



# **ab197746 – DKK1 (Dickkopf-1) Mouse SimpleStep ELISA<sup>®</sup> Kit**

## Instructions for Use

For the quantitative measurement of DKK1 (Dickkopf-1) in mouse serum, plasma, and cell culture supernatants.

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

## INTRODUCTION

|                  |   |
|------------------|---|
| 1. BACKGROUND    | 2 |
| 2. ASSAY SUMMARY | 3 |

## GENERAL INFORMATION

|                                     |   |
|-------------------------------------|---|
| 3. PRECAUTIONS                      | 4 |
| 4. STORAGE AND STABILITY            | 4 |
| 5. MATERIALS SUPPLIED               | 4 |
| 6. MATERIALS REQUIRED, NOT SUPPLIED | 5 |
| 7. LIMITATIONS                      | 5 |
| 8. TECHNICAL HINTS                  | 5 |

## ASSAY PREPARATION

|                          |    |
|--------------------------|----|
| 9. REAGENT PREPARATION   | 7  |
| 10. STANDARD PREPARATION | 8  |
| 11. SAMPLE PREPARATION   | 9  |
| 12. PLATE PREPARATION    | 10 |

## ASSAY PROCEDURE

|                     |    |
|---------------------|----|
| 13. ASSAY PROCEDURE | 11 |
|---------------------|----|

## DATA ANALYSIS

|                           |    |
|---------------------------|----|
| 14. CALCULATIONS          | 13 |
| 15. TYPICAL DATA          | 14 |
| 16. TYPICAL SAMPLE VALUES | 15 |
| 17. SPECIES REACTIVITY    | 19 |

## RESOURCES

|                     |    |
|---------------------|----|
| 18. TROUBLESHOOTING | 20 |
| 19. NOTES           | 21 |

## 1. BACKGROUND

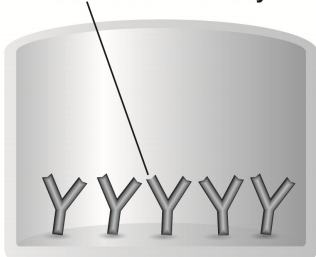
DKK1 (Dickkopf-1) *in vitro* SimpleStep ELISA™ (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of DKK1 protein in mouse serum, plasma, and cell culture supernatants.

The SimpleStep ELISA™ employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB Development Solution is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

Dickkopf-related proteins are secreted proteins with two cysteine-rich domains. They are important in early vertebrate development due to their antagonistic relationship to Wnt signaling, which is involved in embryonic development, limb development, somitogenesis, and eye formation. In adults Dickkopf-related proteins are important to bone formation and bone related diseases, as well as certain cancers and Alzheimer disease. Dickkopf-related protein 1 (DKK1) is one of two secreted inhibitors of canonical Wnt signaling. DKK1 decreases Wnt signaling by inhibiting the interaction between LRP5/6 and Wnt-frizzled complexes. Elevated DKK1 levels are associated with the presence of osteolytic bone lesions in multiple myeloma patients, and may also contribute to the osteopenia and hypercalcemia often associated with myeloma. The inhibition of DKK1 has shown the reversal of the bone deteriorating patterns associated with osteoarthritis.

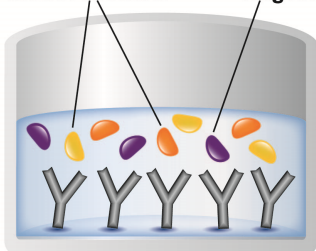
## 2. ASSAY SUMMARY

**Immobilization Antibody**



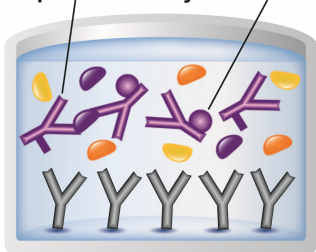
Remove appropriate number of antibody coated well strips. Equilibrate all reagents to room temperature. Prepare all reagents, samples, and standards as instructed.

**Matrix Proteins      Target Analyte**



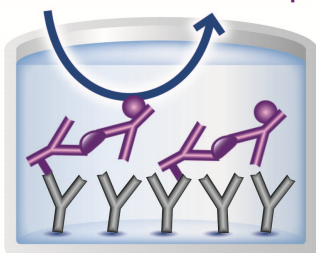
Add standard or sample to appropriate wells.

**Capture Antibody      Detector Antibody**



Add Antibody Cocktail to all wells. Incubate at room temperature.

**Substrate      Color Development**



Aspirate and wash each well. Add TMB Development Solution to each well and incubate. Add Stop Solution at a defined endpoint.

Alternatively, record color development kinetically after TMB substrate addition.

