



# Multi Affinity Removal Column, Human-7

Part Nos. 5188-6409 ( $4.6 \times 50$  mm)

5188-6410 ( $4.6 \times 100$  mm)

5188-6411 ( $10 \times 100$  mm)

**For depletion of seven high-abundance  
proteins from human plasma samples**

## Instructions

Version A1, December 2018

**For Research Use Only. Not for use in diagnostic  
procedures.**



**Agilent Technologies**

## Notices

© Agilent Technologies, Inc. 2016, 2018

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

### Manual Part Number

5990-7228

### Edition

Version A1, December 2018

Printed in USA

Agilent Technologies, Inc.  
5301 Stevens Creek Blvd  
Santa Clara, CA 95051 USA

### Technical Support

#### For US and Canada

Call (800) 227-9770 (option 3,4,4)

Or send an e-mail to:  
[techservices@agilent.com](mailto:techservices@agilent.com)

#### For Europe, Middle East, Africa, and India

Call 00800 345 600 (toll free) or  
+49 69 8679 7730

Or send an e-mail to:  
[genomics\\_tech\\_europe@agilent.com](mailto:genomics_tech_europe@agilent.com)

#### For all other regions

Agilent's world-wide sales and support center telephone numbers can be obtained at [www.agilent.com](http://www.agilent.com) under Contact Us.

Or send an e-mail to:  
[genomics@agilent.com](mailto:genomics@agilent.com)

## Warranty

**The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.**

## Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

## Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## Safety Notices

### CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

### WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

## In this Guide...

This document describes how to use the Multi Affinity Removal Column, Human-7 to chromatographically remove seven interfering high-abundance proteins from human plasma prior to LC/MS or electrophoretic analysis of the samples.

### **1 Before You Begin**

This chapter contains information (such as required reagents and equipment) that you should read and understand before you start an experiment.

### **2 Instructions**

This chapter describes the protocol for chromatographic removal of the targeted proteins from human plasma samples and includes troubleshooting information.

### **3 Reference**

This chapter contains reference information including column specifications and a list of related products.

## What's New in Version A1

- Updates to description of needles included with reagent kits (see [Table 8](#) on page 26)
- Correction to part numbers for spin cartridge Starter Reagent Kit, p/n 5188-5254, and mRP-C18 desalting column p/n 5188-5231 (see [Table 8](#) on page 26)
- Updates to Technical Support contact information (see [page 2](#))

# Contents

|          |                                                  |           |
|----------|--------------------------------------------------|-----------|
| <b>1</b> | <b>Before You Begin</b>                          | <b>6</b>  |
|          | Safety Considerations                            | 7         |
|          | Materials Required                               | 7         |
|          | Storage Conditions                               | 8         |
|          | Overview                                         | 8         |
| <b>2</b> | <b>Instructions</b>                              | <b>10</b> |
|          | Protocol for 4.6 x 50 mm column                  | 11        |
|          | Step 1. Set up the column (4.6 x 50 mm column)   | 11        |
|          | Step 2. Prepare the sample (4.6 x 50 mm column)  | 12        |
|          | Step 3. Run the column (4.6 x 50 mm column)      | 12        |
|          | Protocol for 4.6 x 100 mm column                 | 14        |
|          | Step 1. Set up the column (4.6 x 100 mm column)  | 14        |
|          | Step 2. Prepare the sample (4.6 x 100 mm column) | 15        |
|          | Step 3. Run the column (4.6 x 100 mm column)     | 15        |
|          | Protocol for 10 x 100 mm column                  | 17        |
|          | Step 1. Set up the column (10 x 100 mm column)   | 17        |
|          | Step 2. Prepare the sample (10 x 100 mm column)  | 18        |
|          | Step 3. Run the column (10 x 100 mm column)      | 18        |
|          | Guidelines for flow-through fraction analysis    | 19        |
|          | Recommendations                                  | 20        |
|          | Troubleshooting                                  | 22        |
| <b>3</b> | <b>Reference</b>                                 | <b>23</b> |
|          | Column Specifications                            | 24        |
|          | Related Agilent Products                         | 25        |



# 1

## Before You Begin

Safety Considerations 7

Materials Required 7

Storage Conditions 8

Overview 8

Make sure you read and understand the information in this chapter and have the necessary equipment and reagents listed before you start an experiment.



## Safety Considerations

When preparing biological samples using Agilent Multiple Affinity Removal Columns, follow general guidelines for handling biological materials and wear protective eyewear and gloves.

