



61-8620 1 ml
HRP-Rabbit Anti-
Sheep IgG (H+L)
Lot No.

Peroxidase-Rabbit Anti-Sheep IgG (H+L)

For Research Use Only. Not for use in diagnostic procedures.

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

Form: 1 ml liquid. The product is in peroxidase stabilizing buffer containing 0.1% proclin as preservative.

AFFINITY PURIFIED RABBIT ANTI-SHEEP IgG (H+L) SPECIFIC ANTIBODY is conjugated with highly purified horseradish peroxidase (RZ > 3.0).

RECOMMENDED WORKING DILUTIONS FOR ELISA: 1:2,000 to 1:4,000. Working dilution for specific application should be determined by the investigator to obtain the best conditions.

STORAGE: Store at 4°C.

REMARKS: For research use only. Not for human or drug use. Product contains 0.1% proclin 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

PI618620

(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.



61-8620 1 ml
HRP-Rabbit Anti-
Sheep IgG (H+L)
Lot No.

Peroxidase-Rabbit Anti-Sheep IgG (H+L)

For Research Use Only. Not for use in diagnostic procedures.

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

Form: 1 ml liquid. The product is in peroxidase stabilizing buffer containing 0.1% proclin as preservative.

AFFINITY PURIFIED RABBIT ANTI-SHEEP IgG (H+L) SPECIFIC ANTIBODY is conjugated with highly purified horseradish peroxidase (RZ > 3.0).

RECOMMENDED WORKING DILUTIONS FOR ELISA: 1:2,000 to 1:4,000. Working dilution for specific application should be determined by the investigator to obtain the best conditions.

STORAGE: Store at 4°C.

REMARKS: For research use only. Not for human or drug use. Product contains 0.1% proclin 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

PI618620

(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.