

> TGF- β 1 EIA kit

Catalog # ADI-900-155

96 Well Enzyme Immunoassay Kit

For use with plasma, serum and tissue culture supernatants



All reagents, except standard, should be stored at 4°C. Store standard at -20°C.



Check our website for additional protocols, technical notes and FAQs.



For proper performance, use the insert provided with each individual kit received.

Table of Contents

2 Introduction

2 Principle

3 Materials Supplied

4 Storage

4 Materials Needed but Not Supplied

5 Reagent Preparation

6 Sample Handling

6 Sample Activation

8 Assay Procedure

9 Calculation of Results

9 Calibration

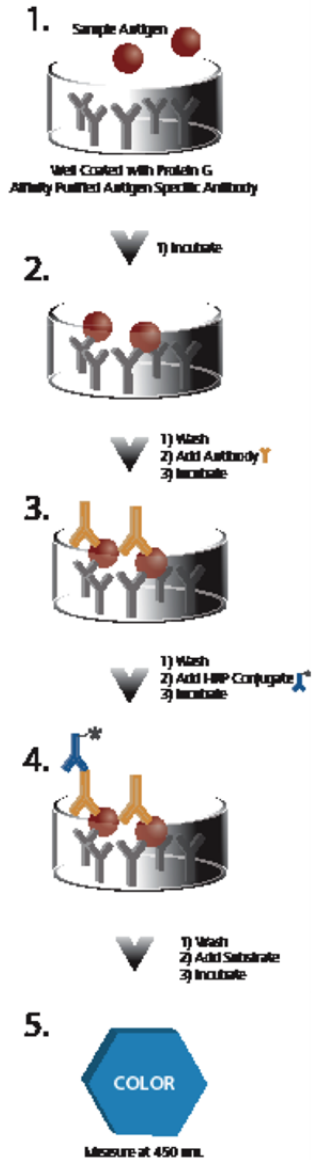
10 Typical Results in Assay Buffer 13

11 Typical Results in Assay Buffer 30

12 Performance Characteristics

14 References

16 Limited Warranty



Introduction

Transforming growth factor beta 1 is also known as TGFβ1, TGF β, and TGFβ. The inactive form consists of a TGF-β1 homodimer that is non-covalently linked to a latency-associated peptide homodimer. The active form is a homodimer of mature TGF-β1 that is disulfide linked. The precursor is cleaved into mature TGF-β1 and the latency-associated peptide^{1, 2}. Many cells can synthesize TGF-β1, which is secreted to function in proliferation, differentiation, transformation, signaling and apoptosis. This cytokine has been implicated in many diseases including diabetes³, renal disease⁴, chronic pulmonary obstructive disease⁵, as well as many cancers including prostate⁶, and colon⁷. This protein has generated tremendous excitement as a potential tumor-specific cancer therapeutic. Specifically, as a stable soluble dimer, it selectively induces apoptosis in many transformed cells but not in normal cells⁸. The literature contains many recent reviews on the biology and function of TGF-β1^{4, 6, 9, 10}.

Principle

1. Samples and standards are added to wells coated with a monoclonal antibody specific for TGF-β1. The plate is then incubated.
2. The plate is washed, leaving only bound TGF-β1 on the plate. A yellow solution of polyclonal antibody to TGF-β1 is then added. This binds the TGF-β1 captured on the plate. The plate is then incubated.
3. The plate is washed to remove excess antibody. A blue solution of HRP conjugate is added to each well, binding to the TGF-β1 polyclonal. The plate is again incubated.
4. The plate is washed to remove excess HRP conjugate. TMB Substrate solution is added. The substrate generates a blue color when catalyzed by the HRP.
5. Stop solution is added to stop the substrate reaction. The resulting yellow color is read at 450 nm. The amount of signal is directly proportional to the level of TGF-β1 in the sample



Do not mix components from different kit lots or use reagents beyond the expiration date of the kit.



The standard should be handled with care due to the known and unknown effects of the molecule.



6.01% urea is harmful if swallowed or inhaled. Causes irritation to skin, eyes, and respiratory tract.



Activity of conjugate is affected by nucleophiles such as azide, cyanide, and hydroxylamine.



Protect substrate from prolonged exposure to light.



Stop solution is caustic. Keep tightly capped.

