



61-0320 1 ml  
HRP-Rabbit Anti-  
Mouse IgG2b  
Lot No.

**Peroxidase-Rabbit Anti-Mouse IgG2b (Gamma 2b Specific)**

**INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY**  
MINIMUM CROSS-REACTIVITY TO HUMAN SERUM PROTEIN

**FORM:** 1 ml liquid. The product is in enzyme conjugate stabilizer and 0.1% Proclin as preservative.

AFINITY PURIFIED RABBIT ANTI-MOUSE IgG2b SPECIFIC ANTIBODY is conjugated with highly purified horseradish peroxidase (RZ > 3.0).

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

**STORAGE:** Store at 4°C.

**Research Use Only**

[www.invitrogen.com](http://www.invitrogen.com)

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

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI610320

(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](http://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.



61-0320 1 ml  
HRP-Rabbit Anti-  
Mouse IgG2b  
Lot No.

**Peroxidase-Rabbit Anti-Mouse IgG2b (Gamma 2b Specific)**

**INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY**  
MINIMUM CROSS-REACTIVITY TO HUMAN SERUM PROTEIN

**FORM:** 1 ml liquid. The product is in enzyme conjugate stabilizer and 0.1% Proclin as preservative.

AFINITY PURIFIED RABBIT ANTI-MOUSE IgG2b SPECIFIC ANTIBODY is conjugated with highly purified horseradish peroxidase (RZ > 3.0).

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

**STORAGE:** Store at 4°C.

**Research Use Only**

[www.invitrogen.com](http://www.invitrogen.com)

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

E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI610320

(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](http://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.