

# Quantikine<sup>®</sup>

## Human Tie-1 Immunoassay

Catalog Number DTE100

**For the quantitative determination of human Tie-1 concentrations in cell culture supernates, cell lysates, serum, and plasma.**

*This package insert must be read in its entirety before using this product.*

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

## TABLE OF CONTENTS

Contents	Page
INTRODUCTION	2
PRINCIPLE OF THE ASSAY . . . . .	3
LIMITATIONS OF THE PROCEDURE	3
MATERIALS PROVIDED . . . . .	4
STORAGE	4
OTHER SUPPLIES REQUIRED . . . . .	5
PRECAUTION	5
SAMPLE COLLECTION AND STORAGE . . . . .	5
CELL LYSIS PROCEDURE	5
SAMPLE PREPARATION . . . . .	6
REAGENT PREPARATION	6
ASSAY PROCEDURE . . . . .	7
ASSAY PROCEDURE SUMMARY	8
CALCULATION OF RESULTS. . . . .	9
TYPICAL DATA	9
TECHNICAL HINTS . . . . .	10
PRECISION	10
RECOVERY . . . . .	11
LINEARITY	11
SENSITIVITY . . . . .	11
CALIBRATION	11
SAMPLE VALUES . . . . .	12
SPECIFICITY	13
REFERENCES . . . . .	14
PLATE LAYOUT	15

### MANUFACTURED AND DISTRIBUTED BY:

R&D Systems, Inc.	TELEPHONE:	(800) 343-7475
614 McKinley Place NE		(612) 379-2956
Minneapolis, MN 55413	FAX:	(612) 656-4400
United States of America	E-MAIL:	info@RnDSystems.com

### DISTRIBUTED BY:

R&D Systems Europe, Ltd.	TELEPHONE:	+44 (0)1235 529449
19 Barton Lane		
Abingdon Science Park	FAX:	+44 (0)1235 533420
Abingdon, OX14 3NB	E-MAIL:	info@RnDSystems.co.uk
United Kingdom		

R&D Systems GmbH	TELEPHONE:	+49 (0)6122 90980
Borsigstrasse 7		
65205 Wiesbaden-Nordenstadt	FAX:	+49 (0)6122 909819
Germany	E-MAIL:	infogmbh@RnDSystems.co.uk

R&D Systems Europe	FREEPHONE:	+0800 90 72 49
77 boulevard Vauban		
59041 LILLE CEDEX	FAX:	+0800 77 16 68
France	E-MAIL:	info@RnDSystems.co.uk

## INTRODUCTION

Tie-1 is a 130 kDa type I transmembrane protein and member of the protein kinase superfamily, the tyrosine protein kinase family, and the Tie subfamily of receptor tyrosine kinases (1). Human Tie-1 is synthesized as an 1138 amino acid (aa) precursor with a 21 aa signal sequence, a 759 aa extracellular domain (ECD), a 25 aa transmembrane sequence, and a 354 aa cytoplasmic domain (SwissProt #: P35590). The ECD contains two immunoglobulin-like (Ig-like C2-type) domains, three EGF-like domains, and three fibronectin type III domains (2). In addition, there are five sites for potential N-linked glycosylation and one  $\text{Ca}^{2+}$  binding site (2). The cytoplasmic domain contains one tyrosine kinase domain (2). Mature human Tie-1 is 93% aa identical to mature mouse Tie-1. The expression of the Tie-1 gene is restricted to endothelial cells and some hematopoietic cell lineages (1, 3).

Targeted Tie-1 disruption results in embryonic lethality due to edema, hemorrhage, and microvascular rupture (4 - 7). Tie-1 deficient embryos also contain hyperactive endothelial cells that exhibit a large number of extensions and filopodia projecting into vessel lumen and have increased capillary densities, suggesting that Tie-1 may act as a brake during embryonic angiogenesis (4, 7). Upregulation of Tie-1 has been observed during wound healing, ovarian follicle maturation, and tumor angiogenesis (3, 8, 9). Tie-1 is expressed in gastrointestinal stromal tumors, leiomyoma, and schwannoma (10), is an independent prognostic marker for gastric cancer (11), and has also been detected in epithelial tumors in breast and thyroid (12). Recently, an alternative splice variant in Tie-1 has been shown to reduce rheumatoid arthritis severity (13).

