



ELISA Kit
Catalog #KHO0281

Human
c-Met
**[pYpYpY1230/
1234/1235]**

www.invitrogen.com
Invitrogen Corporation
542 Flynn Road, Camarillo, CA 93012
Tel: 800-955-6288
E-mail: techsupport@invitrogen.com

TABLE OF CONTENTS

Introduction	4
Principle of the Method.....	6
Reagents Provided.....	7
Supplies Required but Not Provided	8
Procedural Notes/Lab Quality Control	8
Safety	10
Directions for Washing.....	10
Procedure for Extraction of Proteins from Cells.....	11
Reagent Preparation and Storage.....	13
Reconstitution and Dilution of Hu c-Met [pYpYpY1230/1234/1235] Standard.....	13
Dilution of Hu c-Met [pYpYpY1230/1234/1235] Standard	14
Storage and Final Dilution of Anti-Rabbit IgG HRP (100X).....	15
Dilution of Wash Buffer	16
Assay Method.....	16
Typical Data	19
Limitations of the Procedure.....	20
Performance Characteristics	21
Sensitivity	21
Precision	22
Recovery.....	23
Parallelism	23
Linearity of Dilution	24
Specificity.....	25
References	27

INTRODUCTION

c-Met, a member of the tyrosine kinase superfamily, is the receptor for hepatocyte growth factor, also known as scatter factor (HGF/SF). The mature c-Met protein is a disulfide-linked heterodimer with $M_r=190$ kDa composed of a heavily glycosylated α subunit that is completely extracellular in localization, and a β subunit comprised of an extracellular ligand binding domain, a single transmembrane domain, and a cytoplasmic tyrosine kinase domain. c-Met is transcribed from a single open reading frame and translated into a protein precursor that is proteolytically cleaved, yielding the heterodimeric mature protein. Alternative splicing yields several c-Met isoforms, including proteins that remain in the uncleaved, monomeric state or that lack various portions of the c-Met cytoplasmic domain. Cells expressing c-Met include epithelial cells, endothelial cells, blood cells of various types, and glomerular mesenchymal cells.

The ligand for c-Met, HGF/SF, is a member of the plasminogen-related growth factor family, which is synthesized as an inactive pro-form. HGF/SF activation requires cleavage with either urokinase plasminogen activator (uPA), HGF activator, or Coagulation Factor Xa. Sources of HGF/SF include mesenchymal cells, mesangial cells, endothelial cells, macrophages, and tumor cells.

HGF/SF binding to c-Met stimulates receptor dimerization and the phosphorylation of numerous residues within the receptor's cytoplasmic domain, including tyrosines 1230, 1234, and 1235 within the Tyr-X-X-X Tyr-Tyr motif of c-Met's activation loop. This motif is conserved among the activation loops of several receptor tyrosine kinases including insulin receptor, insulin-like growth factor-1 receptor, nerve growth factor receptor/Trks, and RON. Phosphorylation of

tyrosines 1234 and 1235 of c-Met is required for activation of the receptor's tyrosine kinase activity. c-Met phosphorylation also generates docking sites for numerous signaling molecules and stimulates receptor internalization via clathrin-coated vesicles. Signaling proteins that are phosphorylated and/or localized in response to c-Met phosphorylation include: Grb2, Shc, Cbl, Crk, cortactin, paxillin, GAB1, PI-3 K, FAK, Src, Ras, ERK1 and 2, JNK, PLC- γ , AKT, and STAT3.

The Invitrogen c-Met [pYpYpY1230/1234/1235] ELISA is designed to detect and quantify the level of c-Met protein that is phosphorylated at tyrosines 1230, 1234, and 1235. This assay is intended for the detection of c-Met from lysates of human cells. For normalizing c-Met content of the samples, a c-Met (Total) ELISA Kit, which is independent of phosphorylation status, is available from Invitrogen (Cat. # KHO0251).

