

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

<b>Contents</b>	0.1 mg antibody.
<b>Reconstitution</b>	The lyophilized contents of the vial should be reconstituted with 1 ml of sterile, 40% glycerol.
<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. Note that this product contains no preservatives.
<b>Approximate working dilution</b>	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.
<b>Method of preparation</b>	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 <b>Manual of Clinical Immunology</b> , 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

---