Tie-1 shares a high degree of homology with Tie-2, the receptor for three members of the angiopoietin family (Ang-1, Ang-2, and Ang-3/4) (3, 14 - 17). A soluble Ang-1 chimeric protein, COMP-Ang-1, and native Ang-1 and Ang-4 also activate Tie-1 (3). They all induce Tie-1 phosphorylation, although native Ang-1 and Ang-4 are less efficient than COMP-Ang-1 (3). It is reported that Ang-1 induces Tie-1 phosphorylation; however, it is Tie-2 dependent (4). The research shows that Ang-1 fails to induce Tie-1 phosphorylation when Tie-2 is down-regulated in endothelial cells. Tie-1 phosphorylation is induced in the absence of Ang-1 by either a constitutively active form of Tie-2 or a Tie-2 agonistic antibody. In HEK293 cells, Ang-1 phosphorylates a form of Tie-1 without kinase activity when co-expressed with kinase-defective Tie-2 (4). Research also shows that a main role for Tie-1 is to modulate blood vessel morphogenesis by virtue of its ability to down-regulate Tie-2-driven signaling and endothelial survival (4). Both Tie-1 and Tie-2 are also needed for postnatal bone marrow hematopoiesis (3, 18).

The Quantikine Human Tie-1 immunoassay is a 4.5 hour solid phase ELISA designed to measure human Tie-1 in cell culture supernates, cell lysates, serum, and plasma. It contains NS0-expressed recombinant human Tie-1 and has been shown to accurately quantitate the recombinant factor. Results obtained using natural human Tie-1 showed dose-response curves that were parallel to the standard curves obtained using the Quantikine kit standards. These results indicate that the Quantikine Human Tie-1 kit can be used to determine relative mass values for naturally occurring human Tie-1.

## **PRINCIPLE OF THE ASSAY**

This assay employs the quantitative sandwich enzyme immunoassay technique. A monoclonal antibody specific for Tie-1 has been pre-coated onto a microplate. Standards and samples are pipetted into the wells and any Tie-1 present is bound by the immobilized antibody. After washing away any unbound substances, an enzyme-linked polyclonal antibody specific for Tie-1 is added to the wells. Following a wash to remove any unbound antibody-enzyme reagent, a substrate solution is added to the wells and color develops in proportion to the amount of Tie-1 bound in the initial step. The color development is stopped and the intensity of the color is measured.

## **LIMITATIONS OF THE PROCEDURE**

- FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.
- The kit should not be used beyond the expiration date on the kit label.
- Do not mix or substitute reagents with those from other lots or sources.
- If samples generate values higher than the highest standard, dilute samples with the appropriate Calibrator Diluent and repeat the assay.
- Any variation in standard diluent, operator, pipetting technique, washing technique, incubation time or temperature, and kit age can cause variation in binding.
- This assay is designed to eliminate interference by ligands, binding proteins, and other factors present in biological samples. Until all factors have been tested in the Quantikine Immunoassay, the possibility of interference cannot be excluded.

## MATERIALS PROVIDED

**Tie-1 Microplate** (Part 893245) - 96 well polystyrene microplate (12 strips of 8 wells) coated with a mouse monoclonal antibody against Tie-1.

**Tie-1 Conjugate** (Part 893246) - 21 mL of polyclonal antibody against Tie-1 conjugated to horseradish peroxidase with preservatives.

**Tie-1 Standard** (Part 893247) - 100 ng of recombinant human Tie-1 in a buffered protein solution with preservatives; lyophilized.