Materials Supplied

1. Assay Buffer 13
110 mL, Catalog No. 80-1627
Tris buffered saline containing protein and detergents.
2. Assay Buffer 30
110 mL, Catalog No. 80-1610
Tris buffered saline and urea containing detergents.
3. human TGF- β 1 Standard
0.25 mL, Catalog No. 80-1577
One vial containing 10,000 pg/mL of recombinant human TGF- β 1.
4. TGF- β 1 Clear Microtiter Plate
One Plate of 96 Wells, Catalog No. 80-1574
A plate of break-apart strips coated with a mouse monoclonal antibody specific to TGF- β 1.
5. Wash Buffer Concentrate
100 mL, Catalog No. 80-1287
Tris buffered saline containing detergents.
6. human TGF- β 1 EIA Antibody
10 mL, Catalog No. 80-1575
A yellow solution of polyclonal antibody to TGF- β 1.
7. TGF- β 1 EIA Conjugate
10 mL, Catalog No. 80-1576
A blue solution of streptavidin conjugated to horseradish peroxidase.
8. TMB Substrate
10 mL, Catalog No. 80-0350
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide.
9. Stop Solution 2
10 mL, Catalog No. 80-0377
A 1N solution of hydrochloric acid in water.
10. TGF- β 1 Assay Layout Sheet
1 each, Catalog No. 30-0235
11. Plate Sealer
3 each, Catalog No. 30-0012



All reagents, except standard, should be stored at 4°C. Store standard at -20°C.

Storage

All components of this kit, except the TGF- β 1 Standard, are stable at 4°C until the kit's expiration date. The TGF- β 1 Standard must be stored at -20°C .

Materials Needed but Not Supplied

1. Deionized or distilled water.
2. Precision pipets for volumes between 50 μ L and 1,000 μ L.
3. Repeater pipet for dispensing 100 μ L.
4. Disposable beakers for diluting buffer concentrates.
5. Graduated cylinders.
6. A microplate shaker.
7. Lint-free paper for blotting.
8. Microplate reader capable of reading at 450 nm.
9. Graph paper for plotting the standard curve.
10. Glacial acetic acid (A.C.S. Grade, 17.4N).
11. HEPES, free acid (Reagent Grade, M.W. 238.8).
12. Urea (Reagent Grade, M.W. 60.06).
13. Sodium hydroxide (A.C.S. Grade, 10N).
14. Hydrochloric acid (A.C.S. Grade, 12N).



Bring all reagents, except for the standard, to room temperature for at least 30 minutes prior to opening.



Activation reagents may be stored at room temperature for up to one month.



Plastic tubes must be used for standard preparation.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.

Reagent Preparation

1. Wash Buffer

Prepare the wash buffer by diluting 50 mL of the supplied Wash Buffer Concentrate with 950 mL of deionized water. This can be stored at room temperature until the kit expiration, or for 3 months, whichever is earlier.

2. Activation Reagent

For cell culture supernatant

1N HCl, 50 mL

To 25 mL deionized water, slowly add 4.17 mL 12N HCl, mix, adjust volume to 50 mL using deionized water.

1.2N NaOH, 0.5M HEPES, 50 mL

To 30 mL deionized water, slowly add 6 mL 10N NaOH, mix, then add 5.95 g HEPES, stir well until all goes into solution. Bring final volume to 50 mL using deionized water.

For Serum/Plasma

2.5N Acetic Acid, 10M Urea, 50 mL

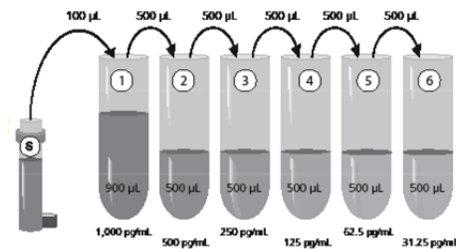
In a glass beaker, add 20 mL deionized water, then add 30.04 g urea, mix until urea dissolves. Slowly add 7.18 mL glacial acetic acid, mix well, bring final volume to 50 mL using deionized water. * The buffer can be stored at room temperature. When urea comes out of solution, place the bottle containing buffer into a 37°C water bath until dissolved.

2.7N NaOH, 1M HEPES, 50 mL

Slowly add 13.5 mL 10N NaOH to 25 mL deionized water, mix well. Add 11.9 g HEPES, mix well until HEPES dissolves. Bring final volume to 50 mL with deionized water.

3. human TGF-β1 Standards

Thaw standard on ice. Use the appropriate assay buffer based on the sample that will be tested. Use Assay Buffer 13 for activated cell culture supernatants and Assay Buffer 30 for activated serum/plasma samples.



