



Product Manual

Dkk-1 (human), ELISA kit

Catalog #: ADI-900-151

96 Well Enzyme-Linked Immunsorbent Assay Kit



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

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

TABLE OF CONTENTS



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.



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

Introduction	2
Principle	3
Materials Supplied	4
Storage	5
Other Materials Needed	5
Reagent Preparation	6
Sample Handling	7
Assay Procedure	8
Calculation of Results	9
Typical Results	10
Typical Standard Curve	10
Performance Characteristics	11
References	13
Contact Information	14

INTRODUCTION

The Dkk-1 (human), EIA kit is a complete kit for the quantitative determination of Dkk-1 in plasma, serum, and culture fluids. Please read the complete kit insert before performing this assay.

Dkk-1 (Dickkopf or Dickkopf-1) is a secreted inhibitor of the Wnt/canonical signaling pathway which mediates a variety of cellular processes. Wnt proteins are extracellular signaling molecules that initiate activation of the canonical pathway by binding low-density lipoprotein receptor-related protein 5&6 (LRP5/6) coreceptors, triggering a cascade of events that stabilize β -catenin¹. Wnts mediate osteoblast development and cell-to-cell interactions during embryogenesis, and are also responsible for regulating bone mass in adults during normal bone remodeling processes. Wnts are also found to be expressed during the stationary phase of the cell cycle and are thought to control expansion of bone marrow stromal cells preventing overpopulation in marrow².

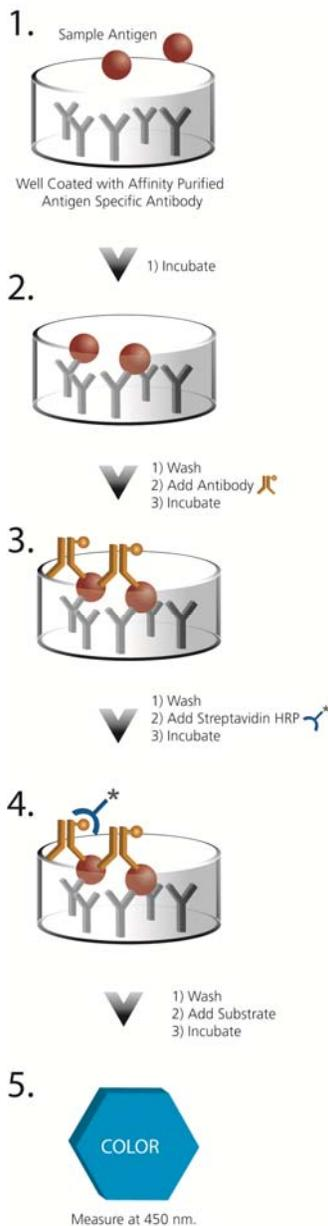
There are at least two families of secreted inhibitors of Wnt signaling: the secreted frizzled-related protein family and the Dickkopf family³. The Dkk-1 protein consists of 266 amino acids including a 31 amino acid N-terminal signal peptide. Dkk-1 has a predicted molecular weight of 26kD and an apparent molecular weight of 35kD by SDS-PAGE and western blot analysis. Dkk-1 disrupts the canonical pathway by binding to the LRP5/6 coreceptors, preventing interaction with Wnt and facilitating degradation of β -catenin, which ultimately results in osteolytic breakdown of bone tissue.

Wnt signaling dysregulation has been implicated in cancer of the colon and stomach⁴. Gonzalez-Sancho reported that the Wnt/ β -catenin pathway is downregulated by the induction of Dkk-1 expression, a mechanism that is lost in colon cancer, suggesting that Dkk-1 acts as a tumor suppressor gene in this cancer type⁵. Dkk-1 proteins are expressed during log phase of the cell cycle, suggesting that dysregulated Dkk-1 could be responsible for the growth of some malignancies. The Dkk-1 gene has also been determined to be a target of p53, as Cappuccio suggests it is a necessary requirement for the execution of cell death⁶. Tian reports that elevated Dkk-1 levels in bone marrow, plasma, and peripheral blood from patients with multiple myeloma correlated with the gene-expression patterns of Dkk-1, and were associated with the presence of focal bone lesions⁷. A recent study reported serum levels of Dkk-1 in patients with multiple myeloma before autologous stem cell transplantation⁸. The range of Dkk-1 serum levels in this group of patients was 6.4-234.8 ng/ml with a mean of

63.6 ng/ml and a standard deviation of 77.3. Hall et al. suggest that Dkk-1 may act like a switch on prostate cancer cell activity, causing cells to change from osteoblastic to a highly osteolytic cell line⁹.

