

Qty: 3 mg/2 ml Streptavidin-AP Conjugate **Catalog No.** 43-4322

Lot No. See product label

AP-Streptavidin Conjugate

FORM

Streptavidin-AP is supplied as 2ml of liquid at 1.5mg/ml in a proprietary enzyme-conjugate stabilizing buffer containing 0.1% sodium azide as a preservative. In the preparation of this product, Zymed uses highly purified streptavidin and alkaline phosphatase.

BACKGROUND

Streptavidin (MW \equiv 66 kDa) binds specifically with biotin (244 Da). It is derived from the bacterium *Streptomyces avidinii* and bears a remarkable similarity to chicken egg-white avidin both in three-dimensional structure and its ability to bind biotin with extremely high affinity (K_d=10⁻¹⁵M). It is a tetrameric protein capable of binding up to 4 biotin molecules. Unlike avidin, Streptavidin is non-glycosylated and is essentially neutral in charge, whereas avidin (pl \cong 10.5) is basic at neutral pH. Because of this, streptavidin has considerably less non-specific binding resulting in less background. It has replaced avidin as the reagent of choice for most applications where protein interactions may cause background.

USAGE

Zymed's Streptavidin-AP is recommended for use in detection systems utilizing biotinylated antibodies, and other biotinylated molecules. This product can be used for a variety of common applications including immunohistochemistry, Western blotting, *In Situ* hybridization, and ELISA. For an excellent, comprehensive review of properties and applications of streptavidin/biotin amplification methods, see Bayer and Wilchek's <u>Avidin-Biotin Technology</u>, Methods in Enzymology, Volume 184, Academic Press (1990).

Working concentrations for specific applications should be determined by the investigator. Appropriate dilutions will be affected by several factors, including primary and secondary antibody affinity, antigen concentration and length of incubations. We recommend the following ranges as starting points.

Immunohistochemistry ⁽¹⁻³⁾ :	1:100 to 1:300
Western blot (chromogenic) ⁽⁴⁾ :	1:500 to 1:1,250
	1:1,000 to 1:2,500
ELISA ⁽⁵⁻⁷⁾ :	1:1,000 to 1:2,500

PROCEDURES

Zymed has general guidelines for ELISA, blotting and other applications available on our Web site at <u>www.invitrogen.com/methods</u>. You may also obtain assistance from our Technical Service department at (800) 874-4494. Another good source of information about general immunoassay procedures is Ed Harlow & David Lane's <u>Antibodies, A</u> <u>Laboratory Manual</u>, Cold Spring Harbor Laboratory (1988). Also see Bayer and Wilchek (referenced above).

(Rev 12/08) DCC-08-1089

PI434322

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, <u>www.invitrogen.com</u>). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

(Rev 12/08) DCC-08-1089

STORAGE AND STABILITY

Store at 2-8°C.

WARRANTY

Zymed products are guaranteed to perform as stated for the recommended applications. This warranty is valid until the expiration date printed on the bottle.

SELECTED PRODUCT REFERENCES

References cite Zymed's standard products.

Immunohistochemistry & Immunocytochemistry

- 1. Greenwald, R.J. et al, J Immunol. 158:4088-4096 (1997).
- 2. Nirenburg, S. and Cepko, C., J Neurosci. 13(8):3238-3251 (1993).
- 3. Yang, X. et al., J Neurosci. 13(7):3006-3017 (1993).

Western blot

4. Zioncheck, T.F. et al, J Biol. Chem. 270(28):16871-16878 (1995).

ELISA

5. Deng, G. et al, *J Cell Biol.* 134(6):1563-1571 (1996).

- 6. Yoshida, A. et al, *J Immunol.* 155:2057-2066 (1995).
- 7. Thomas, A.P. et al, Hybridoma 15(5):359-364(1996).

In Situ Hybridizaion

10. McQuid, S. and Allan, G., J. Histochem. Cytochem. 40:569-574 (1992)

RELATED PRODUCTS

Product	Cat. No.
Goat anti-Mouse IgG (H+L)-Biotin (ZyMAX™ Grade)	81-6540
Goat anti Rabbit IgG (H+L)-Biotin (ZyMAX™ Grade)	81-6140
Rabbit anti-Goat IgG (H+L)-Biotin (ZyMAX™ Grade)	81-1640
Goat anti-Human IgG (H+L)-Biotin (ZyMAX™ Grade)	81-7140
Goat anti-Rat IgG (H+L)-Biotin	62-9540
pNPP Single Solution (chromogen for ELISA) BCIP/NBT Reagent Kit (chromagen for immunoblotting) BCIP/NBT Liquid Reagent Kit (chromogen for IHC, blotting) AP-Fast Red Reagent (chromagen for IHC) AP-Blue Reagent Kit (chromogen for IHC)	00-2212 (100 ml), 00-2213 (500 ml) 00-2210 00-2211 (IHC), 00-2209 (blot) 00-2234 00-2204

TRADEMARKS

PI434322

Zymed[®] is a trademark of Zymed Laboratories, Inc.

For research use only.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com