



Qty: 100 µg/200 µL

Mouse anti-ApoA5

Catalog No. 39-8300

Lot No.

Mouse anti-ApoA5

FORM

This monoclonal antibody is supplied as a 200 µL aliquot at a concentration of 0.5 mg/mL in PBS, pH 7.4, containing 0.1% sodium azide. This antibody is highly purified from mouse ascites by protein A chromatography.

CLONE: ZA010

ISOTYPE: Mouse IgG1-kappa

IMMUNOGEN

Recombinant protein corresponding to the full-length human ApoA5 protein

SPECIFICITY

This antibody is specific for the ApoA5 (apolipoprotein A5, ApoA-V, regeneration associated protein 3 (RAP3)) protein. On Western blots, it identifies the target band at ~41 kDa.

REACTIVITY

Reactivity has been confirmed with human HepG2 cell lysates.

Sample	ELISA	Western Blotting
Human	ND	+++
Immunogen	+++	+++

(Excellent +++, Good++, Poor +, No reactivity 0, Not applicable N/A, Not Determined ND)

USAGE

Working concentrations for specific applications should be determined by the investigator. Appropriate concentrations will be affected by several factors, including secondary antibody affinity, antigen concentration, sensitivity of detection method, temperature and length of incubations, etc. The suitability of this antibody for applications other than those listed below has not been determined. The following concentration ranges are recommended starting points for this product.

ELISA: 0.1-1.0 µg/mL
Western Blotting: 1 µg/mL

STORAGE

Store at 2-8°C for up to one month. Store at -20°C for long-term storage. Avoid repeated freezing and thawing.

(cont'd)

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI398300

(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.

BACKGROUND

Apolipoprotein A5 (ApoA5, ApoA-V, regeneration associated protein 3 (RAP3)) is an apolipoprotein that is an important determinant of plasma triglyceride levels, a major risk factor for coronary artery disease.¹⁻³ It is a component of high density lipoprotein and is highly similar to a rat protein that is upregulated in response to liver injury.¹ ApoA5 is present in human serum although at much lower concentration than other apolipoproteins.¹

REFERENCES

1. van Der Vliet HN, et al. *J Biol Chem* 276:44512-44520, 2001.
2. Weinberg RB, et al. *J Biol Chem* 278:34438-34444, 2003.
3. Pennacchio LA, et al. *Science* 294:169-173, 2001.
4. O'Brien PJ, et al. *Clin Chem* 51:351-359, 2005.

RELATED PRODUCTS

Product	Conjugate	Cat. No.
Protein A	Sepharose® 4B	10-1041
rec-Protein G	Sepharose® 4B	10-1241

Conjugate	ZyMAX™ Goat x Rabbit IgG (H+L)	ZyMAX™ Goat x Mouse IgG (H+L)
Purified	81-6100	81-6500
FITC	81-6111	81-6511
TRITC	81-6114	81-6514
Cy™3	81-6115	81-6515
Cy™5	81-6116	81-6516
HRP	81-6120	81-6520
AP	81-6122	81-6522
Biotin	81-6140	81-6540

Zymed® and ZyMAX™ are trademarks of Zymed Laboratories Inc. Cy™ and Sepharose® are registered trademarks of Amersham Biosciences Ltd.

For Research Use Only

MZ060214

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI398300

(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.