Label six 12 x 75 mm polypropylene tubes #1 through #6. Pipet 900 µL of the appropriate assay buffer into tube #1. Pipet 500 µL of the assay buffer into tubes #2 through #6. Add 100 µL of the 10,000 pg/mL standard stock into tube #1 and vortex thoroughly. Add 500 µL of tube #1 to tube #2 and vortex thoroughly. Add 500 µL of tube #2 to tube #3 and vortex thoroughly. Continue this for tubes #4 through #6.

Diluted standards should be used within 60 minutes of preparation.

The concentration of TGF-β1 in tubes is labeled above.



If buffers other than those provided are used in the assay, the end-user must determine the appropriate dilution and assay validation.



If the end user chooses to vary from the activation procedure noted, it is up to the end user to determine the appropriate dilution of samples and assay validation.



Activated culture supernates must be used immediately. Activated serum and plasma are stable for up to one month at -80°C

Sample Handling

TGF- β 1, *in vivo*, is processed from a latent form to the bioactive form of the protein. Only the bioactive form is immunoreactive and detected in the kit. This process can be performed *in vitro*.

Serum supplements for cell culture contain TGF- β 1. It's recommended that a sample of serum-supplemented cell culture media be analyzed as a control or cell culture can be performed with serum-free media if possible.

To measure circulating levels of TGF- β 1 platelet-poor plasma should be used and care taken to avoid platelet degranulation. Platelets constitute a major source of TGF- β 1.

Sample Activation

Cell Culture Supernates

1. Add 20 μ L of 1N HCl to 100 μ L culture supernate.
2. Mix well. Incubate 10 minutes at room temperature.
3. Neutralize the sample with 20 μ L 1.2N NaOH/0.5M HEPES.
4. Mix well and assay immediately after dilution with assay buffer.

Sample is now at a 1:1.4 dilution and requires an additional 1:2 dilution with **Assay Buffer 13** before running the assay (mix 140 μ L activated sample with 140 μ L Assay Buffer 13). This will bring the samples to a final dilution of 1:2.8. A minimum of a 1:2.8 final dilution is required to eliminate matrix interference in the assay.

Serum or Plasma

1. Add 50 μ L 2.5N acetic acid/10M urea to 50 μ L serum / plasma.
2. Mix well. Incubate 10 minutes at room temperature.
3. Neutralize the sample with 50 μ L 2.7N NaOH/1M HEPES.
4. Add 350 μ L Assay Buffer 13.

Sample is now at a 1:10 dilution and requires an additional 1:12 dilution with **Assay Buffer 30** before running in the assay. This will bring the sample to a final dilution of 1:120. A minimum of a 1:120 final dilution is required to eliminate matrix interference in the assay.

Sample Diluents

Dilute the standards and activated samples in Assay Buffer 13 if activated cell culture supernatants will be tested in the assay. Otherwise use Assay Buffer 30 for the standard and activated serum/plasma samples.

Minimum Recommended Dilution After Sample Activation

Cell Culture supernatant	≥ 1:2
Serum	≥ 1:12
Plasma	≥ 1:12

Sample Recoveries

TGF-β1 concentrations were measured in activated-neutralized human serum samples. TGF-β1 was spiked into the activated samples and diluted with the assay buffer, then assayed in the kit.

The following results were obtained:

Sample	Recommended Dilution	%Recovery
Activated human serum	≥1:12	99.0%

Assay Procedure



Bring all reagents, except for the standard, to room temperature for at least 30 minutes prior to opening.



All standards, controls, and samples should be run in duplicate.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.



Pipet the reagents to the sides of the wells to avoid possible contamination.



Prior to the addition of antibody, conjugate, and substrate, ensure there is no residual wash buffer in the wells. Remaining wash buffer may cause variation in assay results.

Refer to the Assay Layout Sheet to determine the number of wells to be used. Remove the wells not needed for the assay and return them, with the desiccant, to the mylar bag and seal. Store unused wells at 4°C.

