59.6 %C, 10.1 %D<br>2 minutes: 26.0 %A, 5.0 %B, 56.9 %C, 12.1 %D<br>8 minutes: 21.5 %A, 10.0 %B, 54.9 %C, 13.6 %D<br>20 minutes: 13.3 %A, 19.0 %B, 51.9 %C, 15.8 %D<br>35 minutes: 30.3 %A, 0.0 %B, 59.6 %C, 10.1 %D   |
| Post time                           | 10 minutes   |
| First-dimensional pump (Protein A)  |  |
| Solvent A                           | 20 mM sodium phosphate buffer, pH 7.4  |
| Solvent B                           | Acetic acid (500.0 mM)   |
| Flow rate                           | 1 mL/min   |
| Gradient                            | 0 to 0.5 minutes: 0%B (binding)<br>0.6 to 1.7 minutes: 100%B (elution)<br>1.8 to 3.5 minutes: 0%B (regeneration)   |
| Second-dimensional pump             |  |
| Solvent A                           | 0.1% FA  |
| Solvent B                           | 0.1% FA in ACN   |
| Flow rate                           | 0.4 mL/min   |
| Gradient                            | 0 minutes - 5 %B, 0.5 minutes - 5 %B, 3 minutes - 80 %B, 4 minutes - 80 %B, 4.1 minutes - 5 %B, 6 minutes - 5 %B   |
| 2D gradient stop time               |  |
| 2D cycle time                       |  |
| Autosampler                         |  |
| Injection volume                    | 5 µL   |
| Sample temperature                  | 5 °C   |
| Columns (IEX → RP) (Protein A → RP) |  |
| 1st - dimension column              | Agilent Bio MAb NP5, 4.6 x 250 mm, PEEK  |
| 1st - dimension column              | Agilent Bio-Monolith Column Protein A  |
| 2nd-dimension column                | Agilent AdvanceBio RP desalting cartridges , 2.1 x 12.5 mm.  |
| Thermostatted column compartment    |  |
| 1st - dimension column              | Room temperature   |
| 2nd-dimension column                | Room temperature   |
| Multiple heart-cutting              |  |
| Operation mode                      | <ul style="list-style-type: none"> <li>Time-based multiple heart-cutting was performed with the heart-cuts (2D time segments) set according to the first-dimension retention times.</li> <li>Heart-cutting of impurities was conducted with a sampling time of 0.04 minutes (loop filling of &gt; 200 %).</li> </ul> |
| Detection                           |  |
| 1st - dimension                     | UV Wavelength 280 nm/4 nm, Ref.: 360 nm/100 nm   |
| 2nd - dimension                     | MS parameters  |
|                                     | Gas temperature: 350 °C  |
|                                     | Sheath gas temperature: 400 °C   |
|                                     | Gas flow: 8 L/min  |
|                                     | Sheath gas flow: 11 L/min  |
|                                     | Nebulizer: 35 psi  |
|                                     | Vcap: 5,000 V  |
|                                     | Nozzle: 1,000 V  |
|                                     | Fragmentor: 200  |
| Time (min)                          | LC stream to   |
| 0                                   | Waste  |
| 10.8                                | MS   |
| 14.6                                | Waste  |

## Results and Discussion

### Instrument and column configuration

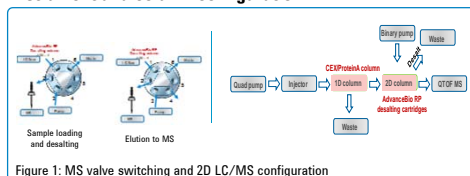


Figure 1: MS valve switching and 2D LC/MS configuration

### High quality MS with AdvanceBio RP desalting cartridges

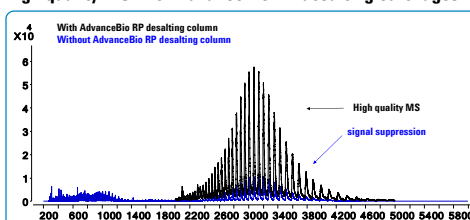


Figure 2: Desalting of mAb1 samples using 2D LC/MS. 2<sup>nd</sup> dimensional mass spectrum profiles with and without AdvanceBio RP desalting cartridges. Use of desalt column provides improvement in signal intensity (6X greater MS signal).