## Materials Required

The Agilent Multiple Affinity Removal Columns and accessories (purchased separately) used in this protocol are shown in [Table 1](#).

**Table 1** Multiple Affinity Removal Column and Accessories

| Part number | Product name                                                | Description                                                                                                                 |
|-------------|-------------------------------------------------------------|-----------------------------------------------------------------------------------------------------------------------------|
| 5188-6409   | Multi Affinity Removal Column Human-7, 4.6 x 50 mm, 1 each  |                                                                                                                             |
| 5188-6410   | Multi Affinity Removal Column Human-7, 4.6 x 100 mm, 1 each | LC column used to remove albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin, and fibrinogen from human plasma samples |
| 5188-6411   | Multi Affinity Removal Column Human-7, 10 x 100 mm, 1 each  |                                                                                                                             |
| 5185-5987   | Buffer A, 1 L                                               | Ready-to-use, optimized buffer for loading, washing, and equilibrating column                                               |
| 5185-5988   | Buffer B, 1 L                                               | Ready-to-use, optimized buffer for elution of bound proteins from column                                                    |
| 5185-5990   | Spin filters, 0.22 µm, 1 pack of 25                         | For sample cleanup before loading column                                                                                    |
| 5185-5991   | Concentrators, 5 kDa MWCO, 1 pack of 25                     | For concentrating flow-through fractions                                                                                    |
| 5185-5989   | Human serum albumin                                         | Dilute standard for checking column capacity (optional)                                                                     |
| 5190-7995   | MARS Column 2 µm Replacement Frit, 2 each                   | One set of 2 frit assemblies for replacement of clogged inlet and outlet column frits                                       |
| 5185-5986   | Starter Reagent Kit                                         | Buffer A: 2 x 1 L<br>Buffer B: 1 L<br>Spin filters 0.22 µm: 2 packs of 25<br>Protein concentrators: 1 pack of 25            |

## CAUTION

Do not expose columns to organic solvents (like alcohols, acetonitrile, etc.), strong oxidizers, acids, reducing agents, or other protein denaturing agents.

Before attaching the column, purge the LC system and run two method blank injections according to protocol to ensure all lines and sample loops are free of organic solvents.

For LC systems shared with other chemical applications, be sure to first purge the LC system, including the sample loop, with isopropyl alcohol, and then extensively with water (approximately 1 hour). After purging, proceed with protocol.

## Storage Conditions

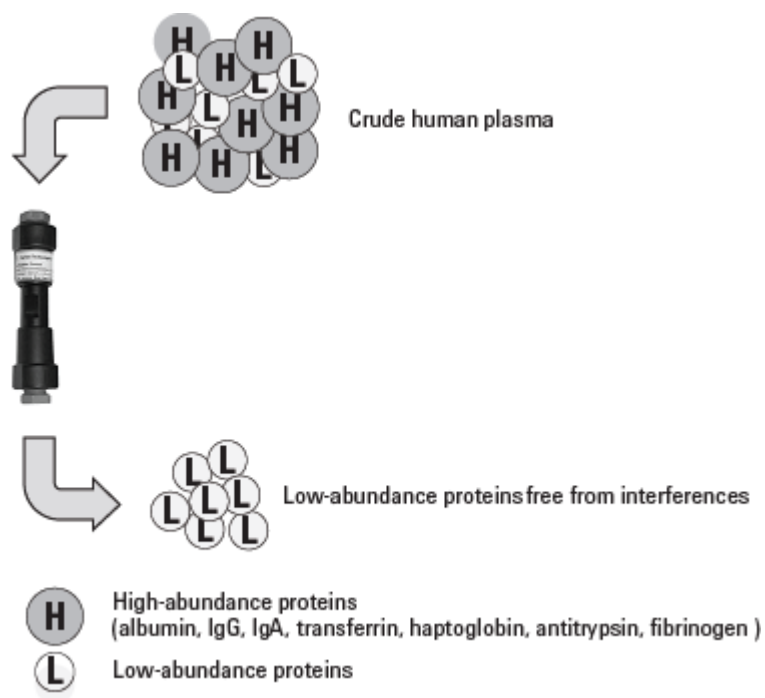
Upon its receipt and when you are not using it, store the column with the end-caps tightly sealed at 2°C to 8°C (35°F to 46°F). **Do not freeze the column.**

## Overview

The Agilent Multiple Affinity Removal System comprises a family of immunodepletion products based on affinity interactions and optimized buffers for sample loading, washing, eluting, and regenerating. This column is specifically designed to remove seven high-abundance proteins from human plasma samples. This technology enables removal of albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin, and fibrinogen with a single device. The targeted high-abundance proteins are simultaneously removed when crude biological samples are passed through the column. Selective immunodepletion provides an enriched pool of low-abundance proteins for downstream proteomics analysis, as depicted in [Figure 1](#) on page 9.

