



~~~~~

## **Immunobead<sup>®</sup> Goat Anti-Rabbit Immunoglobulin**

**Catalog 170-5602**

**Control** 74163

**Contents** 50 mg of Immunobeads with covalently bound goat anti-rabbit immunoglobulin (H+L), lyophilized in 10 mM phosphate buffer, 500 mM sodium chloride, 5 mM EDTA, containing 1.0% BSA, 0.1% Tween-20, and 0.01% sodium azide.

**Capacity** 1.0 µg antibody/mg Immunobead

### **Instructions for Use**

1. Reconstitute Immunobead reagent with 5 ml deionized water to produce a working solution of 10 mg/ml. Further adjustments should be made using assay buffer following the guidelines recommended in step 2. The rehydrated reagent is stable for 6 months when stored at 4 °C. Aliquots may be frozen and stored at -20 °C. Avoid repeated freeze-thaw cycles.
2. This product contains 1.0% BSA as a carrier protein to minimize non-specific binding. Assay buffers should include BSA, be at least 0.15M NaCl, and at a pH between 6 and 8, to minimize non-specific interactions.
3. For most immunoassay procedures, incubation for 2 hours at 37 °C is recommended for complete binding to the solid phase reagent. If the assay is performed at lower temperatures or with dilute reaction volumes, these conditions may need to be extended. If the incubation period exceeds 2 hours, tubes should be mixed briefly at occasional intervals.
4. Centrifuge at 1,000 x g using at least 0.2 mg of Immunobead reagent per 12 x 75 mm tube, to produce a firm pellet. Decant as soon as centrifuge has stopped to avoid loss of the pellet by rapidly inverting the tube over a waste container and then blotting the rim gently on an absorbant towel.

LIT-408 REV B 792