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

READ ENTIRE PROTOCOL BEFORE USE

PRINCIPLE OF THE METHOD

The Invitrogen c-Met [pYpYpY1230/1234/1235] kit is a solid phase sandwich Enzyme Linked-Immuno-Sorbent Assay (ELISA). A monoclonal antibody specific for c-Met has been coated onto the wells of the microtiter strips provided. Samples, including a standard containing c-Met [pYpYpY1230/1234/1235], control specimens, and unknowns, are pipetted into these wells. During the first incubation, the c-Met antigen binds to the immobilized (capture) antibody. After washing, a rabbit antibody specific for c-Met phosphorylated at tyrosines 1230, 1234, and 1235 is added to the wells. During the second incubation, this antibody serves as a detection antibody by binding to the immobilized c-Met protein captured during the first incubation. After removal of excess detection antibody, a horseradish peroxidase-labeled Anti-Rabbit IgG (Anti-Rabbit IgG-HRP) is added. This binds to the detection antibody to complete the four-member sandwich. After a third incubation and washing to remove all the excess anti-rabbit IgG-HRP, a substrate solution is added, which is acted upon by the bound enzyme to produce color. The intensity of this colored product is directly proportional to the concentration of c-Met [pYpYpY1230/1234/1235] present in the original specimen.

REAGENTS PROVIDED

Note: Store all reagents at 2 to 8°C.

<i>Reagent</i>	<i>96 Test Kit</i>
<i>Hu c-Met [pYpYpY1230/1234/1235] Standard:</i> Lyophilized. Contains 0.1% sodium azide. Refer to vial label for quantity and reconstitution volume.	2 vials
<i>Standard Diluent Buffer.</i> Contains 0.1% sodium azide; 25 mL per bottle.	1 bottle
<i>Antibody Coated Wells, 12x8 Well Strips.</i>	1 plate
<i>Hu c-Met [pYpYpY1230/1234/1235] Detection Antibody.</i> Contains 0.1% sodium azide; 11 mL per bottle.	1 bottle
<i>Anti-Rabbit IgG HRP (100X).</i> Contains 3.3 mM thymol; 0.125 mL per vial.	1 vial
<i>HRP Diluent.</i> Contains 3.3 mM thymol; 25 mL per bottle.	1 bottle
<i>Wash Buffer Concentrate (25X);</i> 100 mL per bottle.	1 bottle
<i>Stabilized Chromogen, Tetramethylbenzidine (TMB);</i> 25 mL per bottle.	1 bottle
<i>Stop Solution;</i> 25 mL per bottle.	1 bottle
<i>Plate Covers, adhesive strips.</i>	3

Disposal Note: This kit contains materials with small quantities of sodium azide. Sodium azide reacts with lead and copper plumbing to form explosive metal azides. Upon disposal, flush drains with a large volume of water to prevent azide accumulation. Avoid ingestion and contact with eyes, skin and mucous membranes. In case of contact, rinse affected area with plenty of water. Observe all federal, state and local regulations for disposal.

SUPPLIES REQUIRED BUT NOT PROVIDED

1. Microtiter plate reader capable of measurement at or near 450 nm.
2. Calibrated adjustable precision pipettes, preferably with disposable plastic tips. (A manifold multi-channel pipette is desirable for large assays.)
3. Cell extraction buffer (see Recommended Formulation, p.11).
4. Distilled or deionized water.
5. Plate washer: automated or manual (squirt bottle, manifold dispenser, etc.).
6. Data analysis and graphing software. Graph paper: linear (Cartesian), log-log, or semi-log, as desired.
7. Glass or plastic tubes for diluting and aliquoting standard.
8. Absorbent paper towels.
9. Calibrated beakers and graduated cylinders in various sizes.

PROCEDURAL NOTES/LAB QUALITY CONTROL

1. When not in use, kit components should be refrigerated. All reagents should be warmed to room temperature before use.
2. **Microtiter plates should be allowed to come to room temperature before opening the foil bag.** Once the desired number of strips has been removed, immediately reseal the bag and store at 2 to 8°C to maintain plate integrity.