1. Pipet 100 μL of the appropriate assay buffer into the S0 (0 $\mu\text{g}/\text{mL}$ standard) wells.
2. Pipet 100 μL of Standards #1 through #6 to the bottom of the appropriate wells.
3. Pipet 100 μL of the samples to the bottom of the appropriate wells.
4. Seal the plate. Incubate for 1 hour on a plate shaker (~500 rpm) at room temperature.
5. Empty the contents of the wells and wash by adding 400 μL of wash buffer to every well. Repeat 3 more times for a total of 4 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
6. Pipet 100 μL of yellow antibody into each well except the blank.
7. Seal the plate. Incubate for 2 hours on a plate shaker (~500 rpm) at room temperature.
8. Wash as above (Step 5).
9. Add 100 μL of blue conjugate to each well except the blank.
10. Seal the plate. Incubate for 30 minutes on a plate shaker (~500 rpm) at room temperature.
11. Wash as above (Step 5).
12. Pipet 100 μL of substrate solution into each well.
13. Incubate for 30 minutes on a plate shaker (~500 rpm) at room temperature.
14. Pipet 100 μL of stop solution into each well.
15. After blanking the plate reader against the substrate blank, read optical density at 450 nm. If plate reader is not capable of adjusting for the blank, manually subtract the mean OD of the substrate blank from all readings.



Make sure to multiply sample concentrations by the dilution factor used during sample preparation.

Calculation of Results

Several options are available for the calculation of the concentration of TGF- β 1 in samples. We recommend that the data be handled by an immunoassay software package utilizing a 4-parameter logistic curve fitting program. If data reduction software is not readily available, the concentrations can be calculated as follows:

1. Calculate the average Net OD for each standard and sample by subtracting the average blank OD from the average OD for each standard and sample.

$$\text{Average Net OD} = \text{Average OD} - \text{Average Blank OD}$$

2. Using linear graph paper, plot the average Net OD for each standard versus TGF- β 1 concentration in each standard. Approximate a straight line through the points. The concentration of the unknowns can be determined by interpolation.

Samples with concentrations outside of the standard curve range will need to be re-analyzed using a different dilution.

Calibration

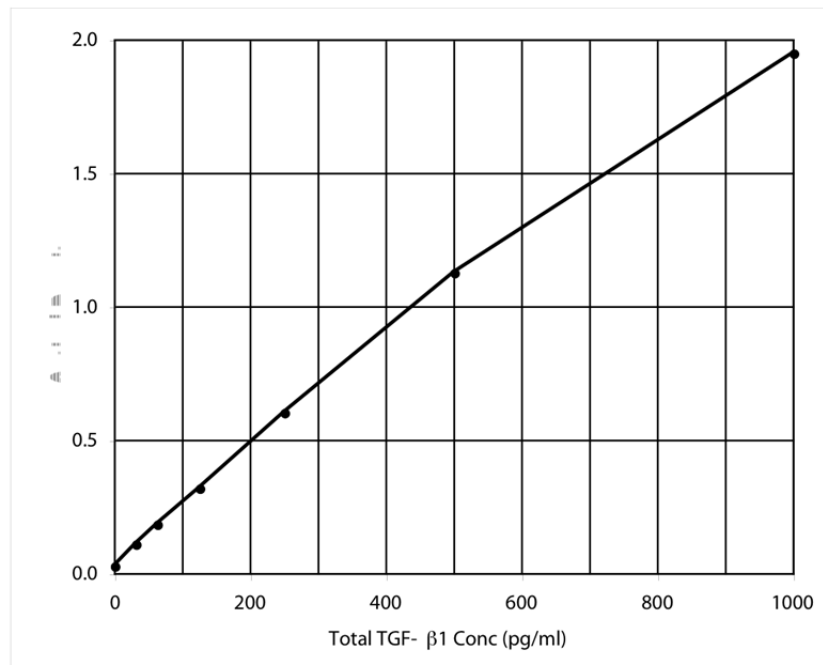
Calibration to the NIBSC/WHO TGF- β 1 International Reference Reagent 87/514 has been determined. To convert sample values obtained in the TGF- β 1 kit to this NIBSC/WHO TGF- β 1 Standard, use the equation below.

$$\text{NIBSC/WHO 87514 value (pg/mL)} = \text{Obtained human TGF-}\beta\text{1 value (pg/mL)} \times 0.915$$

Typical Results in Assay Buffer 13

The results shown below are for illustration only and should not be used to calculate results from another assay.