Specific removal of the seven high-abundance proteins depletes approximately 88–92% of total protein mass from human plasma, facilitating study of the low-abundance proteins in the flow-through fractions. Removal of high-abundance proteins enables improved resolution and dynamic range for one-dimensional gel electrophoresis (1DGE), two-dimensional gel electrophoresis (2DGE) and liquid chromatography/mass spectrometry (LC/MS). The collected flow-through fractions may need to be concentrated dependent upon the downstream applications.





**Figure 1** The Multiple Affinity Removal System.



## 2 Instructions

### Protocol for 4.6 x 50 mm column 11

- Step 1. Set up the column (4.6 × 50 mm column) 11
- Step 2. Prepare the sample (4.6 × 50 mm column) 12
- Step 3. Run the column (4.6 × 50 mm column) 12

### Protocol for 4.6 x 100 mm column 14

- Step 1. Set up the column (4.6 × 100 mm column) 14
- Step 2. Prepare the sample (4.6 × 100 mm column) 15
- Step 3. Run the column (4.6 × 100 mm column) 15

### Protocol for 10 x 100 mm column 17

- Step 1. Set up the column (10 × 100 mm column) 17
- Step 2. Prepare the sample (10 × 100 mm column) 18
- Step 3. Run the column (10 × 100 mm column) 18

### Guidelines for flow-through fraction analysis 19

### Recommendations 20

### Troubleshooting 22



## Protocol for 4.6 x 50 mm column

Column capacity: 30–35 µL human plasma

### Step 1. Set up the column (4.6 × 50 mm column)

- 1 Set up Buffer A and Buffer B as the only mobile phases.
- 2 Purge lines with Buffer A and Buffer B at a flow rate of 1.0 mL/min for 10 min **without a column**.
- 3 Set up LC timetable as specified in [Table 2](#).
- 4 Run two method blanks by injecting 100 µL of Buffer A **without a column**.
- 5 Ensure that you are using the proper sample loop size in the autosampler, and that the sample loop has been flushed with Buffer A.
- 6 Attach the column and equilibrate it in Buffer A for 4 min at a flow rate of 1 mL/min at room temperature.

**Table 2** LC method for 4.6 x 50 mm column \*

| Step | Time (min) | %B     | Flow Rate (mL/min) | Max. Pressure (bar) |
|------|------------|--------|--------------------|---------------------|
| 1    | 0.00       | 0.00   | 0.250              | 120                 |
| 2    | 9.00       | 0.00   | 0.250              | 120                 |
| 3    | 9.01       | 100.00 | 1.000              | 120                 |
| 4    | 12.50      | 100.00 | 1.000              | 120                 |
| 5    | 12.60      | 0.00   | 1.000              | 120                 |
| 6    | 20.00      | 0.00   | 1.000              | 120                 |

\* Solvent A: Buffer A  
Solvent B: Buffer B  
Detection wavelength: 280 nm

## Step 2. Prepare the sample (4.6 × 50 mm column)

Before you begin, consult the Certificate of Analysis for your column to verify the column capacity.

- 1 Dilute the plasma sample four-fold with Buffer A. For example, if the recommended column loading capacity on the Certificate of Analysis is 35 µL of plasma, dilute 35 µL of plasma with 105 µL Buffer A for a final volume of 140 µL.

Addition of protease inhibitors in Buffer A for sample dilution helps prevent protein degradation.

