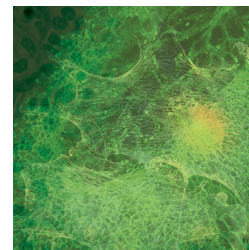


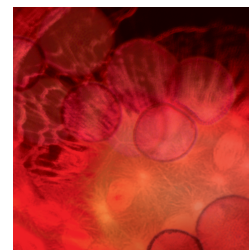
Sample Preparation

Sample preparation prior to 2-D electrophoresis is critical for producing meaningful, reproducible results. Bio-Rad has developed a number of sample preparation products for protein extraction, cleanup (contaminant removal), fractionation, and depletion. These products include chemical reagents, chromatography columns, and preparative electrophoresis tools.



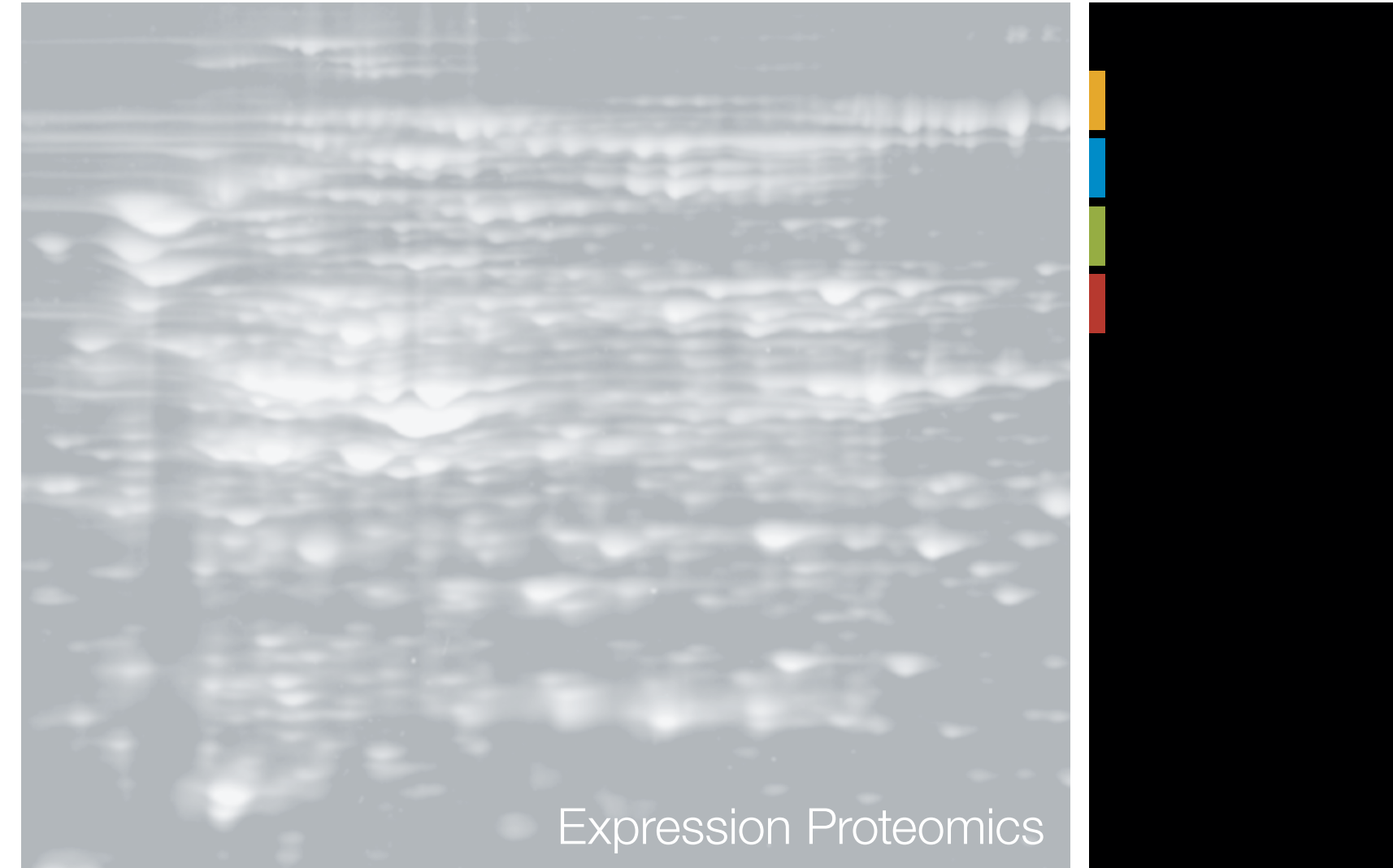
2-D Electrophoresis

One of the most powerful protein separation techniques available, 2-D electrophoresis has the ability to resolve complex mixtures of thousands of proteins simultaneously in a single gel. Depending on the experiment, requirements for resolution, throughput, and speed may vary. Bio-Rad addresses these varying requirements by offering choices for all components of the 2-D process.



Imaging and Analysis

Proteins separated by 2-D gel electrophoresis are stained for visualization, then imaged and analyzed to compare the protein spot patterns of the different samples. Bio-Rad provides a wide assortment of imaging systems and sophisticated PDQuest™ 2-D analysis software to simplify the entire analysis process.



Expression Proteomics

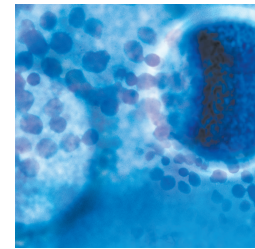
Tools for Protein Sample Preparation,
2-D Electrophoresis,
and Imaging and Analysis



Bio-Rad
Laboratories, Inc.

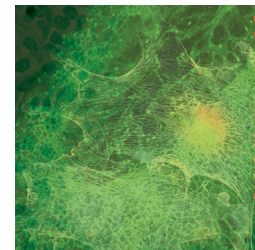
Life Science Group
Web site www.bio-rad.com
USA 800 4BIORAD
Canada 305 364 3435
China 86 21 6426 0808
Czech Republic 420 241 430 532
Denmark 44 52 10 00
Finland 09 804 22 00
France 01 47 95 69 65
Germany 089 316 84 0
Greece 30 210 777 4286
Hong Kong 852 2789 3300
Hungary 36 1 455 8800
India 91 124 4026300
Israel 03 963 0650
Italy 02 216051
Japan 03 6361 7000
Korea 82 2 3473 4460
Mexico 52 555 466 7670
The Netherlands 0316 540666
New Zealand 0608 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 3168
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





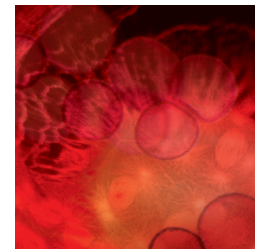
Sample Preparation

Sample preparation prior to 2-D electrophoresis is critical for producing meaningful, reproducible results. Bio-Rad has developed a number of sample preparation products for protein extraction, cleanup (contaminant removal), fractionation, and depletion. These products include chemical reagents, chromatography columns, and preparative electrophoresis tools.



