



**Blotting Grade
Affinity Purified
Goat Anti-Rabbit IgG (H+L)
Alkaline Phosphatase
Conjugate**

**Catalog Number
170-6518**

Specifications

Contents	1.0 ml
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.
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 crossreact 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

Storage conditions

This reagent is shipped frozen on dry ice and can be stored at -20 °C prior to opening. Once thawed, store the reagent at 4 °C. Repeated freeze-thaw cycles will damage the reagent.

Shelf life

One year at 4 °C.

Recommended working dilution

1:3,000

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).