

**EIA Grade
Affinity Purified
Goat anti-Mouse *IgM* (μ)
Alkaline Phosphatase Conjugate
Catalog No. 172-1028**

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 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 IgM from normal mouse sera, were conjugated to alkaline phosphatase using a modified gluteraldehyde procedure [A. Volker, D. Bidwell, and A. Bartlett, pp. 359-371 in Manual of Clinical Immunology , N. Rose and H. Friedman, eds., American Society for Microbiology, Washington, DC. (19&I)].

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