PRINCIPLE

1. Samples and standards are added to wells coated with a monoclonal antibody specific for Dkk-1. The plate is then incubated.
2. The plate is washed, leaving only bound Dkk-1 on the plate. A yellow solution of biotinylated polyclonal antibody to Dkk-1 is then added. This binds the Dkk-1 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 Dkk-1 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 Dkk-1 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 antigen.



Activity of conjugate is affected by concentrations of chelators > 10mM (such as EDTA and EGTA).



Protect substrate from prolonged exposure to light.



Stop solution is caustic. Keep tightly capped.

MATERIALS SUPPLIED

- 1. Assay Buffer 10 Concentrate (10X)**
27ml, Catalog No. 80-0648
Tris buffered saline containing detergents.
- 2. human Dkk-1 Standard**
250pg/vial, Catalog No. 80-1465
Two vials containing 250 pg of recombinant human Dkk-1.
- 3. Dkk-1 Clear Microtiter Plate**
One Plate of 96 Wells, Catalog No. 80-1603
A plate of break-apart strips coated with a mouse monoclonal antibody specific to Dkk-1.
- 4. Wash Buffer Concentrate**
100ml, Catalog No. 80-1287
Tris buffered saline containing detergents.
- 5. Dkk-1 EIA Antibody**
10ml, Catalog No. 80-1461
- 6. human Dkk-1 EIA Conjugate**
10ml, Catalog No. 80-1463
A blue solution of streptavidin conjugated to horseradish peroxidase.
- 7. TMB Substrate**
10ml, Catalog No. 80-0350
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide.
- 8. Stop Solution 2**
10ml, Catalog No. 80-0377
A 1N solution of hydrochloric acid in water.
- 9. Dkk-1 Assay Layout Sheet**
1 each, Catalog No. 30-0230
- 10. Plate Sealer**
3 each, Catalog No. 30-0012
A yellow solution of biotinylated polyclonal antibody to Dkk-1.



Reagents require separate storage conditions.

STORAGE

All components of this kit, except the standard, are stable at 4°C until the kit's expiration date. **The standard should be stored at or below -20°C upon receipt.**

OTHER MATERIALS NEEDED

1. Deionized or distilled water.
2. Phenylmethylsulfonyl fluoride (PMSF), Sigma #P7626 or equivalent.
3. Protease Inhibitor Cocktail (PIC), Sigma #P8340 or equivalent.
4. Precision pipets for volumes between 5µl and 1,000µl.
5. Repeater pipet for dispensing 100µl.
6. Disposable beakers for diluting buffer concentrates.
7. Graduated cylinders.
8. A microplate shaker.
9. Lint-free paper for blotting.
10. Microplate reader capable of reading at 450 nm.
11. Graph paper for plotting the standard curve.



Bring all reagents to room temperature for at least 30 minutes prior to opening.



If inhibitors other than those recommended are used, the end user is responsible for assay validation. In some cases, certain protease inhibitor cocktails may cause performance differences.



Plastic tubes must be used for standard preparation.

