

# ProteinChip® Human Serum Control

Catalog #C10-00006

## Contents

1 tube 500 µl pooled healthy human male serum.

## Storage

Store at -20°C. Avoid freeze-thawing.

## Biohazard Precautions

Appropriate personal protective equipment should always be worn when processing biohazardous samples. Dispose of biohazardous waste in accordance with local laws and/or environmental regulations.

## Example Profile

Figure 1 shows an example profile of the ProteinChip human serum control fractionated using the ProteinChip serum fractionation kit (catalog #K10-00007). The fractions were profiled using the ProteinChip WCX2 array (no longer available).

**Note:** Profiles may vary with instrument settings.

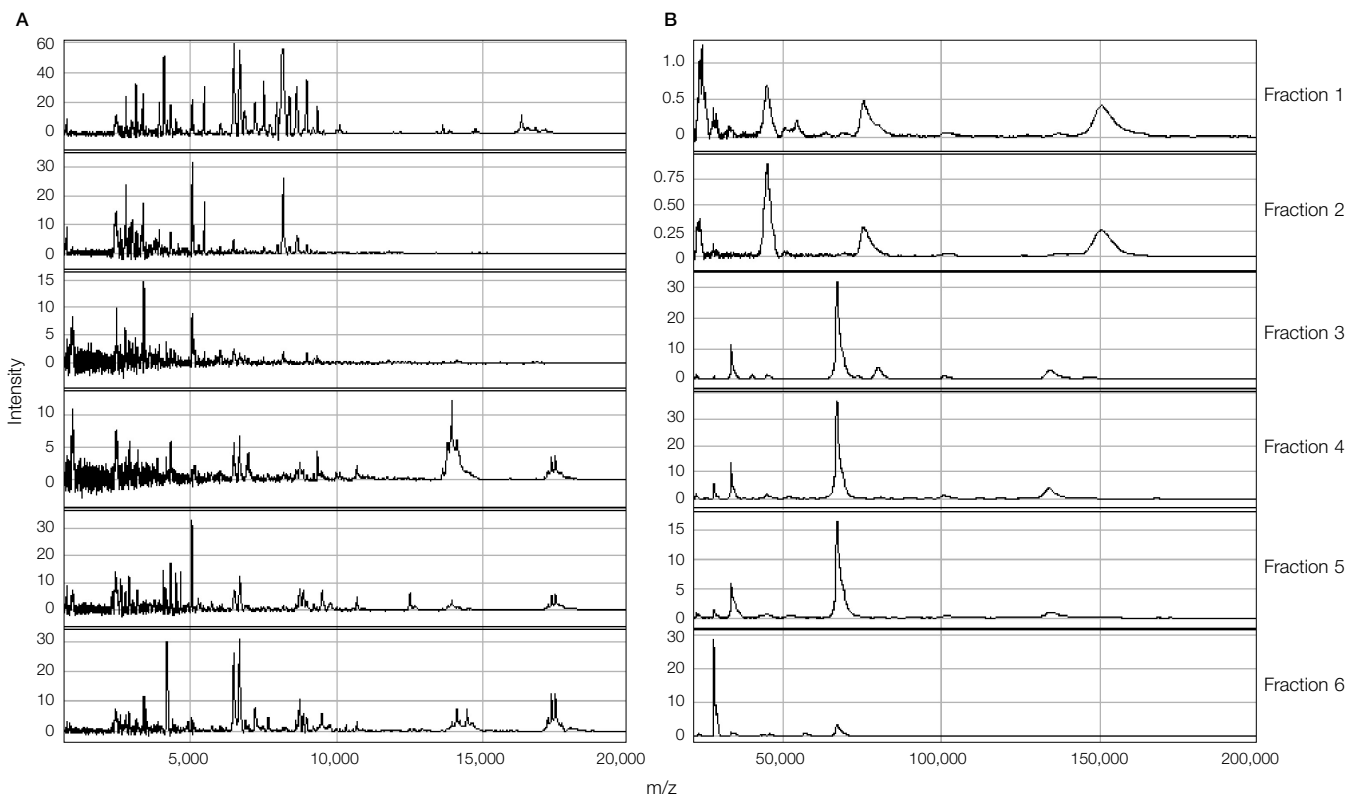


Fig. 1. Profile of fractionated ProteinChip human serum control. A, low molecular mass range, <20 kD; B, high molecular mass range, 20–200 kD.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400  
Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000