

EIA Grade
Protein A - Horseradish Peroxidase Conjugate
Catalog No. 172-1018

Contents	0.5 mg of Protein A.
Reconstitution	The lyophilized contents of the vial should be reconstituted with 1 ml of sterile, distilled/deionized water. When reconstituted in this fashion, the conjugate will be contained in PBS buffer pH 7.4 with 1% BSA and 0.01% Thimerosal as a preservative.
Storage	The unopened vial of lyophilized conjugate, as well as the conjugate when constituted as recommended above, can be stored at 2-5° C.
Approximate working dilution	13,000. Since assay conditions vary, the optimum dilution of this product should be determined for each particular application.
Method of preparation	A 1:1 mixture of Protein A and Horseradish Peroxidase were conjugated by a modification of the method of M. B. Wilson and P. K. Nakane. ¹

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

Reference	1. Wilson, M. B. and Nakane, P. K., [Immunofluorescence and Related Staining Techniques, (Knapp, W., Holubar, K. and Wicks, G., eds.) Elsevier/North Holland, New York, pgs. 215-224 (1978)].
------------------	---