

# Ordering Information

## Bio-Plex™ COOH Beads

### How to Order

Each milliliter of Bio-Plex COOH beads contains enough beads for 10 coupling reactions. This volume is sufficient to produce enough protein-coupled beads for up to twenty 96-well microplates using 5,000 beads per well.



### COOH Beads

Bead		Catalog #		Bead		Catalog #		Bead		Catalog #		Bead		Catalog #		Bead		Catalog #	
1 ml		16 ml		1 ml		16 ml		1 ml		16 ml		1 ml		16 ml		1 ml		16 ml	
1	171-506001	171-606001	21	171-506021	171-606021	41	171-506041	171-606041	61	171-506061	171-606061	81	171-506081	171-606081					
2	171-506002	171-606002	22	171-506022	171-606022	42	171-506042	171-606042	62	171-506062	171-606062	82	171-506082	171-606082					
3	171-506003	171-606003	23	171-506023	171-606023	43	171-506043	171-606043	63	171-506063	171-606063	83	171-506083	171-606083					
4	171-506004	171-606004	24	171-506024	171-606024	44	171-506044	171-606044	64	171-506064	171-606064	84	171-506084	171-606084					
5	171-506005	171-606005	25	171-506025	171-606025	45	171-506045	171-606045	65	171-506065	171-606065	85	171-506085	171-606085					
6	171-506006	171-606006	26	171-506026	171-606026	46	171-506046	171-606046	66	171-506066	171-606066	86	171-506086	171-606086					
7	171-506007	171-606007	27	171-506027	171-606027	47	171-506047	171-606047	67	171-506067	171-606067	87	171-506087	171-606087					
8	171-506008	171-606008	28	171-506028	171-606028	48	171-506048	171-606048	68	171-506068	171-606068	88	171-506088	171-606088					
9	171-506009	171-606009	29	171-506029	171-606029	49	171-506049	171-606049	69	171-506069	171-606069	89	171-506089	171-606089					
10	171-506010	171-606010	30	171-506030	171-606030	50	171-506050	171-606050	70	171-506070	171-606070	90	171-506090	171-606090					
11	171-506011	171-606011	31	171-506031	171-606031	51	171-506051	171-606051	71	171-506071	171-606071	91	171-506091	171-606091					
12	171-506012	171-606012	32	171-506032	171-606032	52	171-506052	171-606052	72	171-506072	171-606072	92	171-506092	171-606092					
13	171-506013	171-606013	33	171-506033	171-606033	53	171-506053	171-606053	73	171-506073	171-606073	93	171-506093	171-606093					
14	171-506014	171-606014	34	171-506034	171-606034	54	171-506054	171-606054	74	171-506074	171-606074	94	171-506094	171-606094					
15	171-506015	171-606015	35	171-506035	171-606035	55	171-506055	171-606055	75	171-506075	171-606075	95	171-506095	171-606095					
16	171-506016	171-606016	36	171-506036	171-606036	56	171-506056	171-606056	76	171-506076	171-606076	96	171-506096	171-606096					
17	171-506017	171-606017	37	171-506037	171-606037	57	171-506057	171-606057	77	171-506077	171-606077	97	171-506097	171-606097					
18	171-506018	171-606018	38	171-506038	171-606038	58	171-506058	171-606058	78	171-506078	171-606078	98	171-506098	171-606098					
19	171-506019	171-606019	39	171-506039	171-606039	59	171-506059	171-606059	79	171-506079	171-606079	99	171-506099	171-606099					
20	171-506020	171-606020	40	171-506040	171-606040	60	171-506060	171-606060	80	171-506080	171-606080	100	171-506000	171-606000					

Beads highlighted in tan are currently available off the shelf and will ship the next day. Orders for any of the remaining beads will ship within 4–6 weeks (minimum volume requirements apply).

Catalog # Description

### Beads

(See table) **Bio-Plex COOH Beads**,  $1.25 \times 10^7$  beads/ml, 1 ml

(See table) **Bio-Plex COOH Beads**,  $1.25 \times 10^7$  beads/ml, 16 ml

### Bio-Plex System and System Performance Validation Tools

- 171-000001 **Bio-Plex Suspension Array System**, 100–240 V, includes array reader, microplate platform, Bio-Plex Manager™ software, PC, monitor, calibration kit, validation kit, maintenance, calibration, and validation (MCV) plate III, 20 L sheath fluid, spare sample needle
- 171-000005 **Bio-Plex Suspension Array System**, 100–240 V, same as 171-000001, with high-throughput fluidics (HTF) system
- 171-000050 **High-Throughput Fluidics (HTF) System**, 100–240 V
- 171-000055 **Sheath Fluid**, 20 L
- 171-001008 **Bio-Plex Manager Version 4.0 Software Upgrade Kit**, for version 3.0 users
- 171-001012 **Bio-Plex Manager Software License**, version 4.0 Security Edition workstation, single user
- 171-203001 **Bio-Plex Validation Kit 4.0**, includes optics validation, reporter validation, classify validation, and fluidics validation bead sets for approximately 50 validation routines using Bio-Plex Manager 4.0 and MCV plate III
- 171-203032 **Bio-Plex MCV Plate III**, for use with Bio-Plex Manager software version 4.0
- 171-203060 **Bio-Plex Calibration Kit**, includes CAL1 and CAL2 calibration beads for approximately 50 daily calibration routines

### Reagent Kits

- 171-304000 **Bio-Plex Cytokine Reagent Kit**, 1 x 96-well, includes assay buffer, wash buffer, detection antibody diluent, streptavidin-PE (100x), 96-well sterile filter plate, sealing tape, instructions
- 171-304004 **Bio-Plex Phosphoprotein Detection Reagent Kit**, 1 x 96-well, includes assay buffer, wash buffer, antibody diluent, resuspension buffer, streptavidin-PE (100x), 96-well sterile filter plate, sealing tape, instructions

### Sample Preparation Kits

- 171-305000 **Bio-Plex Human Serum Diluent Kit**, 1 x 96-well, includes sample diluent, standard diluent
- 171-305004 **Bio-Plex Mouse Serum Diluent Kit**, 1 x 96-well, includes sample diluent, standard diluent
- 171-305008 **Bio-Plex Rat Serum Diluent Kit**, 1 x 96-well, includes sample diluent, standard diluent
- 171-304011 **Bio-Plex Cell Lysis Kit**, 1 x 96-well, includes cell wash buffer, cell lysis buffer, factors, instructions

**BIO-RAD**