

BIO-RAD

**EIA Grade
Affinity Purified
Goat anti-Human IgG (γ)
Horseradish Peroxidase Conjugate
Catalog No. 172-1033**

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	1:1,000 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 goats immunized with purified IgG from normal human sera, were conjugated to horseradish peroxidase by the periodate oxidation method described by P. K. Nakane and A. Kawaoi [<i>J. Histochem. Cytochem.</i> , 22, 1084-1091 (1974)].

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

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

87-0298 487