

BIO-RAD

Affinity Purified
Goat anti-Human IgA + IgG + IgM (H + L)
Alkaline Phosphatase Conjugate

Catalog No. 172-1046

Control No. 26272

- 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, repeated freeze-thaw cycles should be avoided. Note that this product contains no preservatives.
- Approximate Working Dilution:** 1:400 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 immunoglobulins from normal human sera, were conjugated to alkaline phosphatase using a modified gluteraldehyde procedure [A. Volker, D. Bidwell, and A. Bartlett, pp. 359371 in *Manual of Clinical Immunology*, N. R. Rose and H. Friedman, eds., American Society for Microbiology, Washington, D.C. (1980)].

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