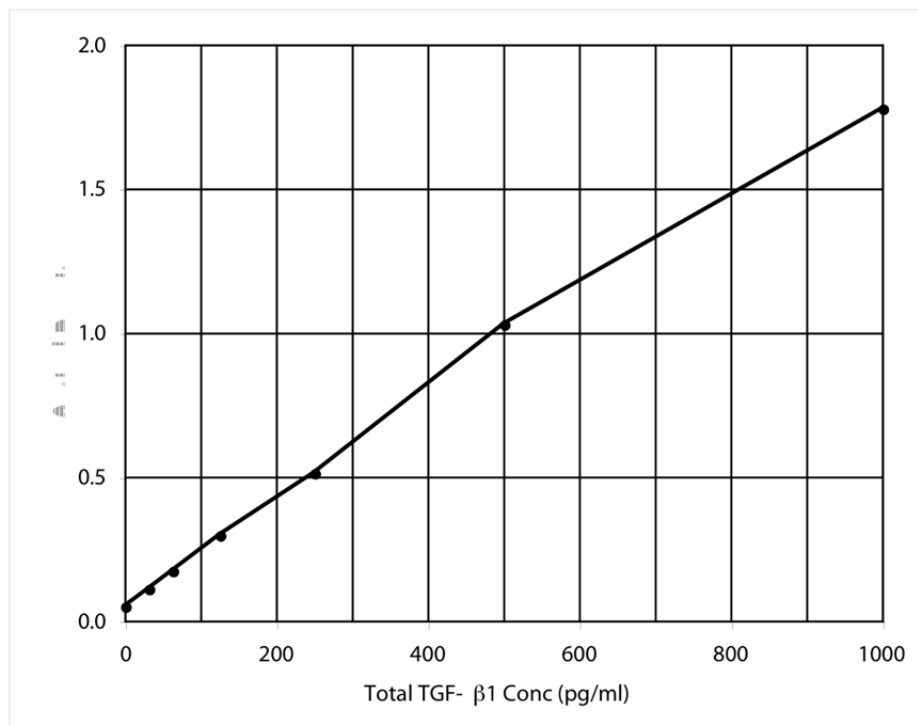
Sample	Net OD	TGF- β 1 (pg/mL)
S0	0.029	0
S1	1.950	1,000
S2	1.128	500
S3	0.604	250
S4	0.321	125
S5	0.186	62.5
S6	0.111	31.25
Unknown 1	1.172	526
Unknown 2	0.326	125



Typical Results in Assay Buffer 30

The results shown below are for illustration only and should not be used to calculate results from another assay.

Sample	Net OD	TGF- β 1 (pg/mL)
S0	0.052	0
S1	1.780	1,000
S2	1.031	500
S3	0.515	250
S4	0.299	125
S5	0.175	62.5
S6	0.113	31.25
Unknown 1	1.003	492
Unknown 2	0.268	115



Performance Characteristics

Specificity

The cross reactivities for a number of related compounds were determined by diluting cross reactants in the assay buffer at three different concentrations of 10,000,000 pg/mL, 100,000 pg/mL and 10,000 pg/mL. These samples were then measured in the assay.

Compound	Cross Reactivity
human TGF- β 1	100% (Assay Buffer 13 & Assay Buffer 30)
human TGF- β 1.2	2.23% (Assay Buffer 13)
	1.21% (Assay Buffer 30)
human TGF- β 2	<0.03% (Assay Buffer 13 & Assay Buffer 30)
human TGF- β 3	<0.03% (Assay Buffer 13 & Assay Buffer 30)
human latent TGF- β 1	<0.71% (Assay Buffer 13)
	<0.38% (Assay Buffer 30)
human LAP (latent associated protein)	<0.03% (Assay Buffer 13 & Assay Buffer 30)

Sensitivity

The sensitivity of the assay, defined as the concentration of TGF- β 1 measured at 2 standard deviations from the mean of 24 zeros along the standard curve, was determined to be 3.3 pg/mL in Assay Buffer 13, and 10.8 pg/mL in Assay Buffer 30.

Linearity

A buffer sample containing TGF- β 1 was serially diluted 1:2 in Assay Buffer 13 and measured in the assay. The results are shown in the table below.

Dilution	Expected (pg/mL)	Observed (pg/mL)	Recovery (%)
Neat	-----	789 pg/mL	-----
1:2	395 pg/mL	410 pg/mL	103.8 %
1:4	205 pg/mL	208 pg/mL	101.4 %
1:8	104 pg/mL	102 pg/mL	98.1 %
1:16	51 pg/mL	54 pg/mL	105.9 %
1:32	27 pg/mL	25 pg/mL	92.6 %

Another buffer sample containing TGF- β 1 was serially diluted 1:2 in Assay Buffer 30 and measured in the assay. The results are shown in the table below.