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. Assay Buffer

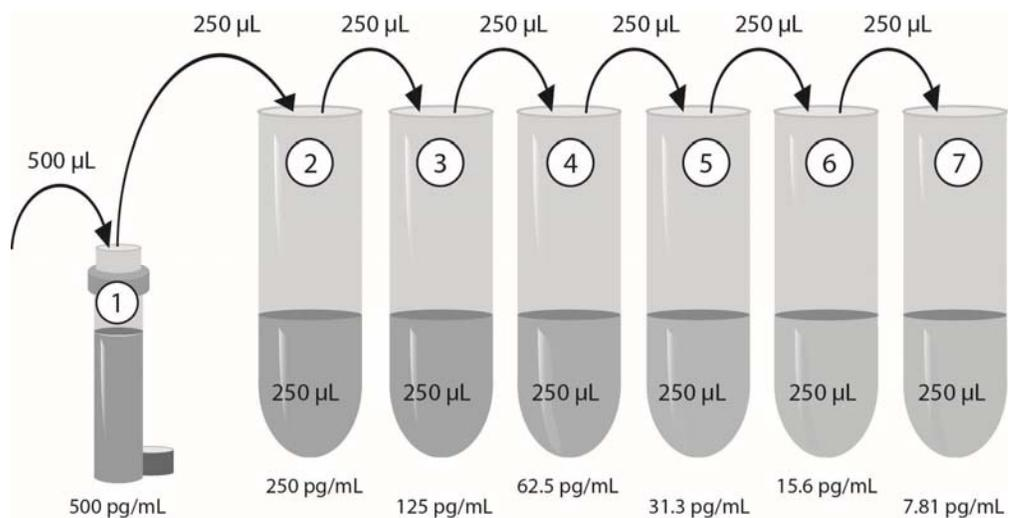
Prepare the assay buffer by diluting 5ml of the supplied Assay Buffer 10 Concentrate with 45ml of deionized water. This can be stored at 4°C until the kit expiration, or for 3 months, whichever is earlier.

3. PIC and PMSF Addition

Immediately prior to use, PIC and PMSF must be added to the assay buffer. If using Sigma Protease Inhibitor Cocktail #P8340, add 0.5 $\mu\text{L}/\text{ml}$ PIC, or equivalent concentration according to alternate vendor's specification sheet. Add PMSF, such as Sigma #P7626, to a final concentration of 1 mM.

Inhibitors must be freshly added to the assay buffer to ensure optimal integrity of the standards and samples. Each inhibitor treated buffer should incubate for 5-10 minutes at room temperature before it is used. Buffers treated with inhibitors should be used within 1 hour of preparation.

4. human Dkk-1 Standards



Add 500 μl of assay buffer plus inhibitors to the lyophilized human Dkk-1 standard vial and vortex. Wait 5 minutes and vortex again prior to use. Label the vial Standard #1. Label six 12 x 75 mm polypropylene tubes #2 through #7. Pipet 250 μl of the assay buffer plus inhibitors into tubes #2 through #7. Add 250 μl of reconstituted Standard #1 to tube #2 and vortex thoroughly. Add 250 μl of tube #2 to tube #3 and vortex thoroughly. Continue this for tubes #4 through #7.

Diluted standards should be used within 30 minutes of preparation.

The concentration of Dkk-1 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.



Samples must be stored at or below -20°C to avoid loss of bioactive analyte. Avoid repeated freeze/thaw cycles.



Add PIC/PMSF to buffers prior to preparing samples.

SAMPLE HANDLING

Human culture fluids, serum, EDTA plasma, and heparin plasma are suitable for use in the assay. Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens.

Dkk-1 concentrations were measured in serum, EDTA plasma, and culture media. Dkk-1 was spiked into the samples and diluted with assay buffer plus inhibitors, then assayed in the kit.

Sample	Recommended Dilution	% Recovery
Human plasma EDTA	≥1:8	109%
MEM tissue culture media	≥1:3	116%
Human serum	≥1:8	109%

Culture media supplemented with 10% fetal bovine serum (FBS) requires a minimum dilution of 1:3 in the assay buffer plus inhibitors. This kit also recognizes bovine Dkk-1.* It is recommended that users evaluate each lot of FBS used for the levels of endogenous bovine Dkk-1 present. The user has the option to either dilute the culture fluids to eliminate the endogenous bovine Dkk-1 or subtract the concentration of bovine Dkk-1 contributing to the total Dkk-1 detected.

