



~~~~~

## **EIA Grade Affinity Purified Rabbit Anti-Goat IgG (H+L) Alkaline Phosphatase Conjugate**

**Catalog 172-1037**

|                                     |                                                                                                                                                                                                                                                                       |
|-------------------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Contents</b>                     | 1 ml                                                                                                                                                                                                                                                                  |
| <b>Reconstitution</b>               | The lyophilized contents of this vial should be reconstituted with 1 ml of sterile, 40% glycerol.                                                                                                                                                                     |
| <b>Buffer</b>                       | When reconstituted according to the instructions above, product contains 10 mM Tris, 250 mM sodium chloride, pH 8.0, 1.5% bovine serum albumin, and 0.05% sodium azide.                                                                                               |
| <b>Storage</b>                      | The unopened vial of lyophilized conjugate, as well as the conjugate when reconstituted as recommended above, can be stored at 4° C. If other methods of reconstitution and storage are employed, the use of azide and repeated freeze-thaw cycles should be avoided. |
| <b>Shelf life</b>                   | One year at 4 °C.                                                                                                                                                                                                                                                     |
| <b>Recommended working dilution</b> | 1:1,000 as determined by a direct ELISA method. Since assay conditions vary, the optimum dilution of this enzyme-antibody conjugate should be determined for each particular application.                                                                             |
| <b>Method of preparation</b>        | Antibody is conjugated to alkaline phosphatase using a modification of the glutaraldehyde procedure described by Engvall and Perlmann. <sup>1</sup>                                                                                                                   |
| <b>Reference</b>                    | 1. Engvall, E. and Perlmann, P., <i>Immunochemistry</i> , 8, 871 (1971).                                                                                                                                                                                              |

This product is intended for research use only. It is not intended for clinical diagnostic purposes.