3. Samples should be frozen if not analyzed shortly after collection. Avoid multiple freeze-thaw cycles of frozen samples. Thaw completely and mix well prior to analysis.
4. If large amounts of particulate matter are present, centrifuge or filter prior to analysis.
5. It is recommended that all standards, controls and samples be run in duplicate.
6. Extracted cell lysate samples containing c-Met [pYpYpY1230/1234/1235] protein should be diluted with *Standard Diluent Buffer* at least 1:10. This dilution is necessary to reduce the matrix effect of the Cell Extraction Buffer.
7. When pipetting reagents, maintain a consistent order of addition from well-to-well. This ensures equal incubation times for all wells.
8. Cover or cap all reagents when not in use.
9. **Do not mix or interchange different reagent lots from various kit lots.**
10. Do not use reagents after the kit expiration date.
11. Read absorbances within 2 hours of assay completion.
12. In-house controls should be run with every assay. If control values fall outside pre-established ranges, the accuracy of the assay is suspect.
13. All residual wash liquid must be drained from the wells by efficient aspiration or by decantation followed by tapping the plate forcefully on absorbent paper. *Never* insert absorbent paper directly into the wells.
14. Because *Stabilized Chromogen* is light sensitive, avoid prolonged exposure to light. Also avoid contact between *Stabilized Chromogen* and metal, or color may develop.

SAFETY

All blood components and biological materials should be handled as potentially hazardous. Follow universal precautions as established by the Centers for Disease Control and Prevention and by the Occupational Safety and Health Administration when handling and disposing of infectious agents.

DIRECTIONS FOR WASHING

Incomplete washing will adversely affect the test outcome. All washing must be performed with *Wash Buffer Concentrate (100X)* provided.

Washing can be performed manually as follows: completely aspirate the liquid from all wells by gently lowering an aspiration tip (aspiration device) into the bottom of each well. Take care not to scratch the inside of the well.

After aspiration, fill the wells with at least 0.4 mL of diluted wash solution. Let soak for 15 to 30 seconds, then aspirate the liquid. Repeat as directed under **ASSAY METHOD**. After the washing procedure, the plate is inverted and tapped dry on absorbent tissue.

Alternatively, the wash solution may be put into a squirt bottle. If a squirt bottle is used, flood the plate with wash buffer, completely filling all wells. After the washing procedure, the plate is inverted and tapped dry on absorbent tissue.

If using an automated washer, the operating instructions for washing equipment should be carefully followed. If your automated washer allows, 30 second soak cycles should be programmed into the wash cycle.

PROCEDURE FOR EXTRACTION OF PROTEINS FROM CELLS

A. Recommended Formulation of Cell Extraction Buffer:

10 mM Tris, pH 7.4

100 mM NaCl

1 mM EDTA

1 mM EGTA

1 mM NaF

20 mM $\text{Na}_4\text{P}_2\text{O}_7$

2 mM Na_3VO_4

1% Triton X-100

10% glycerol

0.1% SDS

0.5% deoxycholate

1 mM PMSF (stock is 0.3 M in DMSO)

Protease inhibitor cocktail (e.g., Sigma Cat. # P-2714, reconstituted according to manufacturer's guideline). Add 250 μL per 5 mL Cell Extraction Buffer.

This buffer is stable for 2 to 3 weeks at 4°C or for up to 6 months when aliquoted (without protease inhibitors and PMSF added) and stored at -20°C. When stored frozen, the Cell Extraction Buffer should be thawed on ice. This buffer (minus protease inhibitor cocktail and PMSF) can be obtained from Invitrogen (Cat. # FNN0011). **Important:** add the protease inhibitors just before using. The stability of protease inhibitor supplemented Cell Extraction Buffer is 24 hours at 4°C. PMSF is very unstable and must be added prior to use, even if added previously.

B. Protocol for Cell Extraction

This protocol has been applied to several cell lines with the Cell Extraction Buffer above. Researchers may optimize the cell extraction procedures that work best in their hands.

