

Affinity Purified
Rabbit anti-Goat IgG (H + L)
Horseradish Peroxidase Conjugate
Catalog No. **172-1034**

- Contents:** 0.1 mg Antibody
- Reconstitution:** The lyophilized contents of the vial should be reconstituted with 1 ml of sterile, 40% glycerol.
- Storage:** The unopened vial of lyophilized conjugate, as well as the conjugate when reconstituted as recommended above, can be stored at 4°C. If other methods of reconstitution and storage are employed, the use of **azide** and repeated freeze-thaw cycles should be avoided. Note that this product contains no preservatives.
- Approximate Working Dilution:** 1:800 as determined by a direct **ELISA** method. Since assay conditions vary, the optimum **dilution of** this antibody-enzyme conjugate should be determined for each particular application.
- Method of Preparation:** Affinity-purified antibodies, isolated from a pool of sera from rabbits immunized with purified **IgG** from normal goat sera, were conjugated to horseradish peroxidase by the periodate oxidation method described by P. K. Nakane and A. Kawaoi [J. *Histochem. Cytochem.*, 22, 1084-1091 (1974)].

This product is intended for research use only. It is not intended for clinical diagnostic purposes.

86-0375 886

Bio-Rad Laboratories, 32nd & Griffin Ave., Richmond, CA 94804