



Qty: 100 µg/400 µL

Rabbit anti-AIDA-1

Catalog No. 36-7000

Lot No.

Rabbit anti-AIDA-1

FORM

This polyclonal antibody is supplied as a 400 µL aliquot at a concentration of 0.25 mg/mL in PBS, pH 7.4, containing 0.1% sodium azide. The antibody is epitope-affinity-purified from rabbit antiserum.

PAD: ZMD.317

IMMUNOGEN

Synthetic peptide derived from the C-terminal region of the human AIDA-1 (EB-1, AT13, E2A-PBX1- associated protein, Cajalin-2) protein.

SPECIFICITY

This antibody is specific for human AIDA-1. On Western blots, it identifies a band at ~150 kDa in pre-B Acute Lymphoblastic Leukemia (ALL) cell line 697.

REACTIVITY

Reactivity has been confirmed with 697 ALL cell lysates. Nalm-6 or Reh Pre-B ALL cell lysates may be used as negative controls. Cross-reactivity with rat AIDA-1 is expected based on amino acid sequence homology.

Sample	Immuno-Precipitation (native)	Western Blotting	ELISA
Human	+++	+++	N/A
Immunogen	N/A	N/A	+++

(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.

Western Blotting: 1-3 µg/mL

ELISA: 0.1-1.0 µg/mL

Immunoprecipitation: 7 µg

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

PI367000

(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

Amyloid-beta precursor protein (β -APP) intracellular domain associated protein -1 (AIDA-1), which was previously named EB-1¹, is a new PTB-containing protein that interacts with the cytoplasmic tail of β -APP. Three isoforms of AIDA-1, AIDA-1a, AIDA-1b and AIDA-1b Δ Ank², were cloned from human fetal brain². Other names include AT13, E2A-PBX1-associated protein, and Cajalin-2. AIDA-1 proteins are expressed at high levels in the brain, and studying their involvement in β -APP processing and AID function may give new insights and may show a possible role for these molecules in normal brain development and in the pathogenesis of Alzheimer's disease². AIDA-1 proteins and β -APP interact *in vitro*, in living cells and, endogenously in leukemia cell lines. The pre-B Acute Lymphoblastic Leukemia (ALL) cell line 697 specifically expresses the longest form (AIDA 1b), while this form is completely absent in other pre-B ALL cell lines such as Nalm-6 and Reh².

REFERENCES

1. Fu X, et al. *Oncogene* 18:4920-4929, 1999.
2. Ghersi E, et al. *J Alzheimers Dis*, 2003. (In press).

RELATED PRODUCTS

Product	Clone or PAD*	Cat. No.
Rb anti-Presenilin 1 (N-term)	PS-1N	71-1300
Rb anti-Presenilin 1 (N-term)	ZMD.97	34-4600
Rb anti-Presenilin 1 (Loop)	PS-1L	51-4200
Rb anti-Presenilin 2 (N-term)	ZMD.96	34-4500
Rb anti-Presenilin 2 (Loop)	ZMD.84	34-4400
Rb anti-Nicastrin	ZMD.242	34-9200
Rb anti-Amyloid- β Precursor Protein	ZMD.316	36-6900
Ms anti-Amyloid- β Precursor Protein	LN27	13-0200
Rb anti-Amyloid- β Peptide	Polyclonal	71-5800
Ms anti-Amyloid- β Peptide	AMY-33	13-0100
Rb anti-BACE	ZMD.116	34-4900
Rb anti-BACE 2 (C-term)	ZMD.117	34-5000
Rb anti-BACE 2 (Mid)	ZMD.167	34-5100
Protein A	Sepharose® 4B	10-1041
rec-Protein G	Sepharose® 4B	10-1241

*PAD: Polyclonal Antibody Designation

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.

Leej120403

For Research Use Only

www.invitrogen.com

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

PI367000

(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