## 3. PRECAUTIONS

**Please read these instructions carefully prior to beginning the assay.**

All kit components have been formulated and quality control tested to function successfully as a kit. Modifications to the kit components or procedures may result in loss of performance.

## 4. STORAGE AND STABILITY

**Store kit at 2-8°C immediately upon receipt.**

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Reagent and Standard Preparation sections.

## 5. MATERIALS SUPPLIED

| Item   | Amount   | Storage Condition (Before Preparation) |
|--|----------|--|
| 10X Mouse DKK1 Capture Antibody                    | 600 µL   | +2-8°C                                 |
| 10X Mouse DKK1 Detector Antibody                   | 600 µL   | +2-8°C                                 |
| DKK1 Mouse Lyophilized Recombinant Protein         | 2 Vials  | +2-8°C                                 |
| Antibody Diluent 5BI                               | 6 mL     | +2-8°C                                 |
| 10X Wash Buffer PT                                 | 20 mL    | +2-8°C                                 |
| TMB Development Solution                           | 12 mL    | +2-8°C                                 |
| Stop Solution                                      | 12 mL    | +2-8°C                                 |
| Sample Diluent NS                                  | 50 mL    | +2-8°C                                 |
| Pre-Coated 96 Well Microplate (12 x 8 well strips) | 96 Wells | +2-8°C                                 |
| Plate Seal   | 1        | +2-8°C                                 |

### 6. MATERIALS REQUIRED, NOT SUPPLIED

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Microplate reader capable of measuring absorbance at 450 or 600 nm
- Method for determining protein concentration (BCA assay recommended)
- Deionized water
- PBS (1.4 mM  $\text{KH}_2\text{PO}_4$ , 8 mM  $\text{Na}_2\text{HPO}_4$ , 140 mM NaCl, 2.7 mM KCl, pH 7.4)
- Multi- and single-channel pipettes
- Tubes for standard dilution
- Plate shaker for all incubation steps
- Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors)

### 7. LIMITATIONS

- Assay kit intended for research use only. Not for use in diagnostic procedures
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted

### 8. TECHNICAL HINTS

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers
- Avoid foaming or bubbles when mixing or reconstituting components
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions

- Ensure plates are properly sealed or covered during incubation steps
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background
- As a guide, typical ranges of sample concentration for commonly used sample types are shown below in Sample Preparation (section 11)
- All samples should be mixed thoroughly and gently
- Avoid multiple freeze/thaw of samples
- Incubate ELISA plates on a plate shaker during all incubation steps
- When generating positive control samples, it is advisable to change pipette tips after each step
- **To avoid high background always add samples or standards to the well before the addition of the antibody cocktail**
- **This kit is sold based on number of tests. A ‘test’ simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions**

## 9. REAGENT PREPARATION

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. **The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.**
- Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

### 9.1 **1X Wash Buffer PT**

Prepare 1X Wash Buffer PT by diluting 10X Wash Buffer PT with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL 10X Wash Buffer PT with 45 mL deionized water. Mix thoroughly and gently.

### 9.2 **Antibody Cocktail**

Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 5BI. To make 3 mL of the Antibody Cocktail combine 300  $\mu$ L 10X Capture Antibody and 300  $\mu$ L 10X Detector Antibody with 2.4 mL Antibody Diluent 5BI. Mix thoroughly and gently.

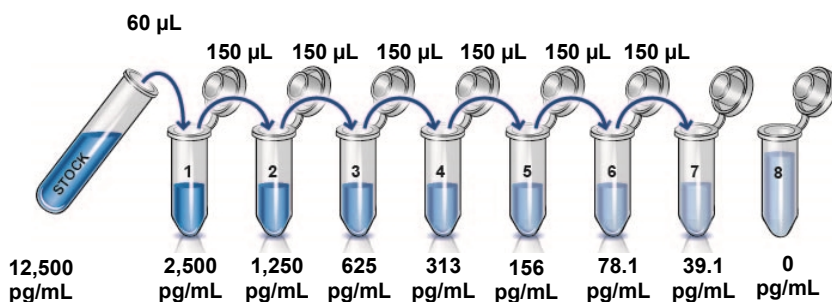


## 10. STANDARD PREPARATION

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of positive controls for every use.

The following table describes the preparation of a standard curve for duplicate measurements (recommended).

