



Bio-Plex™

x-Plex™ Assays

Personal Selection

BIO-RAD

x-Plex Mixed-to-Order Assays

Bio-Plex x-Plex assays are designed for researchers interested in performing multiplex assays with a specific combination of Bio-Plex assay targets that is not available in our preconfigured assay panels. x-Plex assays are made to order with any combination of available Bio-Plex targets. You select the targets specific for your study, and our x-Plex service provides you with a premixed, ready-to-use assay kit. Visit www.bio-rad.com/bioplex/x-plex/ to select your preferred combination of human, mouse, or rat cytokines or phosphoprotein targets. With x-Plex you add the benefit of custom assay configuration to the existing advantages of Bio-Plex assays:

- Premixed conjugated beads
- Premixed detection antibodies
- Premixed standards (cytokines) or controls (phosphoproteins)
- 1-pack and 10-pack options



Take Your Pick — Mixed-to-Order Assays for Any Combination

It's All About You and What You Need

If the particular combination of cytokine or phosphoprotein targets you need for your research is not available in a Bio-Plex preconfigured assay panel, order what you need with premixed x-Plex assays. You select only the analytes you want to test, and we custom make the assay panel for you. The x-Plex assay panel includes premixed conjugated beads, premixed detection antibodies, and premixed standards (for cytokines) or controls (for phosphoprotein assays). Order buffer, cell lysis, and diluent kits to complete your package. This is Bio-Plex x-Plex at your service.

Examine Your x-Plex Assay Options Online

Start by viewing the available selection of x-Plex assay targets online at www.bio-rad.com/bioplex/x-plex/. Visit this site often to get the most up-to-date offering of Bio-Plex assay targets. Use the pull-down menu to choose the species, and the window will display the available assay targets. Simply click the boxes to make your selections, then click “*calculate panel*” to view your custom catalog number as well as the companion buffer and diluent kits. By selecting “*request quote*”, you will be sent a real-time confirmation e-mail with your catalog number; your Bio-Plex representative will contact you with pricing and a ship date (~1 week worldwide). Simply place your order for your x-Plex assay through your normal purchasing process.

The screenshot shows the Bio-Rad website's x-Plex assay configuration page. The page title is "Welcome to Bio-Rad Laboratories". The main heading is "x-Plex Assays". Below this, there are several sections:

- Panel and Package Size:** A dropdown menu for "Human Cytokines" and a package size selector set to "1 x 96".
- Select All/Calculate Panel/View Selection/Request Quote:** A series of buttons for managing the assay panel.
- Assay Targets:** A grid of checkboxes for selecting specific cytokines or phosphoproteins. The visible options include:
 - IL-1beta
 - IL-2
 - IL-4
 - IL-6
 - IL-7
 - IL-10
 - IL-12 (p70)
 - IL-13
 - IL-17
 - G-CSF
- Instructions:** Text explaining that users should select the assay type and click the x-Plex catalog number and request a quote to get more information from a representative.
- Required Items:** A table listing necessary reagents:

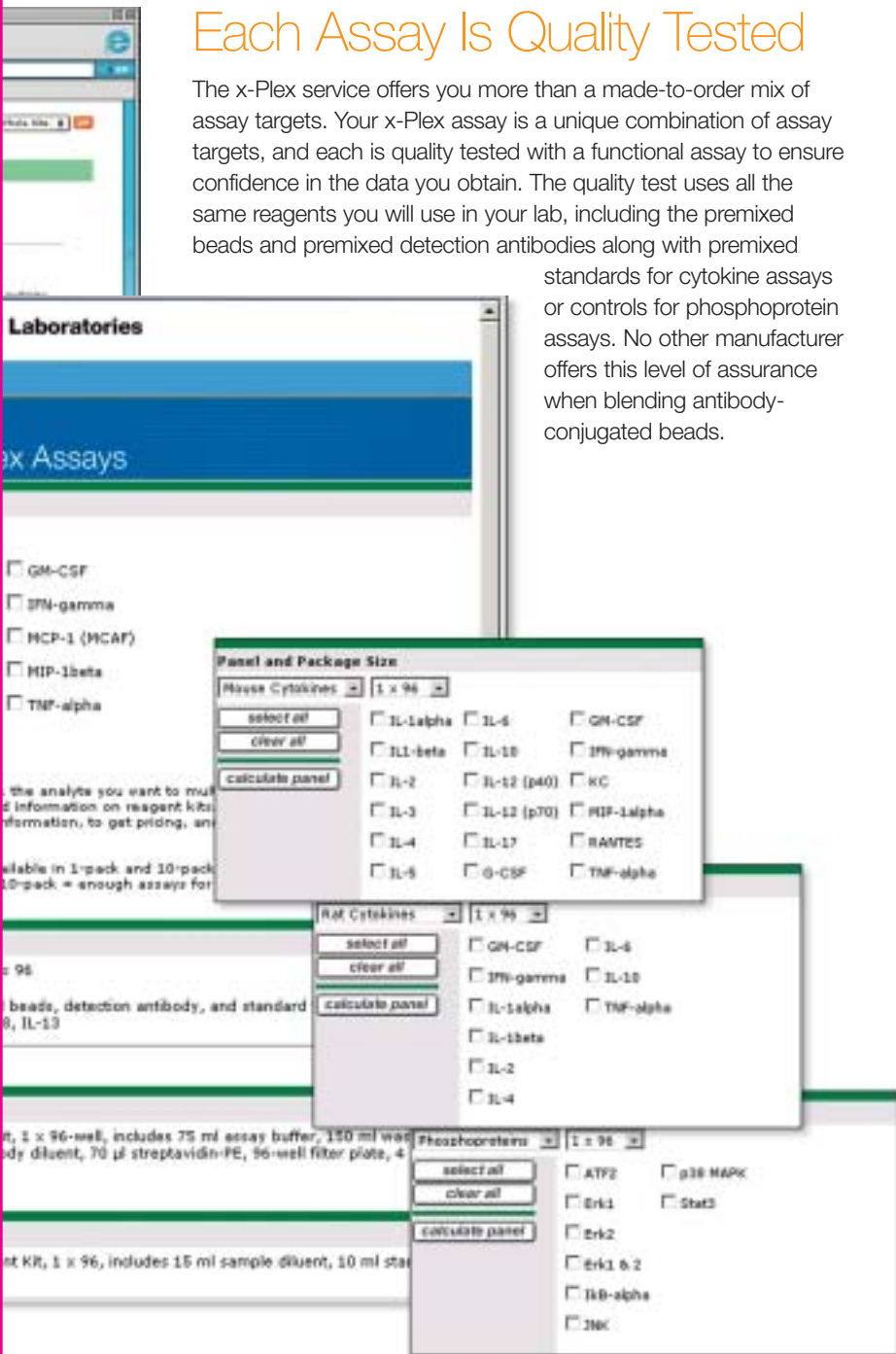
Catalog #	Description
171-304000	Bio-Plex Cytokine Reagent Kit buffer, 15 ml detection antibody sealing tape, instructions
- Optional Items:** A table listing optional reagents:

Catalog #	Description
171-305000	Bio-Plex Human Serum Diluent

Order Assays Combination

Each Assay Is Quality Tested

The x-Plex service offers you more than a made-to-order mix of assay targets. Your x-Plex assay is a unique combination of assay targets, and each is quality tested with a functional assay to ensure confidence in the data you obtain. The quality test uses all the same reagents you will use in your lab, including the premixed beads and premixed detection antibodies along with premixed standards for cytokine assays or controls for phosphoprotein assays. No other manufacturer offers this level of assurance when blending antibody-conjugated beads.

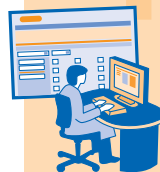


How our delivery system works

1

Make your selection

Configure your x-Plex panel online at www.bio-rad.com/bioplex/x-plex/. Then place your order using your unique x-Plex catalog number.



2

Made to order

Our highly trained specialists will custom mix your assay to your specifications.



3

Quality tested

All custom mixes will be thoroughly tested by our quality control department, then carefully packaged and shipped to you.



4

Get just what you need

Receive the custom assay panel made specifically for your research.



Ordering Information

x-Plex Assays

Once you have obtained a catalog number specific to your x-Plex assay through www.bio-rad.com/bioplex/x-plex/, place your order as with any other Bio-Rad product. For cytokine x-Plex assays, you also need the universal cytokine reagent kit and a serum diluent kit if you are testing serum samples. For phosphoprotein x-Plex assays, you also need the universal phosphoprotein reagent kit and a cell lysis kit for lysate preparation. When reordering your x-Plex panel, simply use the same x-Plex catalog number.

Catalog #	Description
x-Plex Assays (numbers generated online)	
x-Number	Bio-Plex Human Cytokine x-Plex Panel
x-Number	Bio-Plex Mouse Cytokine x-Plex Panel
x-Number	Bio-Plex Rat Cytokine x-Plex Panel
x-Number	Bio-Plex Phosphoproteins x-Plex Panel
Reagent Kits	
171-304000	Bio-Plex Cytokine Reagent Kit , 1 x 96-well
171-304001	Bio-Plex Cytokine Reagent Kit , 10 x 96-well
171-304004	Bio-Plex Phosphoprotein Reagent Kit , 1 x 96-well
171-304005	Bio-Plex Phosphoprotein Reagent Kit , 10 x 96-well
Sample Preparation Kits	
171-304011	Bio-Plex Cell Lysis Kit , 1 x 96-well
171-304012	Bio-Plex Cell Lysis Kit , 10 x 96-well
171-305000	Bio-Plex Human Serum Diluent Kit , 1 x 96-well
171-305001	Bio-Plex Human Serum Diluent Kit , 10 x 96-well
171-305004	Bio-Plex Mouse Serum Diluent Kit , 1 x 96-well
171-305005	Bio-Plex Mouse Serum Diluent Kit , 10 x 96-well
171-305008	Bio-Plex Rat Serum Diluent Kit , 1 x 96-well

Cell Signaling Analysis Just Got Easier

Maximize the benefits of the Bio-Plex system and profile both the extracellular cytokines and intracellular kinases present in the same sample. Use the cell culture supernatant with Bio-Plex cytokine assays, and the cell pellet for activated kinases with phosphoprotein assays.

What You Need For ...

Cytokine Testing	Phosphoprotein Testing
Human, mouse, or rat samples	Mammalian samples
 <p>Bio-Plex Cytokine Assays</p>	 <p>Bio-Plex Phosphoprotein Assays</p>
 <p>Bio-Plex Cytokine Reagent Kit</p>	 <p>Bio-Plex Cell Lysis Kit</p>
 <p>Bio-Plex Diluent Kit (optional)</p>	 <p>Bio-Plex Phosphoprotein Reagent Kit</p>



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191 Canada (905) 712-2771 China (86-21) 63052255 Czech Republic + 420 2 41 43 05 32 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84-1777 Hong Kong 852-2789-3300 India (91-124)-6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20 The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 331 99 99 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-6415 3188 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000