### NOTE

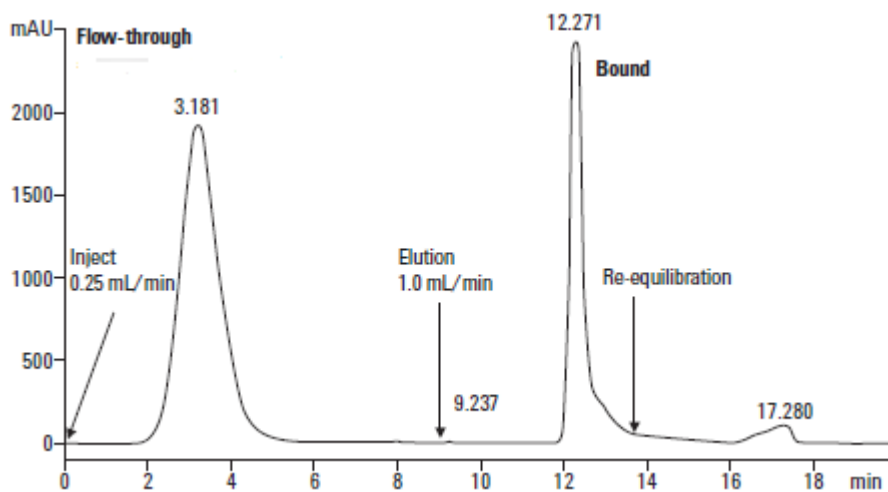
The protocol may be applied to other human biological fluids like serum and CSF with necessary adjustments in sample volume based on albumin concentration.

- 2 Remove particulates with a 0.22 µm spin filter, spinning for 1 min at 16,000 × *g*.

## Step 3. Run the column (4.6 × 50 mm column)

Use the LC timetable (Table 2 on page 11) to complete the following steps.

- 1 Inject the diluted plasma sample at a flow rate of 0.25 mL/min.
- 2 Collect the flow-through fraction (like that which appears from 1.5 to 4.5 min in the typical chromatogram shown in Figure 2 on page 13). Store collected fractions at -20 °C if not analyzed immediately.
- 3 Elute bound proteins from the column with Buffer B (elution buffer) at a flow rate of 1 mL/min for 3.5 min.
- 4 Regenerate column by equilibrating it with Buffer A for an additional 7.4 min at a flow rate of 1 mL/min.
- 5 After equilibration with Buffer A, store the column with ends capped at 2°C to 8°C (35°F to 46°F). **Do not freeze the column.**
- 6 Analyze the flow-through fraction using the guidelines on page 19.

Step 3. Run the column ( $4.6 \times 50$  mm column)

**Figure 2** Representative chromatogram for  $4.6 \times 50$  mm column.

## Protocol for 4.6 x 100 mm column

Column capacity: 60–70 µL human plasma

### Step 1. Set up the column (4.6 × 100 mm column)

- 1 Set up Buffer A and Buffer B as the only mobile phases.
- 2 Purge lines with Buffer A and Buffer B at a flow rate of 1.0 mL/min for 10 min **without a column**.
- 3 Set up LC timetable as specified in [Table 3](#).
- 4 Run two method blanks by injecting 200 µL of Buffer A **without a column**.
- 5 Ensure that you are using the proper sample loop size in the autosampler, and that the sample loop has been flushed with Buffer A.
- 6 Attach the column and equilibrate it in Buffer A for 4 min at a flow rate of 1 mL/min at room temperature.

**Table 3** LC method for 4.6 x 100 mm column\*

| Step | Time (min) | %B     | Flow Rate (mL/min) | Max. Pressure (bar) |
|------|------------|--------|--------------------|---------------------|
| 1    | 0.00       | 0.00   | 0.50               | 120                 |
| 2    | 10.00      | 0.00   | 0.50               | 120                 |
| 3    | 10.01      | 100.00 | 1.000              | 120                 |
| 4    | 17.00      | 100.00 | 1.000              | 120                 |
| 5    | 17.01      | 0.00   | 1.000              | 120                 |
| 6    | 28.00      | 0.00   | 1.000              | 120                 |

\* Solvent A: Buffer A  
 Solvent B: Buffer B  
 Detection wavelength: 280 nm

## Step 2. Prepare the sample (4.6 × 100 mm column)

Before you begin, consult the Certificate of Analysis for your column to verify the column capacity.

- 1 Dilute the plasma sample four-fold with Buffer A. For example, if the recommended column loading capacity on the Certificate of Analysis is 70 µL of plasma, dilute 70 µL of plasma with 210 µL Buffer A for a final volume of 280 µL.

Addition of protease inhibitors in Buffer A for sample dilution helps prevent protein degradation.

