



61-3140 2 ml
Biotin-Rabbit Anti-
Chicken/Turkey IgG (H+L)

Biotin-Rabbit Anti-Chicken/Turkey IgG (H+L)

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

FORM: Liquid. The product is in 10 mM phosphate-0.15 M NaCl buffer, pH 7.4, containing 1% bovine serum albumin, 0.1% sodium azide, and 50% glycerol. The antibody concentration is 0.75 mg/ml.

AFFINITY PURIFIED RABBIT ANTI-CHICKEN/TURKEY IgG (H+L) SPECIFIC ANTIBODY is biotinylated under optimal conditions and is free of unconjugated biotin.

SUGGESTED WORKING DILUTION: 1:8,000 to 1:10,000 in ELISA using peroxidase streptavidin as a probe. Working dilution for specific application should be determined by the investigator to obtain best results.

STORAGE: Store at 4°C.

REMARKS: For research use only. Not for human or drug use. Sodium azide at 0.1% has been added as preservative. This product has not been assayed for species cross-reactivity unless otherwise indicated.

Research Use Only

www.invitrogen.com

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

PI613140

(Rev 12/08) 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.



61-3140 2 ml
Biotin-Rabbit Anti-
Chicken/Turkey IgG (H+L)

Biotin-Rabbit Anti-Chicken/Turkey IgG (H+L)

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

FORM: Liquid. The product is in 10 mM phosphate-0.15 M NaCl buffer, pH 7.4, containing 1% bovine serum albumin, 0.1% sodium azide, and 50% glycerol. The antibody concentration is 0.75 mg/ml.

AFFINITY PURIFIED RABBIT ANTI-CHICKEN/TURKEY IgG (H+L) SPECIFIC ANTIBODY is biotinylated under optimal conditions and is free of unconjugated biotin.

SUGGESTED WORKING DILUTION: 1:8,000 to 1:10,000 in ELISA using peroxidase streptavidin as a probe. Working dilution for specific application should be determined by the investigator to obtain best results.

STORAGE: Store at 4°C.

REMARKS: For research use only. Not for human or drug use. Sodium azide at 0.1% has been added as preservative. This product has not been assayed for species cross-reactivity unless otherwise indicated.

Research Use Only

www.invitrogen.com

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

PI613140

(Rev 12/08) 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.