



Rabbit (polyclonal) Anti-Tie2 [pY¹¹⁰⁸] Phosphospecific Antibody, Unconjugated

PRODUCT ANALYSIS SHEET

Catalog Number:	44-1313G (10 mini-blot size)
Lot Number:	See product label
Volume:	100 μ L
Form of Antibody:	Rabbit polyclonal immunoglobulin in Dulbecco's phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.3 (+/- 0.1), 50% glycerol, with 1.0 mg/mL BSA (IgG, protease free) as a carrier.
Preservative:	0.05% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
Purification:	Purified from rabbit serum by sequential epitope-specific chromatography. The antibody has been negatively pre-adsorbed using a non-phosphopeptide corresponding to the site of phosphorylation to remove antibody that is reactive with non-phosphorylated Tie2. The final product is generated by affinity chromatography using a Tie2-derived peptide that is phosphorylated at tyrosine 1108.
Immunogen:	The antiserum was produced against a chemically synthesized phosphopeptide derived from the region of human Tie2 that contains tyrosine 1108.
Target Summary:	Tie2 (also known as TEK) is a 140 kDa endothelial receptor tyrosine kinase that regulates signaling events mediating embryonic vascular development and tumor angiogenesis. Disruption of Tie2 receptor signaling results in embryonic lethality due to vascular hemorrhage and impaired cardiac development. Tie2 receptor consists of an extracellular region that contains three EGF-like modules flanked by two immunoglobulin (Ig)-like domains, three fibronectin type III (FNIII) repeats, and a cytoplasmic region that contains tyrosine kinase domains including several phosphorylation and binding sites. Phosphorylation of Tie2 at distinct tyrosine residues plays a critical role in vascular development and hematopoiesis including tyrosine 992 in the activation loop which primes the phosphorylation of Tie2 on other tyrosine residues including tyrosine 1102 (also known as tyrosine 1100) and 1108 (also referred to as tyrosine 1106). Tyrosine 1108 is an angiopoietin (Ang 1)-dependent autophosphorylation site that mediates binding and phosphorylation of Dok-R and plays an important role in linking Ang-1 stimulation with downstream cell migration signaling in endothelial cells.
Reactivity:	Human Tie2. Mouse and rat Tie2 (100%) have not been tested but are expected to react.
Applications:	The antibody has been used for Western blotting applications.
Suggested Working Dilutions:	For Western blotting applications, we recommend using the antibody at a 1:1000 dilution. The optimal antibody concentration should be determined empirically for each specific application.
Storage:	Store at -20°C . We recommend a brief centrifugation before opening to settle vial contents. Then, apportion into working aliquots and store at -20°C . For shipment or short-term storage (up to one week), $2-8^{\circ}\text{C}$ is sufficient.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
Positive Controls Used:	Human umbilical vein endothelial (HUVAC) cells plated on fibronectin.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com
This antibody is manufactured under a licensed process covered by Patent # 5, 599, 681.

PI441313G

(Rev 11/08) 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.

Related Products:**Antibodies:**

Tie2 [pY⁹⁹²], Cat. # 44-1310G
 VEGFR2 [pY⁹⁴⁹], Cat # 44- 1041G
 VEGFR2 [pY⁹⁵¹], Cat. # 44-1040G
 VEGFR2 [pY¹⁰⁵⁴], Cat # 44-1046G
 VEGFR2 [pY^{1054/1059}], Cat. # 44-1047G
 FAK [pY³⁹⁷], Cat. # 44-624G
 FAK [pY⁵⁷⁶], Cat. # 44-652G

FAK [pS⁸⁴³], Cat. # 44-594
 Integrin $\alpha 4$ [pS⁹⁸⁸], Cat. # 44-864
 Integrin $\beta 1$ [pS⁷⁸⁵], Cat. # 44-870
 Integrin $\beta 1$ [pTpT^{788/789}], Cat. # 44-872
 Integrin $\beta 3$ [pY⁷⁷³], Cat. # 44-876
 Integrin $\beta 3$ [pY⁷⁸⁵], Cat. # 44-878
 VE-cadherin [pY⁷³¹], Cat. # 44-1145G

References:

Harfouche, R. and S.N. Hussian (2006) Signaling and regulation of endothelial cell survival by angiopoietin-2. *Am. J. Physiol. Heart Circ. Physiol.* 291(4):H1635-H1645.

