

**EIA Grade
Affinity Purified
Goat anti-Mouse *IgM* (μ)
Horseradish Peroxidase Conjugate
Catalog No. 172-1030**

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 IgM from normal mouse sera , were conjugated to horseradish peroxidase by the periodate oxidation method described by I? 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.
