



61-9500 2 ml
Rabbit Anti-
Rat IgG (H+L)
Lot No. See product label

RABBIT ANTI-RAT IgG (H+L)

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

FORM: Lyophilized powder. Reconstitute with 2 ml of distilled water. **AFFINITY PURIFIED RABBIT ANTI-RAT IgG SPECIFIC ANTIBODY** is in 10 mM phosphate buffered saline, pH 7.4, containing 0.1% sodium azide as preservative. The protein concentration is 1 mg per ml determined by $E(1\%) = 14.0$ at 280 nm.

Antibody is produced in rabbits against rat IgG and is affinity purified through rat IgG column. The antibody has been absorbed with human serum proteins.

STORAGE: Store at 4°C. After reconstitution, store frozen in small aliquots. Avoid repeated freezing and thawing.

REMARKS: For research use only. Not for human or drug use. Product contains 0.1% sodium azide as preservative.

Research Use Only

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PI619500

(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.



61-9500 2 ml
Rabbit Anti-
Rat IgG (H+L)
Lot No. See product label

RABBIT ANTI-RAT IgG (H+L)

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

FORM: Lyophilized powder. Reconstitute with 2 ml of distilled water. **AFFINITY PURIFIED RABBIT ANTI-RAT IgG SPECIFIC ANTIBODY** is in 10 mM phosphate buffered saline, pH 7.4, containing 0.1% sodium azide as preservative. The protein concentration is 1 mg per ml determined by $E(1\%) = 14.0$ at 280 nm.

Antibody is produced in rabbits against rat IgG and is affinity purified through rat IgG column. The antibody has been absorbed with human serum proteins.

STORAGE: Store at 4°C. After reconstitution, store frozen in small aliquots. Avoid repeated freezing and thawing.

REMARKS: For research use only. Not for human or drug use. Product contains 0.1% sodium azide as preservative.

Research Use Only

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PI619500

(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.