



62-7540 1 ml
Biotin-Goat Anti-
Human IgM

Biotinylated Goat Anti-Human IgM (μ ch. sp.)

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

FORM: Liquid in 0.01 M phosphate- 0.15 M NaCl, 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 GOAT ANTI-HUMAN IgM (μ chain) SPECIFIC ANTIBODY is biotinylated under optimum conditions and is free of unconjugated biotin.

SUGGESTED DILUTION FOR EIA: 1:8,000 to 1:10,000 with 0.05% Tween 20 in PBS. Working dilution for specific application should be determined by the investigator to obtain the best results.

STORAGE: Store at 4°C.

REMARKS: For research use only. Not for human or drug use. Product contains 0.05% 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

PI627540

(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.



62-7540 1 ml
Biotin-Goat Anti-
Human IgM

Biotinylated Goat Anti-Human IgM (μ ch. sp.)

INVITROGEN AFFINITY PURIFIED ANTIGEN SPECIFIC ANTIBODY

FORM: Liquid in 0.01 M phosphate- 0.15 M NaCl, 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 GOAT ANTI-HUMAN IgM (μ chain) SPECIFIC ANTIBODY is biotinylated under optimum conditions and is free of unconjugated biotin.

SUGGESTED DILUTION FOR EIA: 1:8,000 to 1:10,000 with 0.05% Tween 20 in PBS. Working dilution for specific application should be determined by the investigator to obtain the best results.

STORAGE: Store at 4°C.

REMARKS: For research use only. Not for human or drug use. Product contains 0.05% 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

PI627540

(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.