



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

**Catalog Number
172-1050**

BIO-RAD

Specifications

Contents	2.0 ml
Buffer	10 mM sodium phosphate, 150 mM sodium chloride, pH 7.2. Contains 1.0% bovine serum albumin as a stabilizer and 0.01% thimerosal as a bacteriostat.
Preparation of antibody	Goat anti-human IgG (H + L) is prepared from antisera raised in goats immunized with purified human IgG. Immunoaffinity chromatography procedures are used to isolate antibodies and to remove antibodies which cross-react with bovine immunoglobulin.
Preparation of conjugate	Antibody is conjugated to horseradish peroxidase using a modification of the periodate oxidation procedure described by Wilson and Nakane. 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 product 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. 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).