1. Collect cells in PBS by centrifugation (non-adherent) or scraping from culture flasks (adherent).
2. Wash cells twice with cold PBS.
3. Remove and discard the supernatant and collect the cell pellet. (At this point the cell pellet can be frozen at -80°C and lysed at a later date).
4. Lyse the cell pellet in Cell Extraction Buffer for 30 minutes on ice with vortexing at 10 minute intervals. The volume of Cell Extraction Buffer depends on the cell number in cell pellet and expression of c-Met [pYpYpY1230/1234/1235]. For example, 2×10^7 GTL 16 cells grown in RPMI plus 10% FBS treated with sodium orthovanadate at 1 mM for 16 hours can be extracted in 0.5 mL of Extraction Buffer. Under these conditions, use of 1-10 μL of the clarified cell extract diluted to a volume of 100 μL /well in *Standard Diluent Buffer* (See **Assay Method**) is sufficient for the detection of c-Met [pYpYpY1230/1234/1235].
5. Transfer extracts to microcentrifuge tubes and centrifuge at 13,000 rpm for 10 minutes at 4°C .
6. Aliquot the clear lysate to clean microfuge tubes. These samples are ready for assay. Lysates can be stored at -80°C . Avoid multiple freeze-thaw cycles.

REAGENT PREPARATION AND STORAGE

A. Reconstitution and Dilution of Hu c-Met [pYpYpY1230/1234/1235] Standard

Note: This *Hu c-Met [pYpYpY1230/1234/1235] Standard* (lyophilized cell extract from GTL 16 cells) is designated in Units/mL. One Unit of standard is equivalent to the amount of c-Met [pYpYpY1230/1234/1235] derived from ~21,000 GTL 16 cells treated with 1 mM sodium orthovanadate for 16 hours.

1. Reconstitute *Hu c-Met [pYpYpY1230/1234/1235] Standard* with *Standard Diluent Buffer*. Refer to standard vial label for instructions. Swirl or mix gently and allow to sit for 10 minutes to ensure complete reconstitution. Label as 100 Units/mL c-Met [pYpYpY1230/1234/1235]. Use the standard within 1 hour of reconstitution.
2. Add 0.25 mL of *Standard Diluent Buffer* to each of 6 tubes labeled 50, 25, 12.5, 6.25, 3.13, and 1.56 Units/mL of c-Met [pYpYpY1230/1234/1235].
3. Make serial dilutions of the standard as described in the following dilution table. Mix thoroughly between steps.

B. Dilution of Hu c-Met [pYpYpY1230/1234/1235] Standard

Standard:	Add:	Into:
100 Units/mL	Prepare as described in step 1	
50 Units/mL	0.25 mL of the 100 Units/mL std.	0.25 mL of the Diluent Buffer
25 Units/mL	0.25 mL of the 50 Units/mL std.	0.25 mL of the Diluent Buffer
12.5 Units/mL	0.25 mL of the 25 Units/mL std.	0.25 mL of the Diluent Buffer
6.25 Units/mL	0.25 mL of the 12.5 Units/mL std.	0.25 mL of the Diluent Buffer
3.13 Units/mL	0.25 mL of the 6.25 Units/mL std.	0.25 mL of the Diluent Buffer
1.56 Units/mL	0.25 mL of the 3.13 Units/mL std.	0.25 mL of the Diluent Buffer
0 Units/mL	0.25 mL of the Diluent Buffer	An empty tube

Remaining reconstituted standard should be discarded or frozen in aliquots at -80°C for further use. Do not subject reconstituted standards to more than 1 freeze-thaw cycle.

C. Storage and Final Dilution of Anti-Rabbit IgG HRP (100X)

Please Note: The *Anti-Rabbit IgG HRP (100X)* is in 50% glycerol. This solution is viscous. To ensure accurate dilution, allow *Anti-Rabbit IgG-HRP (100X)* to reach room temperature. Gently mix. Pipette *Anti-Rabbit IgG HRP (100X)* slowly. Remove excess concentrate solution from pipette tip by gently wiping with clean absorbent paper.

1. Dilute 10 μL of this 100X concentrated solution with 1 mL of *HRP Diluent* for each 8-well strip used in the assay. Label as Anti-Rabbit IgG HRP Working Solution.

