

EIA Grade
Affinity Purified
Goat anti-Rabbit IgG (H + L)
Alkaline Phosphatase Conjugate
Catalog No. 172-1016

Contents:	This package contains 1 vial containing 1.0 ml Antibody Solution.
Preparation of antibody	Goat anti-rabbit IgG (H + L) is prepared from antisera raised in goats immunized with purified rabbit IgG. Immunoaffinity chromatography procedures are used to isolate antibodies and to remove antibodies which cross react with human immunoglobulin.
Preparation of conjugate	Antibody is conjugated to alkaline phosphatase using a modification of the one step glutaraldehyde procedure described by Engvall and Perlmann. ¹
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 conditions	4° C for one year.
Recommended working dilution	1:3,000 as determined by direct ELISA. 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., <i>Immunochemistry</i> , 8,871 (1971).
------------------	---