*Bovine Dkk-1 has 90% sequence identity to human Dkk-1. Cross-reactivity to bovine Dkk-1 is observed.

Sample Values

The following human samples were diluted to the minimum recommended dilution and tested in the assay for presence of Dkk-1.

Sample Type	Pathology	# Samples	Concentration (pg/ml)		
			Average	Median	Range
Serum	None	12	1130	891	295-2351
Plasma	None	10	1442	1348	799-2229
Serum	Colon Cancer	10	6354	6513	3137-9791
Plasma	Colon Cancer	10	2633	2354	1026-4604

ASSAY PROCEDURE

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 100µl of the assay buffer plus inhibitors into the S0 (0pg/ml standard) wells.
2. Pipet 100µl of Standards #1 through #7 to the bottom of the appropriate wells.
3. Pipet 100µl of the samples to the bottom of the appropriate wells.
4. Seal the plate. Incubate for 1 hour on a plate shaker (~500rpm) at room temperature.
5. Empty the contents of the wells and wash by adding 400µl 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 100µl of yellow antibody into each well except the blank.
7. Seal the plate. Incubate for 1 hour on a plate shaker (~500rpm) at room temperature.
8. Wash as above (Step 5).
9. Add 100µl of blue conjugate to each well except the blank.
10. Seal the plate. Incubate for 30 minutes on a plate shaker (~500rpm) at room temperature.
11. Wash as above (Step 5).
12. Pipet 100µl of substrate solution into each well.
13. Incubate for 30 minutes on a plate shaker (~500rpm) at room temperature.
14. Pipet 100µl 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.



Bring all reagents 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 reagent to the side of the wells to avoid possible contamination.



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



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 Dkk-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.

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

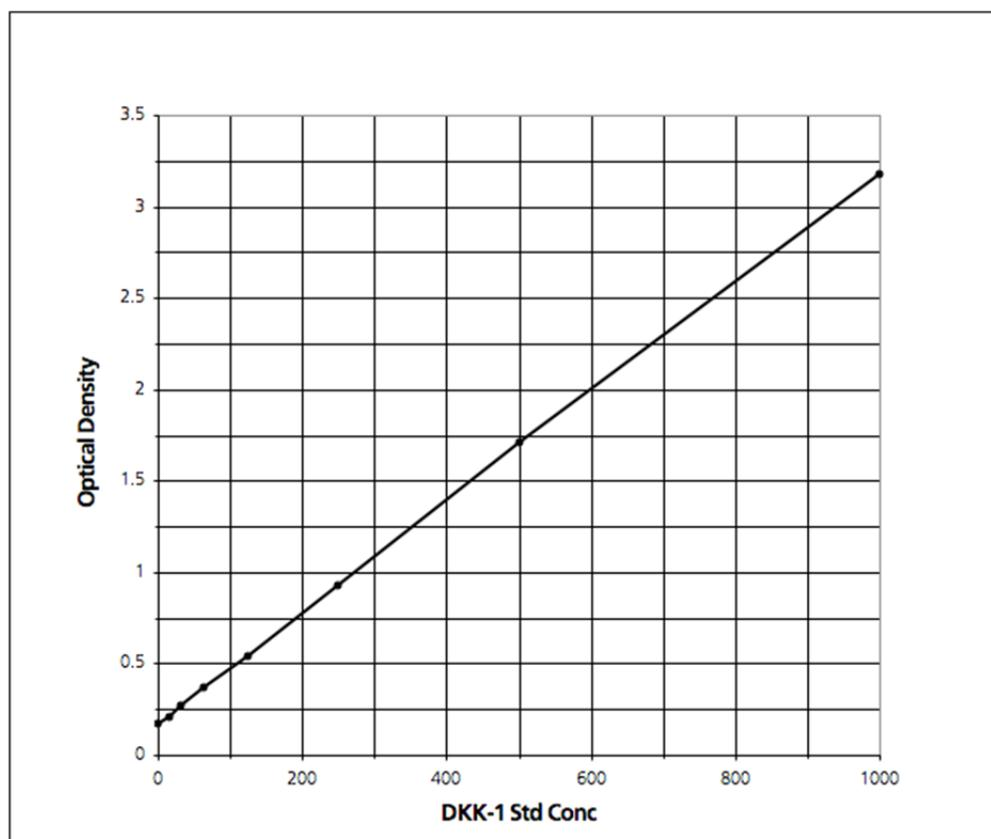
2. Using linear graph paper, plot the average Net OD for each standard versus Dkk-1 concentration in each standard. Approximate a straight line through the points. The concentration of the unknowns can be determined by interpolation.