### NOTE

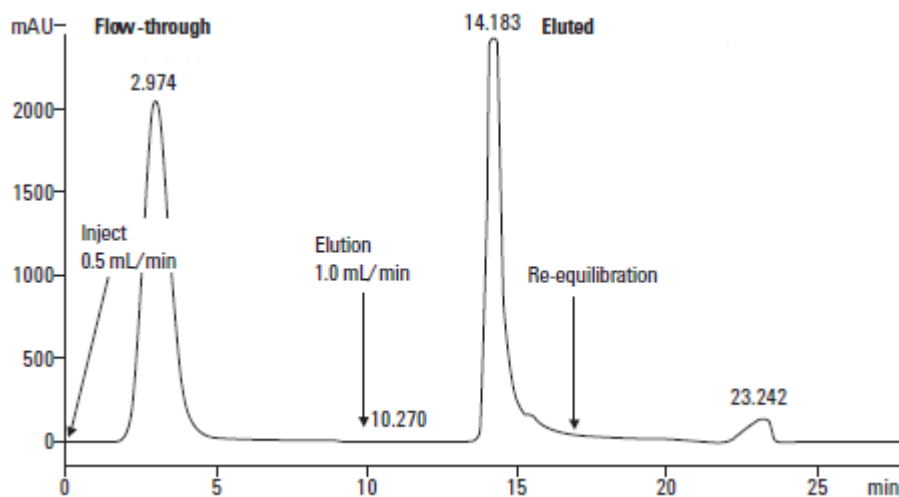
The protocol may be applied to other human biological fluids like serum and CSF with necessary adjustments in sample volume based on albumin concentration.

- 2 Remove particulates with a 0.22 µm spin filter, spinning for 1 min at 16,000 × g.

## Step 3. Run the column (4.6 × 100 mm column)

Use the LC timetable (Table 3 on page 14) to complete the following steps.

- 1 Inject the diluted plasma sample at a flow rate of 0.5 mL/min.
- 2 Collect the flow-through fraction (like that which appears from 2.5 to 6 min in the typical chromatogram shown in Figure 3 on page 16). Store collected fractions at -20 °C if not analyzed immediately.
- 3 Elute bound proteins from the column with Buffer B (elution buffer) at a flow rate of 1 mL/min for 7.0 min.
- 4 Regenerate column by equilibrating it with Buffer A for an additional 11.0 min at a flow rate of 1 mL/min.
- 5 After equilibration with Buffer A, store the column with ends capped at 2°C to 8°C (35°F to 46°F). **Do not freeze the column.**
- 6 Analyze the flow-through fraction using the guidelines on page 19.

Step 3. Run the column ( $4.6 \times 100$  mm column)

**Figure 3** Representative chromatogram for  $4.6 \times 100$  mm column.



## Protocol for 10 × 100 mm column

Column capacity: 250–300 µL of human plasma

### Step 1. Set up the column (10 × 100 mm column)

- 1 Set up Buffer A and Buffer B as the only mobile phases.
- 2 Purge lines with Buffer A and Buffer B at a flow rate of 1.0 mL/min for 10 min **without a column**.
- 3 Set up LC timetable as specified in [Table 4](#).
- 4 Run two method blanks by injecting 200 µL of Buffer A **without a column**.
- 5 Ensure that you are using the proper sample loop size in the autosampler, and that the sample loop has been flushed with Buffer A.
- 6 Attach the column and equilibrate it in Buffer A for 4 min at a flow rate of 1 mL/min at room temperature.

**Table 4** LC method for 10 x 100 mm column \*

| Step | Time (min) | %B     | Flow Rate (mL/min) | Max. Pressure (bar) |
|------|------------|--------|--------------------|---------------------|
| 1    | 0.00       | 0.00   | 0.500              | 120                 |
| 2    | 20.00      | 0.00   | 0.500              | 120                 |
| 3    | 20.01      | 100.00 | 3.000              | 120                 |
| 4    | 27.00      | 100.00 | 3.000              | 120                 |
| 5    | 27.01      | 0.00   | 3.000              | 120                 |
| 6    | 35.00      | 0.00   | 3.000              | 120                 |

\* Solvent A: Buffer A  
 Solvent B: Buffer B  
 Detection wavelength: 280 nm

## Step 2. Prepare the sample (10 × 100 mm column)

Before you begin, consult the Certificate of Analysis for your column to verify the column capacity.

- 1 Dilute the plasma sample four-fold with Buffer A. For example, if the recommended column loading capacity on the Certificate of Analysis is 300 µL of plasma, dilute 300 µL of plasma with 900 µL Buffer A for a final volume of 1200 µL.

