

Bio-Plex Custom Assay Development Reagents

Unlimited Design Capabilities

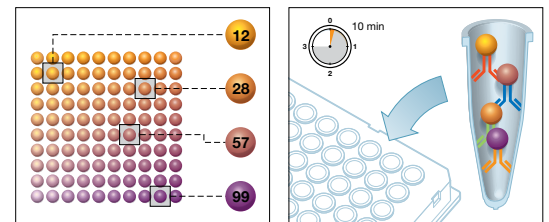
- Choose antibodies, antigens, receptors, ligands, or enzymes as targets
- Covalently couple targets to beads in less than 5 hours
- Multiplex up to 100 targets on the Bio-Plex 200 and up to 500 targets on the Bio-Plex 3D Systems
- Also compatible with the Luminex 200 and FlexMAP 3D Systems

Fulfill your unique research needs by developing custom immunoassays. Bio-Plex COOH Beads and Coupling Kits allow you to create unique multiplex assays specific to your research needs. We recommend Bio-Plex COOH Beads for researchers who:

- Are testing proprietary or unique targets or samples
- Require assay customization
- Require a broader assay selection than is commercially available
- Require flexibility to change targets often
- Have assay development resources
- Have access to a Luminex or Bio-Plex System

Unleash the power of multiplexing by combining beads to gain more information from a single sample of 12.5 μ l or less. Utilizing beads as a substrate enables simultaneous detection and measurement of up to 500 different targets in a single well, saving your precious sample,

hands-on time, and money. The Bio-Plex Reader identifies beads by a signature dye ratio and corresponding bead number. The system uses a reporter dye added just prior to reading to determine the amount of analyte that has bound to the bead.

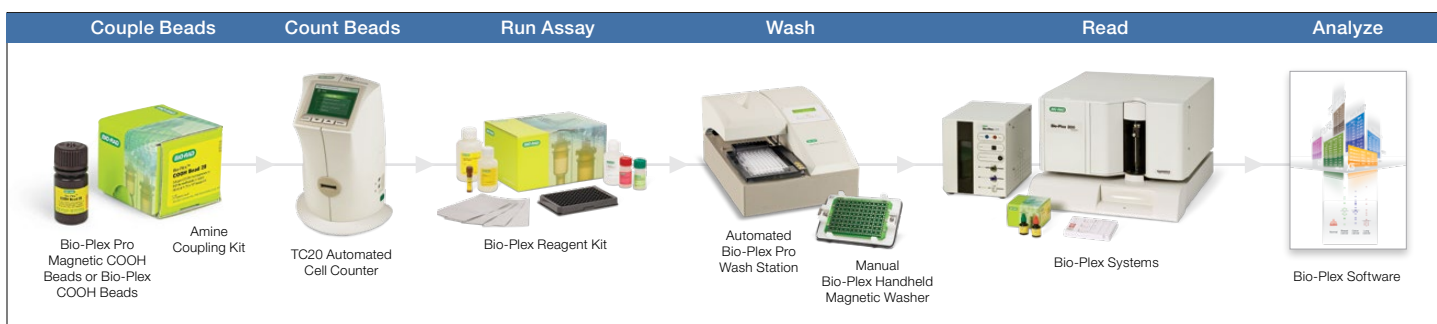


Couple individual proteins (antibodies, antigens, enzymes, ligands, or receptors) to specific bead regions and combine them to create a multiplex assay. Developing custom assays gives you flexibility not only in the targets you choose but also the type of assay, including sandwich, competitive, direct, and indirect assays.

Bead coupling is the first step of developing bead-based multiplex assays. Bio-Rad offers a selection of spectrally distinct COOH beads for use in assay development, including nonmagnetic beads and magnetic beads. The coupling procedure involves a rapid two-step carbodiimide reaction between primary amine groups on the protein and carboxyl functional groups on the bead surfaces. Magnetic beads streamline the coupling reaction and increase assay precision with magnetic separation washes.

Get your data faster while saving money, as you don't have to wait for weeks and pay for commercial custom coupling results. Being able to perform and analyze a multiplex assay reduces need for consumables, buffers, and analysis time versus multiple individual ELISAs.

Custom singleplex and multiplex assays leverage high-performance instruments, industry-leading software, and integrated calibration and validation kits to obtain high-quality data from limited sample volumes.



Bio-Rad has tools to aid you in each step of the process.

1. Couple beads to a protein of your choice with the Bio-Plex Amine Coupling Kit.
2. Get an accurate bead count with the TC20 Automated Cell Counter in less than 30 sec.
3. Run the immunoassays using Bio-Plex Reagent Kits, powered by Luminex xMAP technology.
4. Eliminate manual washes and improve results with the Bio-Plex Pro Wash Station, designed specifically to perform optimal washes for Bio-Plex Assays. Or use the Bio-Plex Handheld Magnetic Washer for manual washes.
5. Measure the multiplexed analytes using the Bio-Plex 200 (up to 100 biomolecules) or Bio-Plex 3D Suspension Array (up to 500 biomolecules) Systems.
6. Analyze with Bio-Plex Data Pro Software, which provides tools to organize and visualize multiple datasets generated from xMAP multiplex experiments.

Ordering Information

Catalog #	Description
171406001	Bio-Plex Amine Coupling Kit
171304501	Bio-Plex Streptavidin-PE , 10 x 96-well reactions
1614916	16-Tube SureBeads Magnetic Rack
1450102	TC20 Automated Cell Counter , includes instrument, power supply, USB flash drive, 30 dual-chamber counting slides (60 counts), 1.5 ml trypan blue
30034376	Bio-Plex Pro Wash Station
171020100	Bio-Plex Handheld Magnetic Washer
171000201	Bio-Plex 200 System , 100–240 V, multiplex array system, includes reader, microplate platform, Bio-Plex Manager 6 Software, PC, calibration kit, validation kit 4.0, MCV plate IV, reservoir, sheath fluid, 2 sample needles
171000205	Bio-Plex 200 System with HTF , same as 171000201 with high-throughput fluidics (HTF)
BIOPLEX3D	Bio-Plex 3D Suspension Array System , includes instrument, xPONENT Software for data acquisition, PC, calibration and verification reagents, Bio-Plex Manager 6 Desktop License

Visit bio-rad.com/BpxBeadCoupling for more information including catalog numbers for COOH beads.

BIO-RAD and BIO-PLEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation. Luminex and xMAP are trademarks of Luminex Corporation. All trademarks used herein are the property of their respective owner.



Bio-Rad Laboratories, Inc.

Life Science Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 6511 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