**Assay Diluent RD1-89** (Part 895881) - 11 mL of a buffered protein solution with preservatives.

**Calibrator Diluent RD5-17** (Part 895512) - 21 mL of a buffered protein solution with preservatives.

**Wash Buffer Concentrate** (Part 895003) - 21 mL of a 25-fold concentrated solution of buffered surfactant with preservatives.

**Color Reagent A** (Part 895000) - 12.5 mL of stabilized hydrogen peroxide.

**Color Reagent B** (Part 895001) - 12.5 mL of stabilized chromogen (tetramethylbenzidine).

**Stop Solution** (Part 895032) - 6 mL of 2 N sulfuric acid.

**Plate Covers** - 4 adhesive strips.

## STORAGE

Unopened Kit	Store at 2 - 8° C. Do not use past kit expiration date.	
Opened/ Reconstituted Reagents	Diluted Wash Buffer	May be stored for up to 1 month at 2 - 8° C.*
	Stop Solution	
	Assay Diluent RD1-89	
	Calibrator Diluent RD5-17	
	Conjugate	
	Unmixed Color Reagent A	
	Unmixed Color Reagent B	
	Standard	
	Microplate Wells	Return unused wells to the foil pouch containing the desiccant pack, reseal along entire edge of zip-seal. May be stored for up to 1 month at 2 - 8° C.*

\*Provided this is within the expiration date of the kit.

## OTHER SUPPLIES REQUIRED

- Microplate reader capable of measuring absorbance at 450 nm, with the correction wavelength set at 540 nm or 570 nm.
- Pipettes and pipette tips.
- Deionized or distilled water.
- Squirt bottle, manifold dispenser, or automated microplate washer.
- 500 mL graduated cylinder.
- Human Tie-1 Controls (optional; available from R&D Systems).

### **If using cell lysate samples, the following supplies are also required:**

- Cell Lysis Buffer 3 (R&D Systems, Catalog # 895366) or equivalent.
- Phosphate-buffered saline (PBS).

## PRECAUTION

The Stop Solution provided with this kit is an acid solution. Wear eye, hand, face, and clothing protection when using this material.

## SAMPLE COLLECTION AND STORAGE

**Cell Culture Supernates** - Remove particulates by centrifugation and assay immediately or aliquot and store samples at  $\leq -20^{\circ}$  C. Avoid repeated freeze-thaw cycles.

**Cell Lysates** - Cells must be lysed before assaying. Refer to the Cell Lysis Procedure below.

**Serum** - Use a serum separator tube (SST) and allow samples to clot for 30 minutes before centrifugation for 15 minutes at 1000 x g. Remove serum and assay immediately or aliquot and store samples at  $\leq -20^{\circ}$  C. Avoid repeated freeze-thaw cycles.

**Plasma** - Collect plasma using heparin or EDTA as an anticoagulant. Centrifuge for 15 minutes at 1000 x g within 30 minutes of collection. Assay immediately or aliquot and store samples at  $\leq -20^{\circ}$  C. Avoid repeated freeze-thaw cycles.

**Note:** *Citrate plasma has not been validated for use in this assay.*

## CELL LYSIS PROCEDURE

Use the following procedure for the preparation of cell lysate samples.

1. Perform a 5-fold dilution of Cell Lysis Buffer 3 (available from R&D Systems) with deionized or distilled water.
2. Wash cells one time in cold PBS.
3. Resuspend cells at  $1 \times 10^6$  cells/mL in diluted lysis buffer.
4. Incubate with gentle agitation for 30 minutes at  $37^{\circ}$  C.
5. Centrifuge to remove cell debris.
6. Assay immediately or aliquot the lysis supernates and store at  $\leq -20^{\circ}$  C until ready for use.

## SAMPLE PREPARATION

Serum and plasma samples require a 10-fold dilution. A suggested 10-fold dilution is 20  $\mu$ L of sample + 180  $\mu$ L of Calibrator Diluent RD5-17.

## REAGENT PREPARATION

**Bring all reagents to room temperature before use.**

**Wash Buffer** - If crystals have formed in the concentrate, warm to room temperature and mix gently until the crystals have completely dissolved. Dilute 20 mL of Wash Buffer Concentrate into deionized or distilled water to prepare 500 mL of Wash Buffer.