- 10.1 **IMPORTANT:** If the protein standard vial has a volume identified on the label, reconstitute the DKK1 standard by adding that volume of Sample Diluent NS indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the DKK1 standard by adding 1 mL Sample Diluent NS. Hold at room temperature for 10 minutes and mix gently. This is the 12,500 pg/mL **Stock Standard Solution**.
- 10.2 Label eight tubes, Standards 1–8.
- 10.3 Add 240  $\mu$ L Sample Diluent NS into tube number 1 and 150  $\mu$ L of Sample Diluent NS into numbers 2-8.
- 10.4 Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



## 11. SAMPLE PREPARATION

| TYPICAL SAMPLE DYNAMIC RANGE |              |
|------------------------------|--------------|
| Sample Type                  | Range (%)    |
| Mouse Serum                  | 6.25 - 0.391 |
| Mouse Plasma - Citrate       | 50 - 0.391   |

### 11.1 Plasma

Collect plasma using citrate. Centrifuge samples at 2,000 x g for 10 minutes. Dilute samples into Sample Diluent NS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

### 11.2 Serum

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Dilute samples into Sample Diluent NS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

### 11.3 Cell Culture Supernatants

Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants and assay. Store samples at -20°C or below. Avoid repeated freeze-thaw cycles.

## 12. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents
- Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C
- For each assay performed, a minimum of two wells must be used as the zero control
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates)
- Differences in well absorbance or “edge effects” have not been observed with this assay

## 13. ASSAY PROCEDURE

- **Equilibrate all materials and prepared reagents to room temperature prior to use.**
- **It is recommended to assay all standards, controls and samples in duplicate.**

- 13.1 Prepare all reagents, working standards, and samples as directed in the previous sections.
- 13.2 Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
- 13.3 Add 50 µL of all sample or standard to appropriate wells.
- 13.4 Add 50 µL of the Antibody Cocktail to each well.
- 13.5 Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
- 13.6 Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and blot it against clean paper towels to remove excess liquid.
- 13.7 Add 100 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm.

*Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.*

*Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.*

- 13.8 Add 100 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.

*Alternative to 13.7 – 13.8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:*

|             |                        |
|-------------|------------------------|
| Mode:       | Kinetic                |
| Wavelength: | 600 nm                 |
| Time:       | up to 20 min           |
| Interval:   | 20 sec - 1 min         |
| Shaking:    | Shake between readings |

*Note that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100  $\mu$ L Stop Solution to each well and recording the OD at 450 nm.*

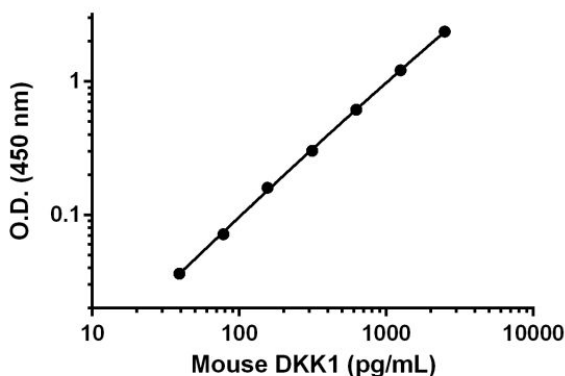
13.9 Analyze the data as described below.

### 14. CALCULATIONS

Subtract average zero standard from all readings. Average the duplicate readings of the positive control dilutions and plot against their concentrations. Draw the best smooth curve through these points to construct a standard curve. Most plate reader software or graphing software can plot these values and curve fit. A four parameter algorithm (4PL) usually provides the best fit, though other equations can be examined to see which provides the most accurate (e.g. linear, semi-log, log/log, 4 parameter logistic). Interpolate protein concentrations for unknown samples from the standard curve plotted. Samples producing signals greater than that of the highest standard should be further diluted and reanalyzed, then multiplying the concentration found by the appropriate dilution factor.

## 15. TYPICAL DATA

**TYPICAL STANDARD CURVE** – Data provided for **demonstration purposes only**. A new standard curve must be generated for each assay performed.



| Standard Curve Measurements |             |       |              |
|-----------------------------|-------------|-------|--------------|
| Conc.<br>(pg/mL)            | O.D. 450 nm |       | Mean<br>O.D. |
|                             | 1           | 2     |              |
| 0                           | 0.069       | 0.070 | 0.069        |
| 39.1                        | 0.087       | 0.090 | 0.089        |
| 78.1                        | 0.123       | 0.126 | 0.124        |
| 156                         | 0.211       | 0.215 | 0.213        |
| 313                         | 0.352       | 0.358 | 0.355        |
| 625                         | 0.655       | 0.676 | 0.666        |
| 1,250                       | 1.266       | 1.257 | 1.261        |
| 2,500                       | 2.341       | 2.473 | 2.407        |

**Figure 1.** Example of DKK1 standard curve. The DKK1 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean  $\pm$  SD) are graphed.

## 16. TYPICAL SAMPLE VALUES

### **SENSITIVITY –**

The calculated minimal detectable dose (MDD) is 6.66 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=8) and adding 2 standard deviations then extrapolating the corresponding concentrations.