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

TYPICAL RESULTS

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

Sample	Net OD	Dkk-1 (pg/ml)
S0	0.168	
S1	3.178	500
S2	1.706	250
S3	0.930	125
S4	0.536	62.5
S5	0.366	31.3
S6	0.268	15.6
S7	0.207	7.81
Unknown 1	2.275	345
Unknown 2	0.373	34.4



PERFORMANCE CHARACTERISTICS

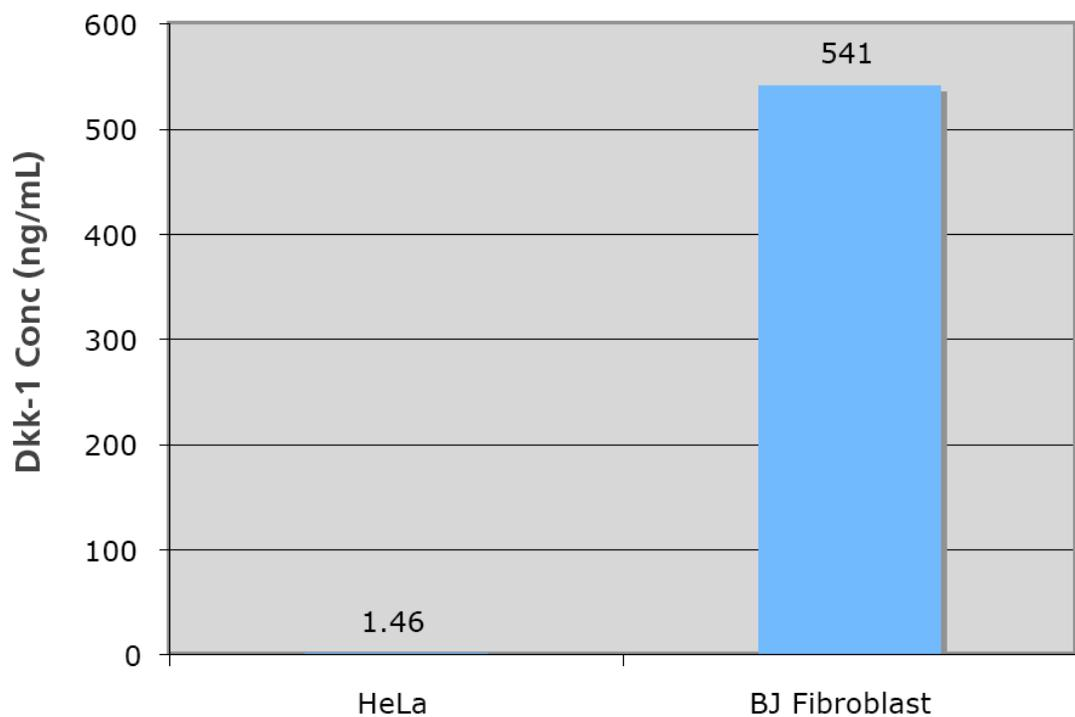
Specificity

The cross reactivities for a number of related compounds were determined by diluting cross reactants in the assay buffer at a concentration of 50,000pg/ml. These samples were then measured in the assay.

Compound	Cross Reactivity
human Dkk-1	100%
Dkk-4	<0.02%
GSK-3 β	<0.02%
β -catenin	<0.02%
PIN 1	<0.02%

Cell Line Experiment

Dkk-1 is expected to be secreted from fibroblast and fibroblast-like cells. HeLa cells are known to be negative for Dkk-1. BJ fibroblast supernatants were serially diluted into assay buffer beginning at a 1:200 dilution. HeLa cell supernatants were serially diluted into assay buffer beginning at a 1:2 dilution. Supernatants came from comparable cell numbers.