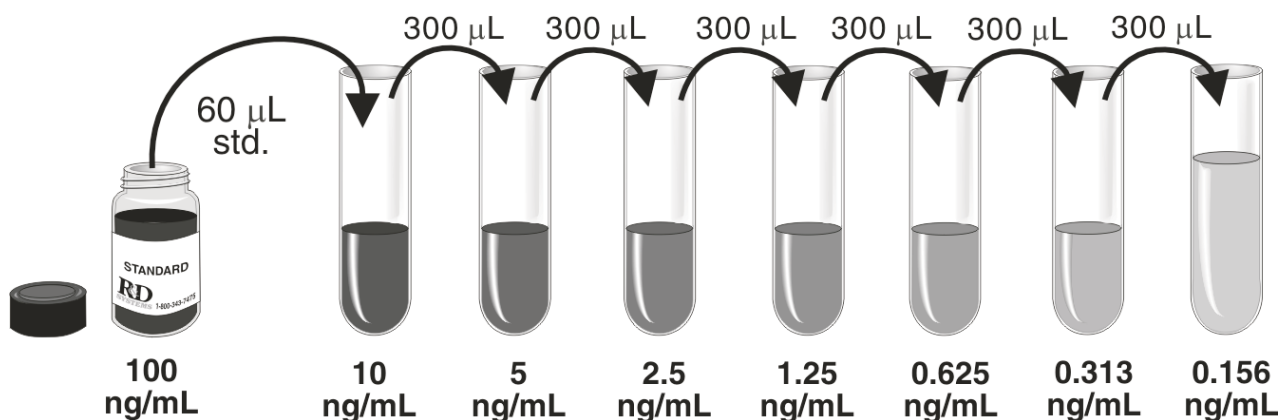
**Substrate Solution** - Color Reagents A and B should be mixed together in equal volumes within 15 minutes of use. Protect from light. 200  $\mu$ L of the resultant mixture is required per well.

**Calibrator Diluent RD5-17 (1:2)** - A 2-fold dilution is required for use with cell culture supernate samples and cell lysates. A suggested 2-fold dilution is 5.0 mL of Calibrator Diluent RD5-17 + 5.0 mL of deionized or distilled water.

**Note:** *Do not dilute Calibrator Diluent RD5-17 for use with serum or plasma samples.*

**Tie-1 Standard** - Reconstitute the Tie-1 Standard with 1.0 mL of deionized or distilled water. This reconstitution produces a stock solution of 100 ng/mL. Mix the standard to ensure complete reconstitution and allow the standard to sit for a minimum of 15 minutes with gentle agitation prior to making dilutions.

Pipette 540  $\mu$ L of Calibrator Diluent RD5-17 (*for serum and plasma samples*) or Calibrator Diluent RD5-17 (1:2) (*for cell culture supernate/cell lysate samples*) into the 10 ng/mL tube. Pipette 300  $\mu$ L of the appropriate Calibrator Diluent RD5-17 into the remaining tubes. Use the stock solution to produce a dilution series (below). Mix each tube thoroughly before the next transfer. The 10 ng/mL standard serves as the high standard. The appropriate Calibrator Diluent serves as the zero standard (0 ng/mL).



## ASSAY PROCEDURE

**Bring all reagents and samples to room temperature before use. It is recommended that all samples, controls, and standards be assayed in duplicate.**

