



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

**Catalog Number  
170-6516**

***BIO-RAD***

# Specifications

---

<b>Contents</b>	2.0 ml
<b>Buffer</b>	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.
<b>Preparation of antibody</b>	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.
<b>Preparation of conjugate</b>	Antibody is conjugated to horseradish peroxidase using a modification of the periodate oxidation procedure described by Wilson and Nakane. <sup>1</sup>

**Storage** 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** 1 year at 4 °C.

**Recommended working dilution** 1:3,000 as determined by Western blotting procedures.

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, pp. 215-224 (1978).

***Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547***

LIT372 Rev D