<u># of 8-Well Strips</u>	<u>Volume of Anti-Rabbit IgG HRP (100X)</u>	<u>Volume of Diluent</u>
2	20 μL solution	2 mL
4	40 μL solution	4 mL
6	60 μL solution	6 mL
8	80 μL solution	8 mL
10	100 μL solution	10 mL
12	120 μL solution	12 mL

For Example:

2. Return the unused *Anti-Rabbit IgG HRP (100X)* to the refrigerator.

D. Dilution of Wash Buffer

Allow the *Wash Buffer Concentrate (25X)* to reach room temperature and mix to ensure that any precipitated salts have redissolved. Dilute 1 volume of the *Wash Buffer Concentrate (25X)* with 24 volumes of deionized water (i.e., 50 mL may be diluted up to 1.25 liters; 100 mL may be diluted up to 2.5 liters). Label as Working Wash Buffer.

Store both the concentrate and the Working Wash Buffer in the refrigerator. The diluted buffer should be used within 14 days.

ASSAY METHOD: PROCEDURE AND CALCULATIONS

Be sure to read the *Procedural Notes/Lab Quality Control* section before carrying out the assay.

Allow all reagents to reach room temperature before use. Gently mix all liquid reagents prior to use.

Note: A standard curve must be run with each assay.

1. Determine the number of 8-well strips needed for the assay. Insert these in the frame(s) for current use. (Re-bag extra strips and frame. Store these in the refrigerator for future use.)
2. Add 100 μL of the *Standard Diluent Buffer* to zero wells. Well(s) reserved for chromogen blank should be left empty.
3. Add 100 μL of standards, samples or controls to the appropriate microtiter wells. Samples prepared in Cell Extraction Buffer must be diluted 1:10 or greater in *Standard Diluent Buffer* (for example, 10 μL sample into 90 μL buffer). While a 1:10 sample dilution has been found to be satisfactory, higher dilutions such as 1:25 or 1:50 may be optimal. The dilution

chosen should be optimized for each experimental system. Tap gently on side of plate to thoroughly mix. (See **REAGENT PREPARATION AND STORAGE**, Section B.)

4. Cover plate with *plate cover* and incubate for **2 hours at room temperature**.
5. Thoroughly aspirate or decant solution from wells and discard the liquid. Wash wells 4 times. See **DIRECTIONS FOR WASHING**.
6. Pipette 100 μ L of *Hu c-Met [pYpYpY1230/1234/1235] Detection Antibody* solution into each well except the chromogen blank(s). Tap gently on the side of the plate to mix.
7. Cover plate with *plate cover* and incubate for **1 hour at room temperature**.
8. Thoroughly aspirate or decant solution from wells and discard the liquid. Wash wells 4 times. See **DIRECTIONS FOR WASHING**.
9. Add 100 μ L *Anti-Rabbit IgG HRP Working Solution* to each well except the chromogen blank(s). (Prepare the working dilution as described in **REAGENT PREPARATION AND STORAGE**, Section C.)
10. Cover plate with the *plate cover* and incubate for **30 minutes at room temperature**.
11. Thoroughly aspirate or decant solution from wells and discard the liquid. Wash wells 4 times. See **DIRECTIONS FOR WASHING**.
12. Add 100 μ L of *Stabilized Chromogen* to each well. The liquid in the wells will begin to turn blue.

13. Incubate for **30 minutes at room temperature and in the dark.** **Please Note: Do not cover the plate with aluminum foil or metalized mylar.** The incubation time for chromogen substrate is often determined by the microtiter plate reader used. Many plate readers have the capacity to record a maximum optical density (O.D.) of 2.0. The O.D. values should be monitored and the substrate reaction stopped before the O.D. of the positive wells exceed the limits of the instrument. The O.D. values at 450 nm can only be read after the *Stop Solution* has been added to each well. If using a reader that records only to 2.0 O.D., stopping the assay after 20 to 25 minutes is suggested.
14. Add 100 μL of *Stop Solution* to each well. Tap side of plate gently to mix. The solution in the wells should change from blue to yellow.
15. Read the absorbance of each well at 450 nm having blanked the plate reader against a chromogen blank composed of 100 μL each of *Stabilized Chromogen* and *Stop Solution*. Read the plate within 2 hours after adding the *Stop Solution*.
16. Plot on graph paper the absorbance of the standards against the standard concentration. (Optimally, the background absorbance may be subtracted from *all* data points, including standards, unknowns and controls, prior to plotting.) Draw the best smooth curve through these points to construct the standard curve. If using curve fitting software, the four parameter algorithm provides the best curve fit.
17. Read the c-Met [pYpYpY1230/1234/1235] concentrations for unknown samples and controls from the standard curve plotted in step 16. **Multiply value(s) obtained for sample(s) by dilution factor to correct for the dilution in step 3.** (Samples still

producing signals higher than the highest standard (100 Units/mL) should be further diluted in *Standard Diluent Buffer* and reanalyzed, multiplying the concentration found by the appropriate dilution factor.)