Sensitivity

Sensitivity was calculated as the ratio of the mean OD plus 2 standard deviations of twenty-four replicates of the 0pg/ml standard to the mean of twenty-four replicates of the lowest standard, multiplied by the concentration of that standard (7.81pg/ml). This value was determined to be 0.979pg/ml.

Linearity

A buffer sample containing Dkk-1, at a concentration of 400pg/ml, was serially diluted 1:2 in the assay buffer and measured in the assay. The results are shown in the table below.

Dilution	Expected (pg/ml)	Observed (pg/ml)	Recovery (%)
Neat	-----	400.44pg/ml	-----
1:2	200.22pg/ml	218.88pg/ml	109.3%
1:4	100.11pg/ml	99.43pg/ml	99.3%
1:8	50.06pg/ml	52.34pg/ml	104.6%
1:16	25.03pg/ml	26.68pg/ml	106.6%
1:32	12.51pg/ml	13.07pg/ml	104.4%

Precision

Intra-assay precision was determined by assaying 20 replicates of three buffer controls containing Dkk-1 in a single assay.

pg/ml	%CV
296.3	7.4
86.2	6.7
35.3	3.7

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

pg/ml	%CV
329.9	7.9
90.9	12.8
30.4	13.3

REFERENCES

1. Venkatesh K, Bryant H, MacDougald O. Regulation of bone mass by Wnt signaling. *J Clin Invest.* 2006 May;166(5):1202-9.
2. Prockop D, Gregory C, Spees J. One strategy for cell and gene therapy: Harnessing the power of adult stem cells to repair tissues. *PNAS* 2003 Sep;100(supp1):11917-23.
3. John Hopkins University. OMIM-Online Mendelian Inheritance in Man. (2000). <http://www.ncbi.nlm.nih.gov/entrez/dispomim.cgi?id=605189>.
4. Byun T, Karimi M, Marsh JL, Milovanovic T, Lin F, Holcombe RF. Expression of secreted Wnt antagonists in gastrointestinal tissues: potential role in stem cell homeostasis. *J Clin Pathol.* 2005 May;58(5):515-9.
5. Gonzalez-Sancho JM, Aguilera O, Garcia JM, Pendas-Franco N, Pena C, Cal S, Garcia de Herreros A, Bonilla F, Munoz A. The Wnt antagonist DICKKOPF-1 gene is a downstream target of β -catenin/TCF and is downregulated in human colon cancer. *Oncogene.* 2005 Feb 3;24(6):1098-103.
6. Cappuccio I, Calderone A, Busceti C, Biagioni F, Pontarelli F, Bruno V, Storto M, Terstappen G, Gaviraghi G, Fornai F, Battaglia G, Melchiorri d, Zukin S, Nicoletti F, Caricasole A. Induction of Dickkopf-1, a negative modulator of the Wnt pathway, is required for the development of ischemic neuronal death. *J Neurosci.* 2005 Mar 9;25(10):2647-57.
7. Tian E, Zhan F, Walker R, Rasmussen E, Ma Y, Barlogie B, Shaughnessy JD Jr. The role of the Wnt-signaling antagonist DKK1 in the development of osteolytic lesions in multiple myeloma. *N Engl J Med.* 2003 Dec 25;349(26):2483-94.
8. Politou MC, Heath DJ, Rahemtulla A, Szydlo R, Anagnostopoulos A, Dimopoulos MA, Croucher PI, Terpos E. Serum concentrations of Dickkopf-1 protein are increased in patients with multiple myeloma and reduced after autologous stem cell transplantation. *Int J Cancer.* 2006 Oct;119(7): 1728-31.
9. Hall, CL, Bafico A, Dai J, Aaronson SA, Keller, ET. Prostate cancer cells use Wnt signaling proteins to promote growth of painful bone tumors. *Cancer Research* 2005 Sep 1;65(17):7554-60.



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com