1. Prepare all reagents, working standards, and samples as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, and reseal.
3. Add 100  $\mu$ L of Assay Diluent RD1-89 to each well.
4. Add 50  $\mu$ L of Standard, control, or sample\* per well. Cover with the adhesive strip provided. Incubate for 2 hours at room temperature. A plate layout is provided to record standards and samples assayed.
5. Aspirate each well and wash, repeating the process three times for a total of four washes. Wash by filling each well with Wash Buffer (400  $\mu$ L) using a squirt bottle, manifold dispenser, or autowasher. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
6. Add 200  $\mu$ L of Tie-1 Conjugate to each well. Cover with a new adhesive strip. Incubate for 2 hours at room temperature.
7. Repeat the aspiration/wash as in step 5.
8. Add 200  $\mu$ L of Substrate Solution to each well. Incubate for 30 minutes at room temperature. **Protect from light.**
9. Add 50  $\mu$ L of Stop Solution to each well. The color in the wells should change from blue to yellow. If the color in the wells is green or the color change does not appear uniform, gently tap the plate to ensure thorough mixing.
10. Determine the optical density of each well within 30 minutes, using a microplate reader set to 450 nm. If wavelength correction is available, set to 540 nm or 570 nm. If wavelength correction is not available, subtract readings at 540 nm or 570 nm from the readings at 450 nm. This subtraction will correct for optical imperfections in the plate. Readings made directly at 450 nm without correction may be higher and less accurate.

\*Serum/Plasma samples require dilution. See the Sample Preparation section.



## ASSAY PROCEDURE SUMMARY

1. Prepare reagents, samples, and standards as instructed.



2. Add 100  $\mu\text{L}$  Assay Diluent RD1-89 to each well.



3. Add 50  $\mu\text{L}$  Standard, control, or sample\* to each well.  
Incubate 2 hours at RT.



4. Aspirate and wash 4 times.



5. Add 200  $\mu\text{L}$  Conjugate to each well.  
Incubate 2 hours at RT.



6. Aspirate and wash 4 times.



7. Add 200  $\mu\text{L}$  Substrate Solution to each well.  
Incubate 30 minutes. **Protect from light.**



8. Add 50  $\mu\text{L}$  Stop Solution to each well.  
Read at 450 nm within 30 min.  
 $\lambda$  correction 540 or 570 nm

\*Serum/Plasma samples require dilution.  
See the Sample Preparation section.

## CALCULATION OF RESULTS

Average the duplicate readings for each standard, control, and sample and subtract the average zero standard optical density.

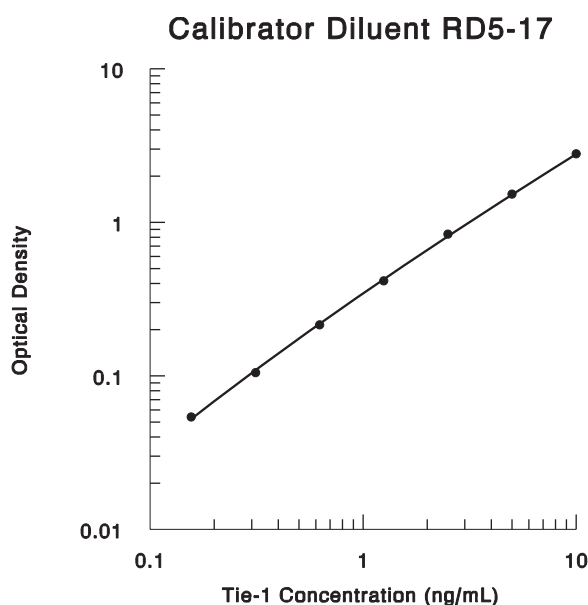
Plot the optical density for the standards versus the concentration of the standards and draw the best fit curve. The data can be linearized by using log/log paper and regression analysis may be applied to the log transformation.

To determine the Tie-1 concentration of each sample, first find the absorbance value on the y-axis and extend a horizontal line to the standard curve. At the point of intersection, extend a vertical line to the x-axis and read the corresponding Tie-1 concentration.

