

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

Contents	1.0 ml
Preparation of antibody	Goat anti-human IgG (H +L) is prepared from antisera raised in goats immunized with purified human IgG . Immunoaffinity chromatography procedures are used to isolate antibodies and to remove antibodies which cross-react with bovine immunoglobulin.
Preparation of conjugate	Antibody is conjugated to alkaline phosphatase using a modification of the one step glutaraldehyde procedure described by Engvall and Perlmann.'
Buffer	10 mM Tris, 150 mM sodium chloride, 1.0 mM magnesium chloride, pH 8.0. Contains 1.0% bovine serum albumin as a stabilizer and 0.1% sodium azide as a bacteriostat.
Storage	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.
Shelf life	One year at 4° C.
Recommended working dilution	1: 3,000 as determined by a direct ELISA 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).