

**EIA Grade Affinity Purified  
Goat Anti-Human IgG (H + L)  
(Bovine Serum Adsorbed)  
Alkaline Phosphatase Conjugate  
Catalog No. 172-1004**

<b>Contents</b>	1.0 ml
<b>Preparation of antibody</b>	Goat anti-human <b>IgG</b> (H +L) is prepared from antisera raised in goats immunized with purified human <b>IgG</b> . Immunoaffinity chromatography procedures are used to isolate antibodies and to remove antibodies which <b>cross-react</b> with bovine immunoglobulin.
<b>Preparation of conjugate</b>	Antibody is conjugated to alkaline phosphatase using a modification of the one step glutaraldehyde procedure described by Engvall and Perlmann.'
<b>Buffer</b>	10 <b>mM</b> Tris, 150 <b>mM</b> sodium chloride, 1.0 <b>mM</b> magnesium chloride, <b>pH</b> 8.0. Contains 1.0% bovine serum albumin as a stabilizer and 0.1% sodium <b>azide</b> as a bacteriostat.
<b>Storage</b>	The product will arrive frozen. It may be stored frozen until needed. Once thawed, store at 4° C. Repeated freeze-thaw cycles should be avoided.
<b>Shelf life</b>	One year at 4° C.
<b>Recommended working dilution</b>	1: <b>3,000</b> as determined by a direct <b>ELISA</b> method. Since assay conditions vary, the optimum dilution should be determined for each particular application.  This product is intended for research use only. It is not intended for clinical diagnostic purposes.

---

**Reference** 1. Engvall, E. and Perlmann, P., *Immunochemistry*: 8, 871 (1971).