

Affinity Purified
*Goat anti-Human **IgA + IgG + IgM (H + L)***
*Horseradish **Peroxidase** Conjugate*

Catalog No. 172-1049

Control No. 26213

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 antibodyenzyme 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 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.