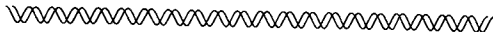


**EIA Grade Affinity
Purified Goat Anti-Mouse
IgG (H+L) Alkaline
Phosphatase Conjugate**

**Catalog Number
172-1015**

BIO-RAD



Bio-Rad Laboratories, 2000 Alfred Nobel Dr., Hercules, CA 94547
LIT-432 91-0808 0692

Specifications

Contents	1.0 ml
Preparation of antibody	Goat Anti-Mouse IgG (H+L) is prepared from antisera raised in goats immunized with purified 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

Goat Anti-Mouse IgG (H+L) AP conjugate 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

$\frac{1}{3,000}$ 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.

References

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