

Bio-Plex Pro Assay Quick Guide

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

- Cover and incubate at 850 ± 50 rpm, as in Step 4, for 30 min at RT.
- Wash the plate three times with 100 μ l 1x assay buffer.
- After the final wash, resuspend the beads in 100 μ l assay buffer. Cover plate as in Step 4 and shake the plate at 850 ± 50 rpm for 30 sec.
- Remove the plate seal and read plate at low PMT (Bio-Plex[®] 200), standard PMT (Bio-Plex 3D), or default settings (Bio-Plex[®] MAGPIX[™]).



BIO-RAD

**Bio-Rad
Laboratories, Inc.**



Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800
Austria 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500
Czech Republic 420 241 430 532 **Denmark** 44 52 10 00
Finland 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0
Greece 30 210 9532 220 **Hong Kong** 852 2789 3300
Hungary 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460
Mexico 52 555 488 7670 **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 7 495 721 14 04 **Singapore** 65 6415 3188
South Africa 27 861 246 723 **Spain** 34 91 590 5200
Sweden 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000

BIO-RAD

Bio-Plex Pro[™] RBM Kidney Toxicity Assays Quick Guide

For Use With	Instruction Manual #
Bio-Plex Pro [™] RBM Human, Rat, and Canine Kidney Toxicity Assays	10028258

This 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 complete instruction manual. New users can download the manual, which includes detailed instructions and a list of kit components, at www.bio-rad.com/bio-plex.

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

A. Reagent Preparation

- Reconstitute the following lyophilized reagents in dH₂O before use according to the table below.

Reagent	Volume, μ l	Reagent	Volume, ml
Standards mix	150	Blocking buffer	1.5
Control 1	100	Standard diluent	1.0
Control 2	100	Detection antibodies	4.8

- Allow vial to sit at room temperature for a minimum of 5 min, not to exceed 30 min.
- Mix by vortexing at a medium setting.
- Bring the 10x assay buffer to room temperature (RT).
 - Mix by inversion to ensure all salts are into solution.
 - Prepare 1x assay buffer — dilute 1 part 10x assay buffer with 9 parts of dH₂O.

B. Dilution of Standard (1:3 Serial Dilution)

1. Label 8 polypropylene tubes **S1** through **S8**.
2. Transfer the reconstituted standard into the tube labeled "**S1**."
3. Add the appropriate amount of the standard diluent into the labeled tubes according to the table below (this will be sufficient for duplicate standard curves and blanks).

Standard	Volume of Standard Diluent, μ l	Volume of Standard, μ l
S2	100	50 of S1
S3	100	50 of S2
S4	100	50 of S3
S5	100	50 of S4
S6	100	50 of S5
S7	100	50 of S6
S8	100	50 of S7
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 pipet tip at each dilution step.

C. Sample Preparation

1. Centrifuge samples at **500 x g** for **5 min** to remove particulates from all samples prior to use.
2. Prepare sample dilutions in **0.5 ml** or **1.0 ml** polypropylene tubes as required for the assay.
3. Dilution scenarios provided below are sufficient to run each sample in duplicate.

Panel	Sample Dilution	Volume of Urine Sample, μ l	Volume of Sample Buffer, μ l
Human Tox 1	1:4	20	60
Human Tox 2	1:50	10	490
Rat Tox 1	1:2	40	40
Rat Tox 2	1:50	10	490
Rat Albumin	1:10,000	5 (A. Prepare 1:100) 5 (B. Prepare 1:100)	495 495
Canine Tox 1	1:15	10	140
Canine Albumin	1:10,000	5 (C. Prepare 1:100) 5 (D. Prepare 1:100)	495 495

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

D. Dispensing of Reagents

1. Add **10 μ l** of blocker to all wells of the plate.
 2. Add **30 μ l** 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 μ l** of the beads to all wells of the plate.
 4. Cover plate with plate seal and protect from light with aluminum foil. Incubate on shaker at **850 \pm 50 rpm** for **1 hr** at RT.
 5. Wash the plate three times with **100 μ l** 1x assay buffer.
 6. **Vortex** the reconstituted detection antibodies at medium speed for **10–20 sec**. Add **40 μ l** to each well.
 7. Cover and incubate at **850 \pm 50 rpm**, as in Step 4, for **1 hr** at RT. **Do not aspirate after incubation.**
 8. Prepare the required dilution of SA-PE as outlined in the following table.
- Note:** Volumes in the table are for an entire 96-well plate. Smaller volumes can be prepared, provided that dilution ratios are maintained.
9. Add **20 μ l** of diluted SA-PE to the required plate wells.