Addition of protease inhibitors in Buffer A for sample dilution helps prevent protein degradation.

### NOTE

The protocol may be applied to other human biological fluids like serum and CSF with necessary adjustments in sample volume based on albumin concentration.

- 2 Remove particulates with a 0.22 µm spin filter, spinning for 1 min at 16,000 × g.

## Step 3. Run the column (10 × 100 mm column)

Use the LC timetable (Table 4 on page 17) to complete the following steps.

- 1 Inject the diluted plasma sample at a flow rate of 0.5 mL/min.
- 2 Collect the flow-through fraction, appearing from 9 to 14 min using the LC method in Table 4 on page 17. Store collected fractions at -20 °C if not analyzed immediately.
- 3 Elute bound proteins from the column with Buffer B (elution buffer) at a flow rate of 3 mL/min for 7.0 min.
- 4 Regenerate column by equilibrating it with Buffer A for an additional 8.0 min at a flow rate of 3 mL/min.
- 5 After equilibration with Buffer A, store the column with ends capped at 2°C to 8°C (35°F to 46°F). **Do not freeze the column.**
- 6 Analyze the flow-through fraction using the guidelines on page 19.

## Guidelines for flow-through fraction analysis

Analyze the flow-through fraction, containing the low-abundance proteins, to verify removal of the targeted high-abundance proteins using the guidelines below:

- For 1D-SDS-PAGE, an aliquot of the flow-through fraction may be used directly.
- For IEF, 2D-GE, and MS analysis of the flow-through fraction, it is necessary to do buffer exchange or desalt to an appropriate buffer. The 5 kDa MWCO spin concentrators (part number 5185-5991) may be used for buffer exchange and concentration. Alternatively, the Agilent mRP-C18 column (part number 5188-5231) may be used for automated desalting and concentration.

## Recommendations

- **Sample dilution using Buffer A**

Do not load crude plasma or other biological samples directly onto the column. Follow instructions for plasma dilution with Buffer A in the protocol for your specific column size.

- **Preventing protein degradation**

Addition of protease inhibitors to Buffer A for sample dilution helps prevent protein degradation.

- **Sample cleanup**

Human plasma may contain particulate materials that can be removed by a quick spin using a 0.22- $\mu$ m spin filter.

- **Variation in column capacity for different samples**

Concentrations of the proteins targeted for depletion can vary among individual plasma samples and in different types of biological samples. Thus column capacity for samples may differ and you may need to adjust the loading volume for a particular plasma sample.

For any samples that require adjustment of the plasma load volume, adhere to the instruction to dilute the samples four-fold with Buffer A; do not vary the proportion of plasma and Buffer A in the diluted sample.

- **Column performance**

Agilent Multiple Affinity Removal Columns should perform reproducibly for greater than 200 runs when handled using the recommended procedures. Buffers A and B are optimized to support column performance and longevity. We cannot guarantee column performance if other buffers are used.

Do not expose columns to organic solvents (like alcohols, acetonitrile, etc.), strong oxidizers, acids, reducing agents, or other protein-denaturing agents.

- **Column storage**

To minimize loss in capacity, equilibrate the column with Buffer A. Cap the ends and store at 2°C to 8°C (35°F to 46°F). **Do not freeze the column.**

- **Analysis of flow-through fractions**

Buffer exchange to an appropriate buffer is recommended for high salt-sensitive applications such as IEF or MS. For 1D-SDS-PAGE, you can load flow-through fractions in Buffer A directly.

- **Fractionation, desalting, or concentration of flow-through fraction**  
Agilent mRP-C18 column (part number 5188-5231) is recommended for fractionation, desalting, or concentration of flow-through fractions with extremely high protein recoveries. Alternatively, spin concentrators with 5 kDa MWCO (part number 5185-5991) can be used to concentrate proteins before analysis.
- **Lyophilization of flow-through fraction**  
If lyophilization of the flow-through fraction (containing the low abundance proteins) is required after recovery from the column, first do buffer exchange to a volatile buffer (such as ammonium bicarbonate). This is recommended due to the high salt concentration of the Buffer A solvent in the flow-through fraction.
- **Bound fraction analysis**  
If you wish to analyze the bound fraction, first do buffer exchange to phosphate-buffered saline (PBS) or to another buffer compatible with your analysis. Buffer B contains compounds that may interfere with some protein assays.

## Troubleshooting

Review the following information for troubleshooting your experiments.

