



03-9820 1 ml
HRP-Mouse Anti-
Rat IgM

PEROXIDASE-MONOCLONAL MOUSE ANTI-RAT IgM

INVITROGEN MOUSE MONOCLONAL ANTIBODY REAGENT

FORM:

Liquid. The product is in enzyme conjugate stabilizer containing 0.1% proclin as preservative.

PURIFIED MONOCLONAL MOUSE ANTI-RAT IgM SPECIFIC ANTIBODY is conjugated with highly purified horseradish peroxidase (RZ > 3.0).

RECOMMENDED WORKING DILUTIONS FOR ELISA:

1:1,000. Working dilution for specific application should be determined by the investigator to obtain the best condition.

STORAGE:

Store at 4°C.



03-9820 1 ml
HRP-Mouse Anti-
Rat IgM

PEROXIDASE-MONOCLONAL MOUSE ANTI-RAT IgM

INVITROGEN MOUSE MONOCLONAL ANTIBODY REAGENT

FORM:

Liquid. The product is in enzyme conjugate stabilizer containing 0.1% proclin as preservative.

PURIFIED MONOCLONAL MOUSE ANTI-RAT IgM SPECIFIC ANTIBODY is conjugated with highly purified horseradish peroxidase (RZ > 3.0).

RECOMMENDED WORKING DILUTIONS FOR ELISA:

1:1,000. Working dilution for specific application should be determined by the investigator to obtain the best condition.

STORAGE:

Store at 4°C.

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PI039820

(Rev 09/09) DCC-09-1334

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.

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PI039820

(Rev 09/09) DCC-09-1334

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.