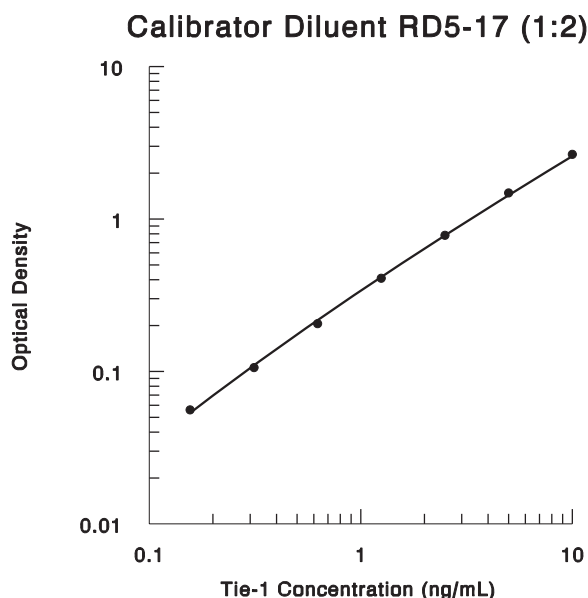
If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

## TYPICAL DATA

This standard curve is provided for demonstration only. A standard curve should be generated for each set of samples assayed.



ng/mL	O.D.	Average	Corrected
0	0.010 0.013 0.064	0.012	—
0.156	0.067 0.116	0.066	0.054
0.313	0.118 0.225	0.117	0.105
0.625	0.228 0.421	0.227	0.215
1.25	0.434 0.826	0.428	0.416
2.5	0.874 1.536	0.850	0.838
5	1.545 2.763	1.541	1.529
10	2.854	2.809	2.797



ng/mL	O.D.	Average	Corrected
0	0.010 0.011 0.054	0.011	—
0.156	0.057 0.101	0.056	0.045
0.313	0.111 0.199	0.106	0.095
0.625	0.212 0.376	0.206	0.195
1.25	0.442 0.773	0.409	0.398
2.5	0.790 1.460	0.782	0.771
5	1.505 2.596	1.483	1.472
10	2.719	2.658	2.647

## TECHNICAL HINTS

- When mixing or reconstituting protein solutions, always avoid foaming.
- To avoid cross-contamination, change pipette tips between additions of each standard level, between sample additions, and between reagent additions. Also, use separate reservoirs for each reagent.
- When using an automated plate washer, adding a 30 second soak period following the addition of wash buffer, and/or rotating the plate 180 degrees between wash steps may improve assay precision.
- To ensure accurate results, proper adhesion of plate sealers during incubation steps is necessary.
- Substrate Solution should remain colorless until added to the plate. Keep Substrate Solution protected from light. Substrate Solution should change from colorless to gradations of blue.
- Stop Solution should be added to the plate in the same order as the Substrate Solution. The color developed in the wells will turn from blue to yellow upon addition of the Stop Solution. Wells that are green in color indicate that the Stop Solution has not mixed thoroughly with the Substrate Solution.

## PRECISION

### Intra-assay Precision (Precision within an assay)

Three samples of known concentration were tested twenty times on one plate to assess intra-assay precision.

### Inter-assay Precision (Precision between assays)

Three samples of known concentration were tested in forty separate assays to assess inter-assay precision.

#### Serum/Plasma Assay

Sample	Intra-assay Precision			Inter-assay Precision		
	1	2	3	1	2	3
n	20	20	20	40	40	40
Mean (ng/mL)	1.17	2.23	4.84	1.21	2.20	4.84
Standard deviation	0.06	0.07	0.16	0.08	0.14	0.17
CV (%)	5.1	3.1	3.3	6.6	6.4	3.5

#### Cell Culture Supernate/Cell Lysate Assay

Sample	Intra-assay Precision			Inter-assay Precision		
	1	2	3	1	2	3
n	20	20	20	40	40	40
Mean (ng/mL)	1.41	2.55	5.73	1.34	2.42	5.27
Standard deviation	0.08	0.09	0.26	0.10	0.15	0.23
CV (%)	5.7	3.5	4.5	7.5	6.2	4.4

## RECOVERY

The recovery of Tie-1 spiked to levels throughout the range of the assay was evaluated.

Sample	Average % Recovery	Range
Cell culture media (n=4)	99	90 - 108%

## LINEARITY

To assess the linearity of the assay, samples containing and/or spiked with high concentrations of Tie-1 were serially diluted with the appropriate Calibrator Diluent to produce samples with values within the dynamic range of the assay.