**Table 5** Troubleshooting suggestions

| Problem                                                                 | Cause                                                                                           | Solution                                                                                                                                               |
|-------------------------------------------------------------------------|-------------------------------------------------------------------------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>High backpressure</b>                                                |                                                                                                 |                                                                                                                                                        |
| <b>Distorted peak shape</b>                                             | Clogged inlet frits                                                                             | Remove particulates from samples with a spin filter before loading and replace plugged frits (part number 5190-7995) on both ends of the column.       |
| <b>Diminished column lifetime</b>                                       |                                                                                                 |                                                                                                                                                        |
| <b>No bound fraction peak</b>                                           | Salt concentration in diluted sample is not appropriate for affinity binding to column          | Sample must be diluted 1:4 in Buffer A to achieve the correct conditions for affinity binding. Follow the recommended sample preparation instructions. |
|                                                                         | Insufficient time for exposure of column to Buffer B (elution buffer)                           | Check LC timetable to ensure enough exposure time to Buffer B for complete removal of bound proteins.                                                  |
| <b>Breakthrough of high-abundance proteins in flow-through fraction</b> | Column plasma capacity exceeded                                                                 | Reduce plasma load per sample.                                                                                                                         |
|                                                                         | Plasma protein levels may be unusually high                                                     | Reduce plasma load per sample.                                                                                                                         |
| <b>Abnormal peak height*</b>                                            | Column may not have been regenerated well enough from previous runs, resulting in lost capacity | Elute bound proteins with Buffer B for an additional 3 min and re-equilibrate the column with Buffer A.                                                |
|                                                                         | Biological growth in the Buffer A reservoir                                                     | Replace with fresh Buffer A.                                                                                                                           |

\* Approximately 90% of plasma proteins will be removed as the bound fraction. The peak height of the bound fraction is expected to be greater than that of the flow-through fraction. If the order is reversed, consider the possible causes in the table above.



### 3 Reference

Column Specifications [24](#)

Related Agilent Products [25](#)

This chapter contains reference information.



## Column Specifications

**Table 6** Column specifications

| Parameter               | Description                                                                                        |                              |                             |
|-------------------------|----------------------------------------------------------------------------------------------------|------------------------------|-----------------------------|
| Type                    | Affinity depletion column                                                                          |                              |                             |
| Part number             | 5188-6409                                                                                          | 5188-6410                    | 5188-6411                   |
| Size                    | 4.6 mm × 50 mm<br>(0.83 mL)                                                                        | 4.6 mm × 100 mm<br>(1.66 mL) | 10 mm × 100 mm<br>(7.85 mL) |
| Column capacity*        | 30–35 µL human<br>plasma                                                                           | 60–70 µL human<br>plasma     | 250–300 µL human<br>plasma  |
| Column body material    | PEEK (polyetheretherketone)                                                                        |                              |                             |
| End-fitting material    | PEEK with 2-µm frits                                                                               |                              |                             |
| Maximum pressure        | 120 bar                                                                                            |                              |                             |
| Operating temperature   | 18–25 °C                                                                                           |                              |                             |
| Column packing material | Affinity ligand-modified resin                                                                     |                              |                             |
| Immobilized ligands     | Affinity ligands to human albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin, and fibrinogen |                              |                             |
| Flow rate range         | 0.25–3.0 mL/min                                                                                    |                              |                             |
| Shipping solution       | Buffer A with 0.02% sodium azide                                                                   |                              |                             |
| Shipping temperature    | 2–8 °C (35–46 °F)                                                                                  |                              |                             |
| Storage temperature     | 2–8 °C (35–46 °F)                                                                                  |                              |                             |

\* Consult the column Certificate of Analysis to verify capacity for plasma samples.



## Related Agilent Products

Agilent Multiple Affinity Removal System spin cartridges and LC columns are listed in [Table 7](#) below.