TYPICAL DATA

The following data were obtained for the various standards over the range of 0 to 100 Units/mL c-Met [pYpYpY1230/1234/1235].

Standard c-Met [pYpYpY1230/1234/1235] (Units/mL)	Average Optical Density (450 nm)
100	2.69
50	1.72
25	1.12
12.5	0.78
6.25	0.55
3.13	0.41
1.56	0.31
0	0.20

LIMITATIONS OF THE PROCEDURE

Do not extrapolate the standard curve beyond the 100 Units/mL standard point; the dose-response is non-linear in this region and accuracy is difficult to obtain. Dilute samples >100 Units/mL with *Standard Diluent Buffer*; reanalyze these and multiply results by the appropriate dilution factor.

The influence of various extraction buffers has not been thoroughly investigated. The rate of degradation of native c-Met or dephosphorylation of c-Met [pYpYpY1230/1234/1235] in various matrices has not been investigated. Although c-Met degradation or dephosphorylation of c-Met [pYpYpY1230/1234/1235] in the Cell Extraction Buffer described in this protocol has not been seen to date, the possibility of this occurrence cannot be excluded.

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

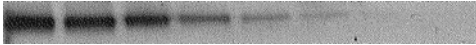
PERFORMANCE CHARACTERISTICS

SENSITIVITY

The analytical sensitivity of this assay is <0.25 Units/mL of c-Met [pYpYpY1230/1234/1235]. This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 30 times. This level of sensitivity is approximately equivalent to the c-Met [pYpYpY1230/1234/1235] content of 5,000 cells, lysed as described above.

The sensitivity of this ELISA was compared to Western blotting using known quantities of c-Met [pYpYpY1230/1234/1235]. The data presented in Figure 1 show that the sensitivity of the ELISA is approximately 2x greater than that of Western blotting. The bands shown in the Western blot data were developed using rabbit anti-c-Met [pYpYpY1230/1234/1235] and an alkaline phosphatase conjugated anti-rabbit IgG followed by chemiluminescent substrate and autoradiography.

Figure 1: Detection of c-Met [pYpYpY1230/1234/1235] by ELISA vs Western Blot:

Western Blot (140 kDa)								
ELISA (O.D. 450 nm)	2.690	1.719	1.179	0.784	0.552	0.414	0.306	0.204
c-Met [pYpYpY 1230/1234/1235] (Units/test)	10	5	2.5	1.25	0.625	0.313	0.156	0

PRECISION

1. Intra-Assay Precision

Samples of known c-Met [pYpYpY1230/1234/1235] concentration were assayed in replicates of 16 to determine precision within an assay.

	Sample 1	Sample 2	Sample 3
Mean (Units/mL)	68.6	25.3	10.8
SD	2.8	2.0	0.6
%CV	4.1	7.8	5.4

SD = Standard Deviation

CV = Coefficient of Variation

2. Inter-Assay Precision

Samples were assayed 48 times in multiple assays to determine precision between assays.

	Sample 1	Sample 2	Sample 3
Mean (Units/mL)	67.8	25.7	10.3
SD	3.3	2.0	0.9
%CV	4.8	7.9	8.8

SD = Standard Deviation

CV = Coefficient of Variation

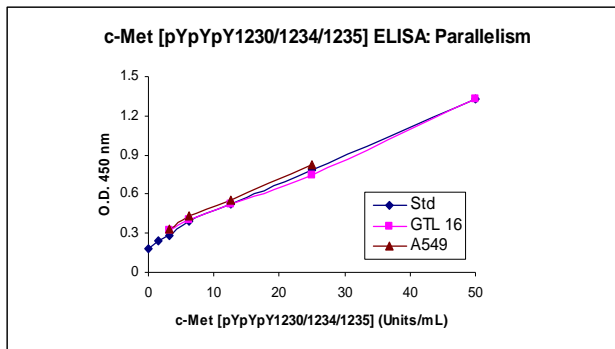
RECOVERY

To evaluate recovery, extraction buffer was diluted 1:10 with *Standard Diluent Buffer* to bring the SDS concentration to <0.01%. c-Met [pYpYpY1230/1234/1235] Standard was spiked into the cell extraction buffer. The average recovery was 107%.

