



Qty: 100 µg/400 µl

Rabbit anti-Angiopoietin 2

Catalog No. 51-4100

Lot No. See product label

## Rabbit anti-Angiopoietin 2

### FORM

This polyclonal antibody is supplied as a 400 µl aliquot at 0.25 mg/ml in phosphate buffered saline (PBS), pH 7.4, containing 0.1% sodium azide (NaN<sub>3</sub>). The antibody is epitope-affinity purified from rabbit antiserum.

**ISOTYPE:** Rabbit Ig

**POLYCLONAL ANTIBODY DESIGNATION (PAD):** NE8

### IMMUNOGEN

A synthetic peptide derived from the amino-terminus of human angiopoietin 2.

### SPECIFICITY

This antibody is specific for the ~ 70 kDa Angiopoietin 2 protein. Reactivity of the antibody was confirmed by Western blotting using lysates from bovine aortic endothelial (BAEC) cells. This antibody does not appear to cross-react with any other related endogenous proteins. Positive control: Lysates from BAEC and HuVEC cells.

### REACTIVITY

Sample	ELISA	Western Blot
Human		+++
Bovine		+++
Immunogen	+++	

### 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:** 1-3 µg/ml

The suitability of this antibody for applications other than those listed above has not been determined.

### 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](http://www.invitrogen.com)

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

PI514100

(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](http://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**

Angiopoietin-1 (Ang1) and Angiopoietin-2 (Ang2) are ligands that regulate tyrosine phosphorylation of the Tie2/Tek receptor on endothelial cells. Proper regulation of Tie2/Tek is required for normal vascular development, by regulating vascular remodeling and endothelial cell interactions.<sup>(3)</sup> Angiopoietin-2 (Ang2) is an antagonist of Ang1 that competes for binding of the Tie2 receptor and blocks Ang-1 induced Tie2 phosphorylation during vasculogenesis.<sup>(1,2)</sup>

**REFERENCES**

1. Maisonpierre, P.C. et. al. (1997) Science 277; 55-60.
2. Kim, I. et. al. (2000) J Biol Chem 275; 18550-6.
3. Zagzag D. et. al. (1999) Exp Neurol 159; 391-400.

**RELATED PRODUCTS**

<b>Primary antibodies</b>	<b>PAD</b>	<b>Cat. No.</b>
Rabbit anti-CXCR4 + Peptide	CXCR4	52-1107
Rabbit anti-MetAP-2	CM33	71-7200
Rabbit anti-Neuropilin-1 + Peptide	Z-CN1	52-0107
Rabbit anti-Neuropilin-2 + Peptide	Z-CN2	52-0207
Rabbit anti-PIGF-1 + Peptide	Z-CP1	52-0307
Rabbit anti-PIGF-2 + Peptide	Z-CP2	52-0407
Rabbit anti-VEGF (NT) + Peptide	Z-NVF8	52-0507
Rabbit anti-VEGF (CT)	Z-CVF3	18-0254
Rabbit anti-VEGF-B + Peptide	Z-CVB2	52-0707
Rabbit anti-VEGF-C	Z-CVC7	18-0255
Rabbit anti-VEG1 + Peptide	Z-CV12	52-0907
Rabbit anti-Vezf1 + Peptide	Z-CVZ1	52-2007

PAD: Polyclonal Antibody Designation

<b>Immunoassay reagents</b>	<b>Conjugate</b>	<b>Cat. No.</b>
Goat anti-Rabbit IgG (H+L) (ZyMAX™ Grade)	Purified	81-6100
	FITC	81-6111
	TRITC	81-6114
	Cy™3	81-6115
	Cy™5	81-6116
	HRP	81-6120
	AP	81-6122
	Biotin	81-6140
Protein A	Sepharose® 4B	10-1041
rec-Protein G	Sepharose® 4B	10-1241

ZyMed® and ZyMAX™ are trademarks of Zymed Laboratories Inc. Cy™ is a trademark of Amersham Life Sciences, Inc. Sepharose® is a registered trademark of Pharmacia LKB.

**For Research Use Only**

SS000919

[www.invitrogen.com](http://www.invitrogen.com)

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

PI514100

(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](http://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.