Tsigkos, S., et al. (2006) Regulation of Ang2 release by PTEN/PI3-kinase/Akt in lung microvascular endothelial cells. *J. Cell Physiol.* 207(2):506-511.

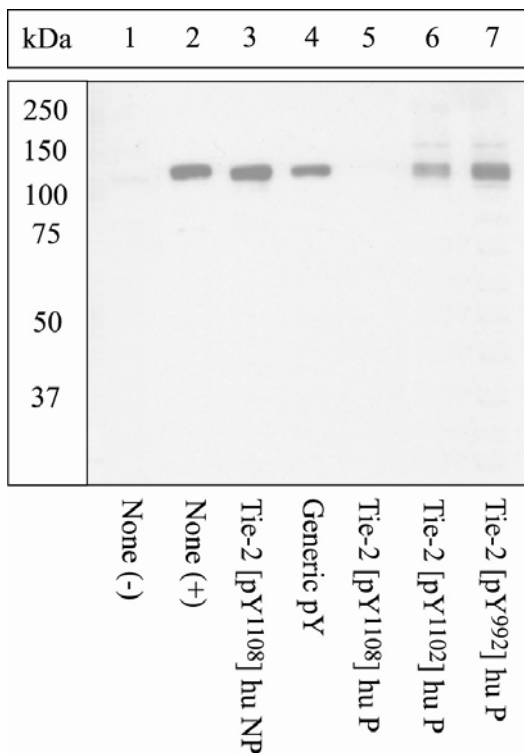
Cascone, I., et al. (2005) Stable interaction between alpha5beta1 integrin and Tie2 tyrosine kinase receptor regulates endothelial cell response to Ang-1. *J. Cell Biol.* 170(6):993-1004.

Valable, S., et al. (2003) Angiopoietin-1-induced PI3-kinase activation prevents neuronal apoptosis. *FASEB J.* 17(3):443-445.

Niu, X.L., et al. (2002) Deletion of the carboxyl terminus of Tie2 enhances kinase activity, signaling, and function. Evidence for an autoinhibitory mechanism. *J. Biol. Chem.* 277(35):31768-31773.

Murray, B.W., et al. (2001) Mechanistic effects of autophosphorylation on receptor tyrosine kinase catalysis: enzymatic characterization of Tie2 and phospho-Tie2. *Biochemistry* 40(34):10243-10253.

Teichert-Kuliszewska, K., et al. (2001) Biological action of angiopoietin-2 in a fibrin matrix model of angiogenesis is associated with activation of Tie2. *Cardiovasc. Res.* 49(3):659-670.

**Antibody Specificity**

Lysates prepared from HUVAC cells left alone (1) or allowed to grow on fibronectin-coated plates for 90 min (2-7) were resolved on a 10% polyacrylamide gel and transferred to PVDF. The membrane was blocked with a 3% BSA-TBST buffer for one hour at room temperature, and then incubated with the Tie-2 [pY¹¹⁰⁸] antibody for two hours at room temperature in 3% BSA-TBST buffer, following prior incubation with: no peptide (2), the non-phosphopeptide corresponding to the phosphopeptide immunogen (3) a generic phosphotyrosine-containing peptide (4) or the phosphopeptide immunogen corresponding to Tie2 [pY¹¹⁰⁸] (5), a phosphopeptide immunogen corresponding to Tie2 [pY¹¹⁰²] (6) or phosphopeptide corresponding to Tie2 [pY⁹⁹²] (7). After washing, the membrane was incubated with goat F(ab')₂ anti-rabbit IgG HRP conjugate (Cat. # ALI4404) and signals were detected using the Pierce SuperSignal™ reagent.

The data show that the signal was induced by fibronectin and selectively blocked by the phosphopeptide corresponding to Tie2 [pY¹¹⁰⁸] indicating that the signal is phosphorylation site-specific.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

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

This antibody is manufactured under a licensed process covered by Patent # 5, 599, 681.