### On-line desalting 2D LC/MS (IEX → RP): mAb1

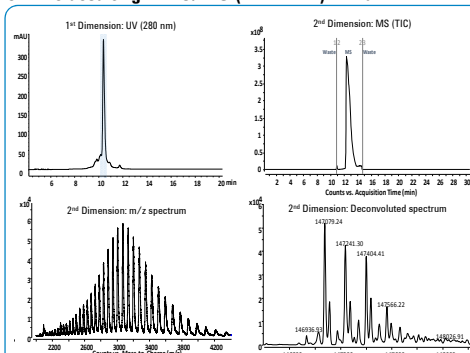


Figure 3: 2D LC/MS of mAb1 profiles. 1D column: IEX; 2D column: Desalt cartridges. The selected main peak in the 1D chromatogram is captured and transferred to the AdvanceBio desalting cartridges that effectively remove salts and provide high quality MS results.

### AdvanceBio RP desalting cartridges: reproducibility

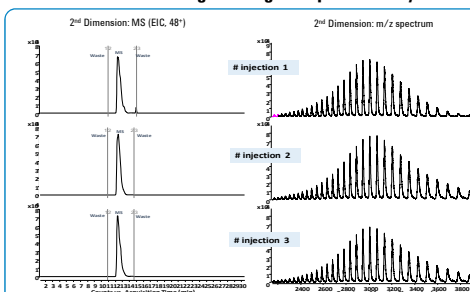


Figure 4: 2D LC/MS of mAb1 profiles. 1D column: IEX; 2D column: Desalting cartridges. 2<sup>nd</sup> dimension EIC (left) and m/z (right) profiles. Injection to injection repeatability shows consistent performance of AdvanceBio RP desalting cartridges.

## Results and Discussion

### Desalting of reduced mAb with different concentrations of NaCl (1D experiments)

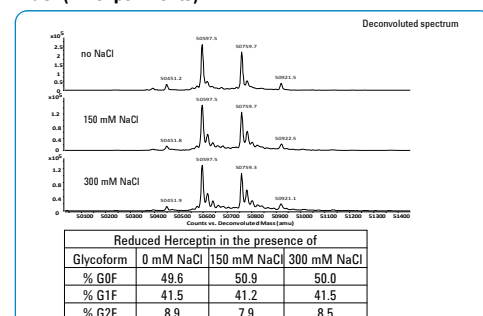


Figure 5: Deconvoluted spectrum of Herceptin heavy chain. NaCl containing mAb samples were analyzed on the AdvanceBio RP desalting cartridges following reduction using DTT. The signal decreases slightly when samples contain more salt, but relative quantification is similar between samples in the presence or absence of NaCl.

### On-line desalting 2D LC/MS (IEX → RP): mAb2

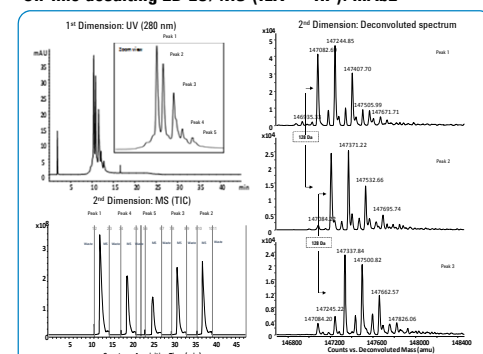


Figure 6: 2D LC/MS of mAb2 profiles. 1D column: IEX; 2D column: Desalt cartridges. Selected peaks (multiple heart-cutting) in the 1D chromatogram was captured and transferred to the AdvanceBio desalting cartridges that effectively removed salts and provided high quality MS results. The 3 main peaks (peaks 1,2,3) corresponds to C-terminal lysine variants.

### On-line desalting 2D LC/MS (Protein A → RP): mAb1

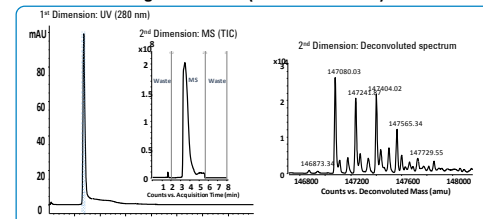


Figure 7: 2D LC/MS of mAb1 profiles. 1D column: Protein A; 2D column: Desalt cartridges. Selected peaks in the 1D chromatogram was captured and transferred to the AdvanceBio desalt cartridges that effectively removed salts and provided high quality MS results.

## Conclusions

- AdvanceBio desalting cartridges effectively remove salts for better MS sensitivity.
- AdvanceBio desalting cartridges provide reproducible MS results.
- AdvanceBio desalting cartridges allow automated 2D LC in Multiple Heart-Cutting 2D-LC/MS methods for characterization of mAbs.