



Blotting Grade Protein A - Horseradish Peroxidase Conjugate

**Catalog Numbers
170-6522**

BIO-RAD

Specifications

Contents	1.0 ml
Buffer	10 mM sodium phosphate, 150 mM sodium chloride, pH 7.2, containing 1.0% bovine serum albumin as a stabilizer and 0.01% thimerosal as a bacteriostat.
Preparation of conjugate	A 1:1 mixture of protein A and horseradish peroxidase was conjugated using a modification of the periodate procedure described by Wilson and Nakane. ¹
Storage	The product will arrive frozen. It may be stored frozen until needed. Once thawed, store at 4 °C. Repeated freeze-thaw cycles should be avoided.
Shelf life	One year at 4 °C.

**Recommended 1/3,000
working
dilution**

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 Wick, G., eds., Elsevier/North Holland Biomedical Press, Amsterdam, pp. 215-244 (1978).