		Cell culture media (n=4)	Cell lysates (n=3)	Serum* (n=4)	Heparin plasma* (n=4)	EDTA plasma* (n=4)
1:2	Average % of Expected	104	95	101	106	102
	Range (%)	99 - 110	85 - 103	98 - 105	100 - 109	99 - 104
1:4	Average % of Expected	105	92	102	106	102
	Range (%)	94 - 114	85 - 97	99 - 105	98 - 111	97 - 106
1:8	Average % of Expected	102	98	101	103	99
	Range (%)	92 - 109	—	95 - 105	97 - 109	93 - 104
1:16	Average % of Expected	98	97	101	102	96
	Range (%)	87 - 108	—	91 - 110	89 - 111	90 - 100

\*Samples were diluted prior to assay, as directed in the Sample Preparation section.

## SENSITIVITY

One hundred ten assays were evaluated and the minimum detectable dose (MDD) of Tie-1 ranged from 0.003 - 0.021 ng/mL. The mean MDD was 0.010 ng/mL.

The MDD was determined by adding two standard deviations to the mean optical density value of twenty zero standard replicates and calculating the corresponding concentration.

## CALIBRATION

This immunoassay is calibrated against a highly purified NS0-expressed recombinant human Tie-1 produced at R&D Systems.

## SAMPLE VALUES

**Serum/Plasma** - Samples drawn from apparently healthy volunteers were evaluated for the presence of Tie-1 in this assay. No medical histories were available for the donors used in this study.

Sample Type	Mean (ng/mL)	Range (ng/mL)	Standard Deviation (ng/mL)
Serum* (n=35)	61.2	30.2 - 88.0	12.1
EDTA plasma* (n=35)	58.8	32.1 - 92.3	11.0
Heparin plasma* (n=35)	59.4	28.4 - 96.3	11.9

\*Samples were diluted prior to assay, as directed in the Sample Preparation section.

### Cell Culture Supernates -

Human peripheral blood lymphocytes ( $1 \times 10^6$  cells/mL) were cultured in DMEM supplemented with 5% fetal calf serum, 5  $\mu$ M  $\beta$ -mercaptoethanol, 2 mM L-glutamine, 100 U/mL penicillin, and 100  $\mu$ g/mL streptomycin sulfate. Cells were cultured unstimulated or stimulated with 10  $\mu$ g/mL PHA for 1 and 6 days. Aliquots of the cell culture supernates were removed and assayed for levels of natural Tie-1.

Condition	Day 1 (ng/mL)	Day 6 (ng/mL)
Unstimulated	ND	0.340
Stimulated	ND	0.226

ND = Non-detectable

Human umbilical vein endothelial cells (HUVEC) were grown in EGM-2 media at 37° C, 95% humidity, and 5% CO<sub>2</sub> for 3 days. Cells were cultured unstimulated or stimulated with 40 ng/mL of recombinant human TNF- $\alpha$  and 50 ng/mL of recombinant human IFN- $\gamma$ . Aliquots of the cell culture supernates were removed and assayed for levels of natural Tie-1.

Condition	Day 3 (ng/mL)
Unstimulated	6.16
Stimulated	7.92

Human lung microvascular endothelial cells (HMVEC) were grown in EGM-2 media at 37° C, 95% humidity, and 5% CO<sub>2</sub>. An aliquot of the cell culture supernate was removed, assayed for levels of natural Tie-1, and measured 1.190 ng/mL.

## Cell Lysates -

Human umbilical vein endothelial cells (HUVEC) were grown in EGM-2 media at 37° C, 95% humidity, and 5% CO<sub>2</sub> for 3 days. Cells were cultured unstimulated or stimulated with 40 ng/mL of recombinant human TNF- $\alpha$  and 50 ng/mL of recombinant human IFN- $\gamma$  for 1 day. The cell culture supernate was removed, and cells were lysed as described in the Cell Lysis Procedure. The cell lysates were assayed for levels of natural Tie-1.

