

## SAMPLE PROTOCOL FOR NO WASH CAPTURE SANDWICH IMMUNOASSAY

*Microspheres should be protected from prolonged exposure to light throughout this procedure.*

1. Select the appropriate antibody-coupled microsphere sets.
2. Resuspend the microspheres by vortex and sonication for approximately 20 seconds.
3. Prepare a Working Microsphere Mixture by diluting the coupled microsphere stocks to a final concentration of 200 microspheres of each set/ $\mu\text{L}$  in PBS-1% BSA. (Note: 25  $\mu\text{L}$  of Working Microsphere Mixture is required for each reaction.) See **Technical Note 1**.
4. Aliquot 25  $\mu\text{L}$  of the Working Microsphere Mixture to the appropriate wells of a round-bottom microtiter plate.
5. Add 25  $\mu\text{L}$  of PBS-1%BSA to each background well.
6. Add 25  $\mu\text{L}$  of standard or sample to the appropriate wells.
7. Mix the reactions gently by pipetting up and down several times with a multi-channel pipettor.
8. Cover the plate and incubate for 30 minutes at room temperature.
9. Dilute the biotinylated detection antibody to the appropriate concentration in PBS-1% BSA. (Note: 25  $\mu\text{L}$  of diluted detection antibody is required for each reaction). See **Technical Note 2**.
10. Add 25  $\mu\text{L}$  of the diluted detection antibody to each well.
11. Mix the reactions gently by pipetting up and down several times with a multi-channel pipettor.
12. Cover the plate and incubate for 30 minutes at room temperature.
13. Dilute streptavidin-R-phycoerythrin to the appropriate concentration in PBS-1% BSA. (Note: 25  $\mu\text{L}$  of diluted streptavidin-R-phycoerythrin is required for each reaction). See **Technical Note 2**.
14. Add 25  $\mu\text{L}$  of the diluted streptavidin-R-phycoerythrin to each well.
15. Mix the reactions gently by pipetting up and down several times with a multi-channel pipettor.
16. Cover the plate and incubate for 15-30 minutes at room temperature.
17. OPTIONAL – Include the following steps if high backgrounds occur:

- a. Pre-wet a 1.2  $\mu\text{m}$  Millipore filter plate with 100  $\mu\text{L}$ /well of PBS-1% BSA and aspirate by vacuum manifold.
- b. Add 50  $\mu\text{L}$  of PBS-1% BSA to the appropriate wells of the filter plate.
- c. Transfer the contents of round-bottom plate to the filter plate.
- d. Aspirate the supernatant by vacuum manifold.
- e. Wash each well twice with 100  $\mu\text{L}$  of PBS-1% BSA and aspirate by vacuum manifold.
- f. Resuspend the microspheres in 100  $\mu\text{L}$  of PBS-1% BSA by gently pipetting up and down five times with a multi-channel pipettor.

18. Analyze 50-75  $\mu\text{L}$  on the Luminex analyzer according to the system manual.

**Technical Note 1:** Either PBS-1% BSA or PBS-BN (PBS, 1% BSA, 0.05% Azide, pH 7.4) may be used as Assay Buffer.

**Technical Note 2:** For a no-wash assay, the concentration of the detection antibody and the streptavidin-R-phycoerythrin should be optimized. The optimal concentrations tend to be higher than in a washed assay.