Dilution	Expected (pg/mL)	Observed (pg/mL)	Recovery (%)
Neat	-----	778 pg/mL	-----
1:2	389 pg/mL	368 pg/mL	94.6 %
1:4	184 pg/mL	185 pg/mL	100.5 %
1:8	93 pg/mL	105 pg/mL	112.9 %
1:16	52 pg/mL	52 pg/mL	100.0 %
1:32	26 pg/mL	28 pg/mL	107.7 %

Precision

Intra-assay precision was determined by assaying 24 replicates of three buffer controls containing TGF- β 1 in a single assay.

In Assay Buffer 13

pg/mL	%CV
98	8.5
251	3.4
475	3.9

In Assay Buffer 30

Inter-assay precision was determined by measuring buffer controls of varying TGF- β 1 concentrations in multiple assays over several days.

In Assay Buffer 13

pg/mL	%CV
110	9.4
250	5.9
481	5.3

In Assay Buffer 30

References

- 1 Russette M. Lyons, Jorma Keski-Oja, and Harold L. Moses.
Proteolytic Activation of Latent Transforming Growth Factor-beta from Fibroblast-conditioned Medium. *The journal of Cell Biology*. Volume 106, May 1998 1659-1665.
- 2 David A. Lawrence, Renee Pircher and Pierre Jullien.
Conversion of a high molecular weight latent beta-TGF from chicken embryo fibroblasts into a low molecular weight active beta-TGF under acidic conditons. *Biochemical and biophysical research communications*. Volume 133, No. 5, December 31, 1985.
- 3 Green EA, Gorelik L, McGregor CM, Tran EH, Flavell RA.
CD4+CD25+ T regulatory cells control anti-islet CD8+ T cells through TGF-beta-TGF-beta receptor interactions in type 1 diabetes. *Proc Natl Acad Sci U S A*. 2003 Sep 16;100(19):10878-83. Epub 2003 Aug 29.
- 4 Ziyadeh FN. Mediators of diabetic renal disease: the case for tgf-Beta as the major mediator. *J Am Soc Nephrol*. 2004 Jan;15 Suppl 1:S55-7. Review.
- 5 Pons AR, Sauleda J, Noguera A, Pons J, Barcelo B, Fuster A, Agusti AG.
Decreased macrophage release of TGF-{beta} and TIMP-1 in chronic obstructive pulmonary disease. *Eur Respir J*. 2005 Jul;26(1):60-6.
- 6 Danielpour D. Functions and regulation of transforming growth factor-beta (TGF-beta) in the prostate. *Eur J Cancer*. 2005 Apr;41(6):846-57. Review.
- 7 Li F, Cao Y, Townsend CM Jr, Ko TC. TGF-beta signaling in colon cancer cells. *World J Surg*. 2005 Mar;29(3):306-11.
- 8 Almasan A, Ashkenazi A. Apo2L/TRAIL: apoptosis signaling, biology, and potential for cancer therapy. *Cytokine Growth Factor Rev*. 2003 Jun-Aug;14(3-4):337-48.
- 9 Lebrin F, Deckers M, Bertolino P, Ten Dijke P. TGF-beta receptor function in the endothelium. *Cardiovasc Res*. 2005 Feb 15;65(3):599-608. Review.
- 10 Bachman KE, Park BH. Duel nature of TGF-beta signaling: tumor suppressor vs. tumor promoter. *Curr Opin Oncol*. 2005 Jan;17(1):49-54. Review.

Notes





Use of Product

This product contains research chemicals. As such, they should be used and handled only by or under the supervision of technically qualified individuals. This product is not intended for diagnostic or human use.

Warranty

Enzo Life Sciences International, Inc. makes no warranty of any kind, expressed or implied, which extends beyond the description of the product in this brochure, except that the material will meet our specifications at the time of delivery. Enzo Life Sciences International, Inc. makes no guarantee of results and assumes no liability for injuries, damages or penalties resulting from product use, since the conditions of handling and use are beyond our control.

www.enzolifesciences.com
Enabling Discovery in Life Science®

Global Headquarters

Enzo Life Sciences Inc.

10 Executive Blvd
Farmingdale, NY 11735

(p) 1-800-942-0430

(f) 1-631-694-7501

(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach
CH-4415 Lause / Switzerland

(p) +41/0 61 926 89 89

(f) +41/0 61 926 89 79

(e) info-ch@enzolifesciences.com

Please visit our website at www.enzolifesciences.com for additional contact information.