

EIA Grade
Affinity Purified
Goat anti-Chicken IgG (H + L)
Alkaline Phosphatase Conjugate
Catalog No. 172-1025

| | |
|-------------------------------------|---|
| Contents | 0.1 mg antibody. |
| Reconstitution | The lyophilized contents of the vial should be reconstituted with 1 ml of sterile, 40% glycerol. |
| Storage | 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. Note that this product contains no preservatives. |
| Approximate working dilution | 1:1,000 as determined by a direct ELISA method. Since assay conditions vary, the optimum dilution of this antibody-enzyme conjugate should be determined for each particular application. |
| Method of preparation | Affinity purified antibodies isolated from a pool of sera from goats immunized with purified IgG from normal chicken sera, were conjugated to alkaline phosphatase using a modified glutaraldehyde procedure [A. Volker, D. Bidwell, and A. Bartlett, pp. 359-371 in Manual of Clinical Immunology , N. Rose and H. Friedman, eds., American Society for Microbiology, Washington, D.C. (1980)]. |

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