

# Bio-Plex Pro™ Human Apolipoprotein 10-Plex Assay

## Quick Guide

For Use with	Instruction Manual #
Bio-Plex Pro Human Apolipoprotein 10-Plex Assay	10000077901

This quick guide can be used to prepare and run a full 1 x 96-well assay plate. For more information on a given step, refer to the corresponding section of the Bio-Plex Pro Human Apolipoprotein 10-Plex Assay Instruction Manual (#10000077901). Go to [bio-rad.com/web/bio-plex](http://bio-rad.com/web/bio-plex) to download the instruction manual, which includes detailed instructions and a list of kit components.

**IMPORTANT!** Pay close attention to vortexing, shaking, and incubation instructions. Deviation from the protocol may result in low assay signal and assay variability.

## Sample Preparation

1. Centrifuge serum or plasma samples at **1,000 x g** for **15 min** at 4°C to remove particulates from all samples prior to use.
2. Prepare sample dilutions in **0.5** or **1.0 ml** polypropylene tubes as required for the assay.
3. Prepare the sample dilution as outlined in Table 1. The dilution example in Table 1 provides a sufficient amount to run each sample in duplicate.

**Table 1. Preparation of the sample dilutions.**

Sample Dilution	Volume of Sample, $\mu$ l	Volume of Sample Dilution Buffer, $\mu$ l
1:50,000	(a) Prepare 1:10 10	90
	(b) Prepare 1:50 10 (a)	490
	Prepare 1:100 5 (b)	495

**Note:** Controls are ready to use after reconstitution. No dilution is needed.

## Bio-Plex Pro Human Apolipoprotein 10-Plex Assay

### Reagent Preparation

1. Reconstitute the following lyophilized reagents in distilled water (dH<sub>2</sub>O) before use, as outlined in Table 2.

**Table 2. Reconstitution of the lyophilized reagents.**

Reagent	Volume dH <sub>2</sub> O
Standards mix	150 µl
Control 1	100 µl
Control 2	100 µl
Blocking buffer	1.5 ml
Standard diluent	1.0 ml
Detection antibodies	4.8 ml

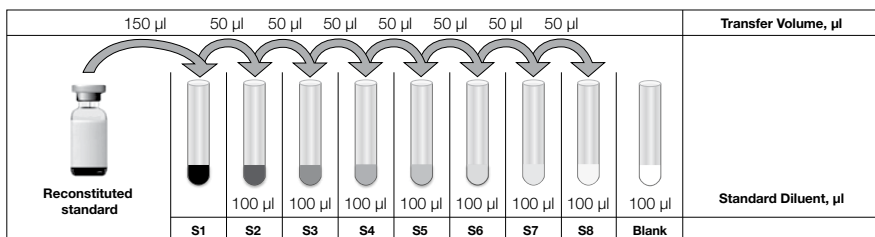
- a. Allow vial to sit at room temperature for **5 min**.
  - b. Mix by **vortexing** at a medium setting.
2. Bring the 10x assay buffer to ambient temperature (RT).
    - a. Mix by inversion to ensure all salts are in solution.
    - b. Prepare 1x assay buffer by diluting **1 part** 10x assay buffer (60 ml) with **9 parts** of dH<sub>2</sub>O (540 ml).

### Dilution of Standard (1:3 serial dilution)

1. Label 9 polypropylene tubes S1 through S8 and Blank.
2. Transfer the reconstituted standard into the tube labeled **S1**.
3. Add the appropriate amount of standard diluent into the labeled tubes as outlined in Table 3 (this will be sufficient for duplicate standard curves and blanks).

**Table 3. Dilution of the standards.**

Standard	Volume of Standard Diluent, µl	Volume of Standard, µl
S1	—	150 from reconstituted vial
S2	100	50 of <b>S1</b>
S3	100	50 of <b>S2</b>
S4	100	50 of <b>S3</b>
S5	100	50 of <b>S4</b>
S6	100	50 of <b>S5</b>
S7	100	50 of <b>S6</b>
S8	100	50 of <b>S7</b>
Blank	100	—



4. Prepare working standards (**S2–S8**) by serial dilution. Transfer the appropriate volume of standard into each of the labeled tubes with standard diluent as outlined above.
5. **Vortex** each standard at a medium setting before proceeding with the next serial dilution. Change the pipet tip at each dilution step.
6. The **Blank** tube consists of standard diluent alone.

**Dispensing Reagents**

---

1. Add **10 µl** of blocker to all wells of the plate.
2. Add **30 µl** of the standard, control, sample, or blank to the appropriate well of the plate.
3. **Vortex** the capture beads at medium speed for **10–20 sec**. Add **10 µl** of the beads to each well of the plate.
4. Cover the plate with a plate seal and protect it from light with aluminum foil. Incubate on a shaker at **850 ± 50 rpm** for **1 hr** at RT.
5. Wash the plate three times with **100 µl** per well of 1x assay buffer.
6. **Vortex** the reconstituted detection antibodies at medium speed for **10–20 sec**. Add **40 µl** to each well.
7. Cover the plate as in step 4 and incubate on a shaker at **850 ± 50 rpm** for **1 hr** at RT. **Do not aspirate after incubation.**
8. Prepare the required dilution of streptavidin-phycoerythrin (SA-PE) as outlined in Table 4.

**Note:** Volumes in Table 4 are for an entire 96-well plate. Different volumes can be prepared, provided that the dilution ratios are maintained.

**Table 4. Dilution of SA-PE.**

SA-PE Dilution	Volume of SA-PE, µl	Volume of 1x Assay	
		Buffer, µl	Total Volume, µl
1:10	225	2,025	2,250

9. Add **20 µl** of diluted SA-PE to each plate well.
10. Cover the plate as in step 4 and incubate on a shaker at **850 ± 50 rpm** for **30 min** at RT.
11. Wash the plate three times with **100 µl** per well of 1x assay buffer.
12. After the final wash, resuspend the beads in each assay well with **100 µl** 1x assay buffer. Cover the plate as in step 4 and shake at **850 ± 50 rpm** for **30 sec**.
13. Remove the plate seal and read the plate at low photomultiplier tube (PMT) (Bio-Plex® 200 System), standard PMT (Bio-Plex 3D Suspension Array System), or default setting (Bio-Plex® MAGPIX™ Multiplex Reader).

MAGPIX is a trademark of Luminex Corporation.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800  
**Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532  
**Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00 **France** 33 01 47 95 69 65  
**Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100  
**India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000  
**Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666  
**New Zealand** 64 9 415 2280 **Norway** 47 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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700  
**Switzerland** 41 026 674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311  
**United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

