



Qty: 100 µg/200 µL

Mouse anti-AIB1

Catalog No. 39-7900

Lot No.

Mouse anti-AIB1

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: ZA008

ISOTYPE: Mouse IgG₁-kappa

IMMUNOGEN

Recombinant protein derived from the N-terminal region of the human AIB1 protein, which is 91% homologous to mouse and rat

SPECIFICITY

This antibody is specific for the AIB1 (amplified in breast cancer 1, nuclear receptor coactivator 3 (NCOA3), ACTR, CAGH16, CBP-interacting protein, CTG26, pCIP, RAC3) protein. On Western blots, it identifies the target band at ~110-150 kDa, depending on the experimental sample.

REACTIVITY

Reactivity has been confirmed with human MCF-7 cell lysates and with mouse lymph node and rat skeletal muscle homogenates.

Sample	ELISA	Western Blotting
Human	ND	+++
Mouse	ND	+++
Rat	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

PI397900

(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

AIB1 (amplified in breast cancer 1, nuclear receptor coactivator 3 (NCOA3), ACTR, CAGH16, CBP-interacting protein, CTG26, pCIP, RAC3) is a nuclear receptor coactivator that interacts with nuclear hormone receptors to enhance their transcriptional activator functions.¹ The encoded protein has histone acetyltransferase activity and recruits p300/CBP-associated factor and CREB binding protein as part of a multisubunit coactivation complex. AIB1 is amplified and overexpressed in breast and ovarian cancer cell lines as well as in breast cancer biopsies, and it also interacts with estrogen receptors in a ligand-dependent fashion.^{2,3} AIB1 also functions to enhance estrogen-dependent transcription (2-3).^{2,3} This protein is initially found in the cytoplasm but is translocated into the nucleus upon phosphorylation. Two transcript variants encoding different isoforms have been found for the *AIB1* gene.

REFERENCES

1. Anzick SL, et al. *Science* 277:965-968, 1997.
2. Bai J, et al. *Cell* 103: 1047-1058, 2000.
3. Shirazi SK, et al. *Clin Genet* 54: 102-103, 1998.

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

MZ060111

www.invitrogen.com

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

PI397900

(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.