### **RECOVERY –**

Three concentrations of DKK1 were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

| Sample Type            | Average % Recovery | Range (%) |
|------------------------|--------------------|-----------|
| Mouse Serum            | 114                | 109 - 120 |
| Mouse Plasma - Citrate | 112                | 99 - 121  |
| Cell Culture Media     | 99                 | 93 - 105  |



## LINEARITY OF DILUTION –

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Native DKK1 was measured in mouse serum and plasma samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

DKK1 protein was spiked into cell culture media and diluted in a 2-fold dilution series in Sample Diluent NS.

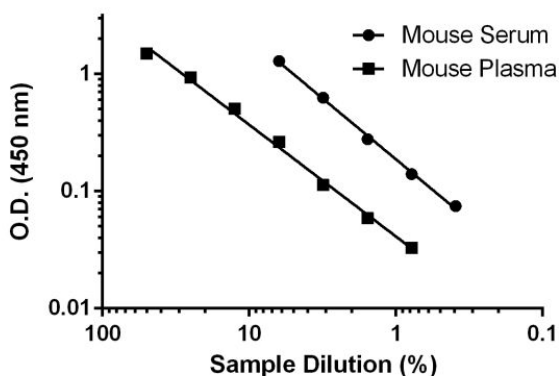
| Dilution Factor | Interpolated value | 6.25% Mouse Serum | 50% Mouse Plasma (Citrate) | 100% Cell Culture Media |
|-----------------|--------------------|-------------------|----------------------------|-------------------------|
| Undiluted       | pg/mL              | 1,357             | 1,604                      | 2,961                   |
|                 | % Expected value   | 100               | 100                        | 100                     |
| 2               | pg/mL              | 622               | 952                        | 1,394                   |
|                 | % Expected value   | 92                | 119                        | 94                      |
| 4               | pg/mL              | 266               | 495                        | 614                     |
|                 | % Expected value   | 78                | 124                        | 83                      |
| 8               | pg/mL              | 132               | 243                        | 301                     |
|                 | % Expected value   | 78                | 121                        | 81                      |
| 16              | pg/mL              | 69.6              | 106                        | 152                     |
|                 | % Expected value   | 82                | 106                        | 82                      |
| 32              | pg/mL              | NL                | 54.6                       | 76.8                    |
|                 | % Expected value   | NL                | 109                        | 83                      |
| 64              | pg/mL              | NL                | 29.2                       | 40.6                    |
|                 | % Expected value   | NL                | 117                        | 88                      |

NL – Non-Linear

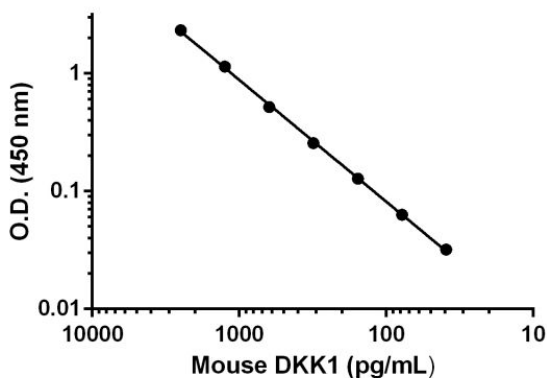
## PRECISION –

Mean coefficient of variations of interpolated values from 3 concentrations of DKK1 within the working range of the assay.

|        | Intra-Assay | Inter-Assay |
|--------|-------------|-------------|
| n=     | 8           | 3           |
| CV (%) | 5.20        | 5.26        |



**Figure 2.** Titration of mouse Serum and mouse plasma within the working range of the assay. Background-subtracted data values (mean  $\pm$  SD,  $n = 2$ ) are graphed.



**Figure 3.** Titration of protein spiked cell culture media within the working range of the assay. Background-subtracted data values (mean  $\pm$  SD,  $n = 2$ ) are graphed.

### 17. SPECIES REACTIVITY

This kit recognizes mouse DKK1 protein.

Other species reactivity was determined by measuring 6.25% (1:16) serum samples of various species, interpolating the protein concentrations from the Mouse standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in mouse serum assayed at the same dilution.

Reactivity < 3% was determined for the following species:

Human

Rat

Hamster

Guinea Pig

Rabbit

Dog

Goat

Pig

Cow

Please contact our Technical Support team for more information.

## 18. TROUBLESHOOTING

| Problem                                      | Cause   | Solution   |
|--|---|--|
| Difficulty pipetting lysate; viscous lysate. | Genomic DNA solubilized                         | Prepare 1X Cell Extraction Buffer PTR (without enhancer). Add enhancer to lysate after extraction.                               |
| Poor standard curve                          | Inaccurate Pipetting                            | Check pipettes   |
|  | Improper standard dilution                      | Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing                       |
| