



# All you need for custom fills and formulas

Bio-Rad *Solutions*™ Program

We will meet your specifications for:  
Ready-to-use powder or liquid  
Packaging options  
Lot size  
In-process and final product QC

**BIO-RAD**

# bulk electrophoresis & chromatography reagents available:

Listed below are examples of ready-to-use chemical reagents and mixtures which are now available in either liquid or dry powder format (neat or concentrates), based on your specifications. Packaging options range from milliliters to liters and from grams to kilograms.



## electrophoresis buffers and reagents

Glycine  
TBE  
Tris  
SDS  
Urea

## stains

Coomassie® stains  
Silver  
Zinc

## chromatography buffers

Hepes  
MAPS® buffers  
MES  
Tris

## solvents

Acetonitrile  
Ethanol  
Methanol

## ELISA reagents

Antibody conjugates  
Enzyme conjugates  
Blockers

*Coomassie is a trademark of ICI Organics.*

Bio-Rad has been your trusted source for research-scale electrophoresis and chromatography reagents and supplies for over 40 years. Our Quality Management Systems are ISO 9001 approved, and many of our products have Drug Master Files (DMFs). As a special service, Bio-Rad can prepare and record DMFs for your made-to-order products. For more information, contact your local Bio-Rad sales representative.

**BIO-RAD**

Bio-Rad  
Laboratories

Life Science  
Group

Web Site [www.bio-rad.com](http://www.bio-rad.com) U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771  
China 86-10-62051850/51 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France 01 43 90 46 90 Germany 089 318 84-0  
Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 39-02-216091 Japan 03-5811-6270 Korea 82-2-3473-4460  
Latin America 305-894-5950 Mexico 514-2210 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 22-74-18-70 Russia 7 095 979 98 00  
Singapore 65-2729877 Spain 34-91-661-7085 Sweden 46 (0)8-55 51 27 00 Switzerland 01-809 55 55 United Kingdom 0800-181134