

Qty: 1.5 mg/1 mL
ZyMax™
Goat anti-Rabbit IgG (H+L)
Biotin Conjugate



Catalog No. 81-6140
Lot No. See product label

Biotin-Goat Anti-Rabbit IgG (H+L) Conjugate (ZyMax™ Grade)

FORM

This ZyMax™ antibody conjugate is supplied in liquid form and contains 1.5 mg/mL of conjugate in 10 mM phosphate-0.15 M NaCl buffer, pH 7.4, containing 1% bovine serum albumin and 0.1% sodium azide. Purified antibody is biotinylated under optimal conditions and free biotin is removed from the preparation.

SPECIFICITY

Prior to conjugation, this polyclonal antibody is raised in goats against the whole rabbit IgG molecule and purified by antigen-affinity chromatography. It reacts strongly with rabbit IgG heavy and light chains. This antibody has been absorbed with human serum proteins so reactivity with human IgG is minimal or absent.

USAGE

ZyMax™ grade reagents offer enhanced assay sensitivity, higher titers, and greater flexibility in assay design, than standard grade reagents. This biotin conjugate is recommended as a secondary reagent for ELISA, Western blotting, immunohistochemistry, and other applications. Working concentrations for specific applications should be determined by the investigator. Optimal dilutions will be affected by several factors, including: primary antibody affinity, sensitivity of detection method, antigen concentration, temperature and length of incubations, etc. We recommend the following ranges as starting points.

ELISA 1:10,000 to 1:30,000 (in PBS containing 0.05% Tween 20)
Western blot: 1:5,000 to 1:10,000
Immunohistochemistry: 1:100 to 1:300

PROCEDURES

Invitrogen has recommended procedures for ELISA, immunoprecipitation and other applications available on our Web site at www.invitrogen.com. You may also obtain assistance from our Technical Service department at (800) 955.6288. Another good source of information about immunoassay procedures is Ed Harlow & David Lane's Antibodies, A Laboratory Manual, Cold Spring Harbor Laboratory (1988).

STORAGE

Store at 2-8°C. This product is guaranteed to perform as stated if used before the expiration date printed on the bottle.

SELECTED INVITROGEN PRODUCT REFERENCES

References cite Invitrogen's standard grade products--ZyMax™ antibodies may be used at considerably higher dilutions.

Immunohistochemistry

1. Boag, A.H. & Young, I.D.: *Amer. J. Pathology* 144(3):585 (1994).
2. Grone, A. et al: *Vet. Pathology* 31:308-315 (1994).
3. Boag, A.H. & Young, I.D.: *Modern Pathology* 6:65 (1993).

ELISA

4. Seiffert, D., et al. *J. Biol. Chem.* 269(31): 19836-19842 (1994)

(81-6140 continued)

www.invitrogen.com

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

PI816140

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). 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.

RELATED PRODUCTS

Product	Conjugate	Cat. No.
Goat anti-Rabbit IgG (H+L) (ZyMax™ Grade)	Purified	81-6100
	FITC	81-6111
	TRITC	81-6114
	Cy™3	81-6115
	Cy™5	81-6116
	HRP	81-6120
	AP	81-6122
Goat F(ab') ₂ anti-Rabbit IgG (H+L)	Purified	62-1800
	FITC	62-1811
	HRP	62-1820
	AP	62-1822
	Biotin	62-1840
Streptavidin (ZyMax™ Grade)	HRP	43-8123
	AP	43-8122
Streptavidin	FITC	43-4311
	TRITC	43-4314
	Cy3	43-4315
	Cy5	43-4316
	Texas Red®	43-4317
	Phycoerythrin	15-4301
	Sepharose® 4B	43-4341

NOTES

Zymed® and ZyMax™ are trademarks of Zymed Laboratories, Inc. Sepharose® and Cy™ are trademarks of Amersham Biosciences Limited.

For Research Use Only

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: techsupport@invitrogen.com

PI816140

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). 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.