



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

**Catalog Number  
170-6515**

***BIO-RAD***

# Specifications

<b>Contents</b>	2.0 ml
<b>Buffer</b>	10 mM sodium phosphate, 150 mM sodium chloride, pH 7.2. Contains 1.0% bovine serum albumin and 0.05% Proclin 300.
<b>Preparation of antibody</b>	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 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(1978).

**Storage**

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**

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

Wilson MB and Nakane PK(1978). *Immuno-fluorescence and Related Staining Techniques*, Knapp W, Holubar K, and Wick G, eds. (Amsterdam: Elsevier/North Holland Biomedical Press), pp. 215-244.

For further information please contact your local Bio-Rad representative, or in the U.S. dial 1-800-424-6723.