



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

**Catalog Number
172-1011**



Specifications

Contents	2.0 ml antibody solution.
Preparation of antibody	Goat Anti-Mouse IgG (H + L) is prepared from antisera raised in goats immunized with purified mouse 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 horseradish peroxidase using a modification of the periodate oxidation procedure described by Wilson and Nakane. ¹
Buffer	10 mM sodium phosphate, 150 mM sodium chloride, pH 7.2 containing 1.0% bovine serum albumin as a stabilizer and 0.01% thimerosal as a bacteriostat.

Storage conditions 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 1 year at 4 °C.

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. Wilson, M. B. and Nakane, P. K., Immunofluorescence and Related Staining Techniques, Knapp, W., Holubar, K. and Wick, G. eds. Elsevier/North Holland Biomedical Press, Amsterdam, 215-224 (1978).