Condition	Day 1 (ng/mL)
Unstimulated	1.13
Stimulated	0.186

Human umbilical vein endothelial cells (HUVEC) were grown in EGM-2 media at 37° C, 95% humidity, and 5% CO<sub>2</sub> for 5 days. Cells were then stimulated with 10 ng/mL of recombinant human IL-1 $\beta$  for 1 day. The cell culture supernate was removed, and cells were lysed as described in the Cell Lysis Procedure. The cell lysate was assayed for levels of natural Tie-1 and measured 0.886 ng/mL.

Human lung microvascular endothelial cells (HMVEC) were grown in EGM-2 media at 37° C, 95% humidity, and 5% CO<sub>2</sub>. The cell culture supernate was removed, and cells were lysed as described in the Cell Lysis Procedure. The cell lysate was assayed for levels of natural Tie-1 and measured 2.60 ng/mL.

## SPECIFICITY

This assay recognizes recombinant and natural human Tie-1. The factors listed below were prepared at 100 ng/mL in the appropriate Calibrator Diluent and assayed for cross-reactivity. Preparations of the following factors at 100 ng/mL in a mid-range recombinant human Tie-1 control were assayed for interference. No significant cross-reactivity or interference was observed.

### Recombinant human:

Ang-1  
Ang-2  
Ang-4  
ANGPT-L1  
ANGPT-L7  
EGF  
EG-VEGF  
HB-EGF  
Tie-2  
VEGF<sub>121</sub>  
VEGF<sub>162</sub>  
VEGF<sub>165</sub>  
VEGF-C  
VEGF/P/IGF

### Recombinant mouse:

Ang-3  
EGF  
EG-VEGF  
Tie-1  
Tie-2  
VEGF<sub>120</sub>  
VEGF<sub>164</sub>  
VEGF-B<sub>167</sub>  
VEGF-B<sub>186</sub>

### Recombinant rat:

EGF  
VEGF<sub>164</sub>

### Recombinant zebrafish:

Tie-2  
VEGF<sub>165</sub>

## REFERENCES

1. Hashiyama, M. *et al.* (1996) Blood **87**:93.
2. Partanen, J. *et al.* (1992) Mol. Cell. Biol. **12**:1698.
3. Saharinen, P. *et al.* (2005) J. Cell Biol. **169**:239.
4. Yuan, H.T. *et al.* (2007) FASEB J. **21**:3171.
5. Puri, M.C. *et al.* (1995) EMBO J. **14**:5884.
6. Sato, T.N. *et al.* (1995) Nature **376**:70
7. Patan, S. *et al.* (1998) Microvasc. Res. **56**:1.
8. Korhonen, J. *et al.* (1992) Blood **80**:2548.
9. Kaipainen, A. *et al.* (1994) Cancer Res. **54**:6571.
10. Nakayama, T. *et al.* (2007) World J. Gastroenterol. **13**:4473.
11. Lin, W-C. *et al.* (1999) Clin. Cancer Res. **5**:1745.
12. Rees, K.A. *et al.* (2007) Int. J. Oncol. **31**:893.
13. Jin, P. *et al.* (2008) Arthritis Res. Ther. **10**:R73.
14. Peters, K.G. *et al.* (2004) Recent Prog. Hormone. Res. **59**:51.
15. Yancopoulos, G.D. *et al.* (2000) Nature **407**:242.
16. Jones, N. *et al.* (2001) Nat. Rev. Mol. Cell. Biol. **2**:257.
17. Lee, H.J. *et al.* (2004) FASEB J. **18**:1200.
18. Puri, M.C. and A. Bernstein (2003) Proc. Natl. Acad. Sci. USA **100**:12753.

**PLATE LAYOUT**

Use this plate layout as a record of standards and samples assayed.

1								
2								
3								
4								
5								
6								
7								
8								
9								
10								
11								
12								
	A	B	C	D	E	F	G	H