**Table 7** Agilent Multiple Affinity Removal System spin cartridges and LC columns

| Product Group        | Proteins Removed                                                                                                                                                                               | Format                 | Capacity            | Part No.  |
|----------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|------------------------|---------------------|-----------|
| <b>Human-14</b>      | albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin, fibrinogen, alpha2-macroglobulin, alpha1-acid glycoprotein, IgM, apolipoprotein AI, apolipoprotein AII, complement C3, transthyretin | spin cartridge         | 8–10 µL plasma      | 5188-6560 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | up to 20 µL plasma  | 5188-6557 |
|                      |                                                                                                                                                                                                | 4.6 x 100 mm LC column | up to 40 µL plasma  | 5188-6558 |
|                      |                                                                                                                                                                                                | 10 x 100 mm LC column  | up to 250 µL plasma | 5188-6559 |
| <b>Human-7</b>       | albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin, fibrinogen                                                                                                                           | spin cartridge         | 12–14 µL plasma     | 5188-6408 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | 30–35 µL plasma     | 5188-6409 |
|                      |                                                                                                                                                                                                | 4.6 x 100 mm LC column | 60–70 µL plasma     | 5188-6410 |
|                      |                                                                                                                                                                                                | 10 x 100 mm LC column  | 250–300 µL plasma   | 5188-6411 |
| <b>Human-6HC</b>     | albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin                                                                                                                                       | spin cartridge         | 14–16 µL serum      | 5188-5341 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | 30–40 µL serum      | 5188-5332 |
|                      |                                                                                                                                                                                                | 4.6 x 100 mm LC column | 60–80 µL serum      | 5188-5333 |
|                      |                                                                                                                                                                                                | 10 x 100 mm LC column  | up to 340 µL serum  | 5188-5336 |
| <b>Human-6</b>       | albumin, IgG, IgA, transferrin, haptoglobin, antitrypsin                                                                                                                                       | spin cartridge         | 7–10 µL serum       | 5188-5230 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | 15–20 µL serum      | 5185-5984 |
|                      |                                                                                                                                                                                                | 4.6 x 100 mm LC column | 30–40 µL serum      | 5185-5985 |
| <b>Human-HSA/IgG</b> | albumin, IgG                                                                                                                                                                                   | spin cartridge         | up to 50 µL serum   | 5188-8825 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | up to 100 µL serum  | 5188-8826 |
| <b>Human-HSA</b>     | albumin                                                                                                                                                                                        | spin cartridge         | up to 75 µL serum   | 5188-5334 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | up to 175 µL serum  | 5188-6562 |
| <b>Mouse-3</b>       | albumin, IgG, transferrin                                                                                                                                                                      | spin cartridge         | 25–30 µL serum      | 5188-5289 |
|                      |                                                                                                                                                                                                | 4.6 x 50 mm LC column  | 37–50 µL serum      | 5188-5217 |
|                      |                                                                                                                                                                                                | 4.6 x 100 mm LC column | 75–100 µL serum     | 5188-5218 |

Additional related products for use with the Agilent Multiple Affinity Removal System are listed in [Table 8](#) below.

**Table 8** Additional related products

| Part Number | Description                                                                     | Content                                                                        |
|-------------|---------------------------------------------------------------------------------|--------------------------------------------------------------------------------|
| 5185-5986   | <b>Starter Reagent Kit for Multiple Affinity Removal System LC columns</b>      | Buffer A: 2 x 1 L                                                              |
|             |                                                                                 | Buffer B: 1 L                                                                  |
|             |                                                                                 | Spin filters 0.22 µm: 2 packs of 25                                            |
|             |                                                                                 | Protein concentrators: 1 pack of 25                                            |
| 5188-5254   | <b>Starter Reagent Kit for Multiple Affinity Removal System spin cartridges</b> | Buffer A: 1 L                                                                  |
|             |                                                                                 | Buffer B: 1 L                                                                  |
|             |                                                                                 | Spin filters 0.22 µm: 2 packs of 25                                            |
|             |                                                                                 | Protein concentrators: 1 pack of 25                                            |
|             |                                                                                 | Luer-Lok adapters: 1 pack of 2                                                 |
|             |                                                                                 | 5-mL plastic Luer-Lok syringes: 1 pack of                                      |
|             |                                                                                 | 1.5-mL microtubes: 6 packs of 100                                              |
|             |                                                                                 | Spin cartridge extra caps and plugs, 1 pack of 6 each                          |
| 5188-5231   | <b>mRP-C18 High Recovery Protein Fractionation and Desalting Column</b>         | 1 Column                                                                       |
|             |                                                                                 | (see <a href="http://www.agilent.com">www.agilent.com</a> for product details) |

## **In This Book**

This document describes how to use the Multi Affinity Removal Column, Human-7 to chromatographically remove interfering high-abundance proteins from human biological samples.

© Agilent Technologies, Inc. 2016, 2018

Version A1, December 2018



5990-7228



**Agilent Technologies**