PI441313G

(Rev 11/08) 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.

Western Blotting Procedure

1. Lyse approximately 10^7 cells in 0.5 mL of ice cold Cell Lysis Buffer (formulation provided below). This buffer, a modified RIPA buffer, is suitable for recovery of most proteins, including membrane receptors, cytoskeletal-associated proteins, and soluble proteins. This cell lysis buffer formulation is available as a separate product which requires supplementation with protease inhibitors immediately prior to use (Cat. # FNN0011). Other cell lysis buffer formulations, such as Laemmli sample buffer and Triton-X 100 buffer, are also compatible with this procedure. Additional optimization of the cell stimulation protocol and cell lysis procedure may be required for each specific application.
2. Remove the cellular debris by centrifuging the lysates at 14,000 x g for 10 minutes. Alternatively, lysates may be ultracentrifuged at 100,000 x g for 30 minutes for greater clarification.
3. Carefully decant the clarified cell lysates into clean tubes and determine the protein concentration using a suitable method, such as the Bradford assay. Polypropylene tubes are recommended for storing cell lysates.
4. React an aliquot of the lysate with an equal volume of 2x Laemmli Sample Buffer (125 mM Tris, pH 6.8, 10% glycerol, 10% SDS, 0.006% bromophenol blue, and 130 mM dithiothreitol [DTT]) and boil the mixture for 90 seconds at 100°C.
5. Load 10-30 µg of the cell lysate into the wells of an appropriate single percentage or gradient minigel and resolve the proteins by SDS-PAGE.
6. In preparation for the Western transfer, cut a piece of PVDF membrane slightly larger than the gel. Soak the membrane in methanol for 1 minute, then rinse with ddH₂O for 5 minutes. Alternatively, nitrocellulose may be used.
7. Soak the membrane, 2 pieces of Whatman paper, and Western apparatus sponges in transfer buffer (formulation provided below) for 2 minutes.
8. Assemble the gel and membrane into the sandwich apparatus.
9. Transfer the proteins at 140 mA for 60-90 minutes at room temperature.
10. Following the transfer, rinse the membrane with Tris buffered saline for 2 minutes.
11. Block the membrane with blocking buffer (formulation provided below) for one hour at room temperature or overnight at 4°C.
12. Incubate the blocked blot with primary antibody at a 1:1000 dilution in Tris buffered saline supplemented with 3% Ig-free BSA and 0.1% Tween 20 overnight at 4°C or for one hour at room temperature.
13. Wash the blot with several changes of Tris buffered saline supplemented with 0.1% Tween 20.
14. Detect the antibody band using an appropriate secondary antibody, such as goat F(ab')₂ anti-rabbit IgG alkaline phosphatase conjugate (Cat. # ALI4405) or goat F(ab')₂ anti-rabbit IgG horseradish peroxidase conjugate (Cat. # ALI4404) in conjunction with your chemiluminescence reagents and instrumentation.

Cell Lysis Buffer

Formulation:

10 mM Tris, pH 7.4
100 mM NaCl
1 mM EDTA
1 mM EGTA
1 mM NaF
20 mM Na₄P₂O₇
2 mM Na₃VO₄
0.1% SDS
0.5% sodium deoxycholate
1% Triton-X 100
10% glycerol
1 mM PMSF (made from a 0.3 M stock in DMSO) or 1 mM AEBSF (water soluble version of PMSF)
60 µg/mL aprotinin
10 µg/mL leupeptin
1 µg/mL pepstatin
(alternatively, protease inhibitor cocktail such as Sigma Cat. # P2714 may be used)

Transfer Buffer

Formulation:

2.4 gm Tris base
14.2 gm glycine
200 mL methanol
Q.S. to 1 liter, then add
1 mL 10% SDS.
Cool to 4°C prior to use.

Tris Buffered Saline

Formulation:

20 mM Tris-HCl, pH 7.4
0.9% NaCl

Blocking Buffer

Formulation:

100 mL Tris buffered saline
3 gm Ig-free BSA
0.1 mL Tween 20

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

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

This antibody is manufactured under a licensed process covered by Patent # 5, 599, 681.

PI441313G

(Rev 11/08) 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.