PARALLELISM

Natural c-Met [pYpYpY1230/1234/1235] from sodium orthovanadate-treated GTL 16 and A549 cell lysates was serially diluted in *Standard Diluent Buffer*. The optical density of each dilution was plotted against the c-Met [pYpYpY1230/1234/1235] standard curve. Parallelism was demonstrated by the figure below and indicated that the standard accurately reflects c-Met [pYpYpY1230/1234/1235] content in samples.

Figure 2



LINEARITY OF DILUTION

GTL 16 cells grown in tissue culture medium containing 10% fetal calf serum were treated with sodium orthovanadate at 1 mM for 16 hours and lysed with Cell Extraction Buffer. This lysate was diluted in *Standard Diluent Buffer* over the range of the assay and measured for c-Met [pYpYpY1230/1234/1235] content. Linear regression analysis of samples versus the expected concentration yielded a correlation coefficient of 0.99.

Dilution	Cell Lysate		
	Measured (Units/mL)	Expected (Units/mL)	% Expected
Neat	12.8	12.8	100
1/2	6.9	6.4	108
1/4	3.4	3.2	106
1/8	2.0	1.6	120

SPECIFICITY

The specificity of this assay for c-Met phosphorylated at tyrosines 1230, 1234, and 1235 was confirmed by peptide competition. The data presented in Figure 3 show that the phospho-peptide containing the phosphorylated tyrosines could block the ELISA signal. The same c-Met sequences without phosphate group could not block the ELISA signal.

Figure 3

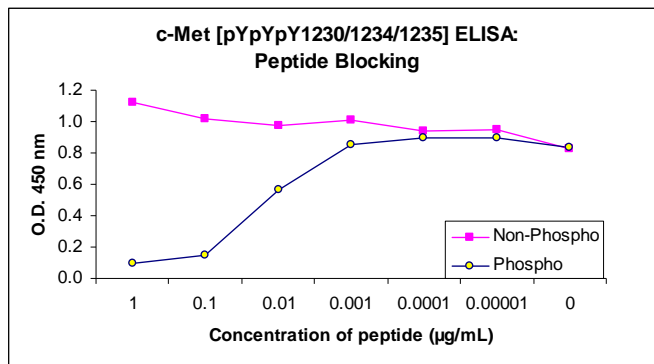


Figure 4

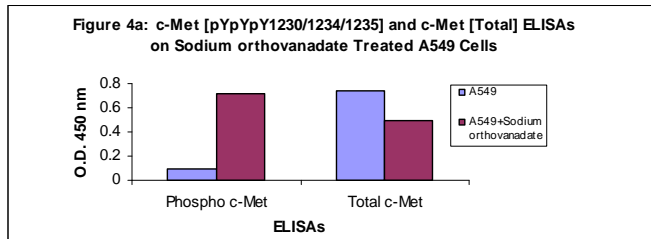
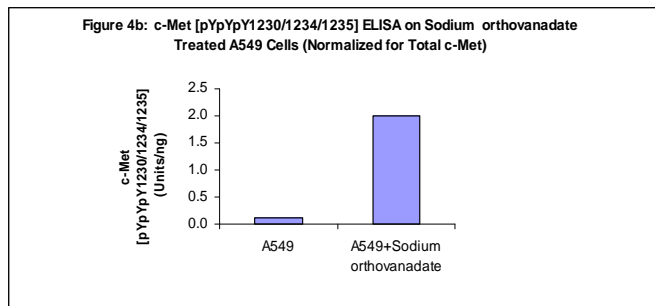


Figure 4a shows that treatment with sodium orthovanadate, a tyrosine phosphatase inhibitor, increased phosphorylation at 1230, 1234, and 1235 of c-Met in A549 cells. A549 cells were treated with sodium orthovanadate at 1 mM for 16 hours, lysed, and assayed in parallel for both c-Met [total] amount and c-Met [pYpYpY1230/1234/1235]. The amount of c-Met remained comparable, while the levels of phosphorylation at tyrosines 1230, 1234, and 1235 increased with sodium orthovanadate treatment. The level of c-Met phosphorylation has been normalized for total c-Met content (Figure 4b).