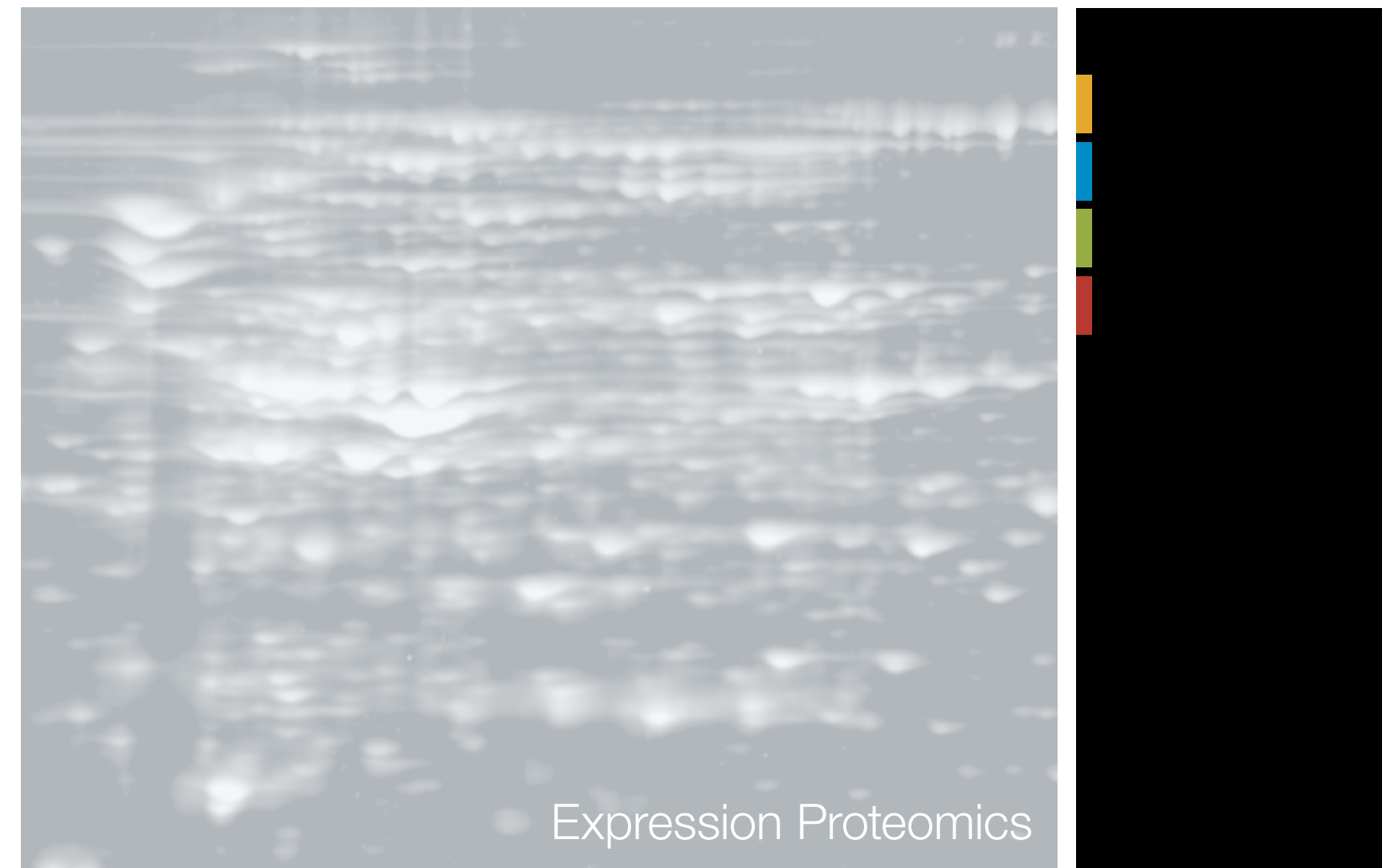
2-D Electrophoresis

One of the most powerful protein separation techniques available, 2-D electrophoresis has the ability to resolve complex mixtures of thousands of proteins simultaneously in a single gel. Depending on the experiment, requirements for resolution, throughput, and speed may vary. Bio-Rad addresses these varying requirements by offering choices for all components of the 2-D process.



Imaging and Analysis

Proteins separated by 2-D gel electrophoresis are stained for visualization, then imaged and analyzed to compare the protein spot patterns of the different samples. Bio-Rad provides a wide assortment of imaging systems and sophisticated PDQuest™ 2-D analysis software to simplify the entire analysis process.



Expression Proteomics

Tools for Protein Sample Preparation,
2-D Electrophoresis,
and Imaging and Analysis

www.expressionproteomics.com



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9814 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3337 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 03 Finland 09 804 22 00 France 01 47 95 89 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 01 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 498 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