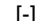
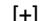



REFERENCES

1. Crostella, L., et al. (2001) Hepatocyte Growth Factor/scatter factor-induces phosphorylation of cortactin in A431 cells in a Src kinase-independent manner. *Oncogene* 20(28):3735-3745.
2. Fan, S., et al. (2001) The multisubstrate adapter Gab1 regulates hepatocyte growth factor (scatter factor)-c-Met signaling for cell survival and DNA repair. *Mol. Cell. Biol.* 21(15):4968-4984.
3. Follenzi, A., et al. (2000) Cross-talk between the proto-oncogenes Met and Ron. *Oncogene* 19(27):3041-3049.
4. Furge, K.A., et al. (2000) Met receptor tyrosine kinase: enhanced signaling through adapter proteins. *Oncogene* 19(49):5582-5589.
5. Guiton, M., et al. (1994) Identification of in vivo brain-derived neurotrophic factor-stimulated autophosphorylation sites on the TrkB receptor tyrosine kinase by site-directed mutagenesis. *J. Biol. Chem.* 269(48):30370-30377.

6. Longati, P., et al. (2001) Receptor tyrosine kinases as therapeutic targets: the model of the MET oncogene. *Curr. Drug Targets* 2(1):41-55.
7. Kang, J.Y, et al. (2003) Tissue microarray analysis of hepatocyte growth factor/Met pathway components reveals a role from Met, matrilin, and hepatocyte growth factor activator inhibitor 1 in the progression of node-negative breast cancer. *Cancer Res.* 63:1101-1105.
8. Matsumoto, K. and T. Nakamura (2001) Hepatocyte growth factor: renoprotective role and potential therapeutics for renal diseases. *Kidney Int.* 59(6):2023-2038.
9. Maulik, G., et al. (2002) Role of the hepatocyte growth factor receptor, c-Met, in oncogenesis and potential for therapeutic inhibition. *Cytokine Growth Factor Rev.* 13(1):41-59.
10. Park, M., et al. (1987) Sequence of MET protooncogene cDNA has features characteristic of the tyrosine kinase family of growth-factor receptors. *Proc. Nat'l. Acad. Sci.* 84:6379-6383.
11. Parr, C., et al. (2001) The HGF/SF-induced phosphorylation of paxillin, matrix adhesion, and invasion of prostate cancer cells were suppressed by NK4, an HGF/SF variant. *Biochem. Biophys. Res. Commun.* 285(5):1330-1337.
12. Wei, L., et al. (1995) Expression, characterization, and crystallization of the catalytic core of the human insulin receptor protein-tyrosine kinase domain. *J. Biol. Chem.* 270(14):8122-8130.
13. Zhang, X., et al. (2003) Sp1 and Sp3 transcription factors synergistically regulate HGF receptor gene expression in kidney. *Am. J. Renal Physiol.* 284:F82-F94.

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.

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

Copyright © Invitrogen Corporation, 15 January 2010

NOTES

Hu c-Met [pYpYpY1230/1234/1235] Assay Summary

Incubate 100 μ L Standard or Cell Extract (>1:10)
for 2 hours at RT



↓ aspirate and wash 4x

Incubate 100 μ L of Detection Antibody
for 1 hour at RT



↓ aspirate and wash 4x

Incubate 100 μ L of HRP Anti-Rabbit Antibody
for 30 minutes at RT



↓ aspirate and wash 4x

Incubate 100 μ L of Stabilized Chromogen
for 30 minutes at RT



↓
Add 100 μ L of Stop Solution and read at 450 nm



Total time: 4 hours

 c-Met
[pYpYpY1230/1234/1235]

 HRP Anti-Rabbit
Antibody

 Detection
Antibody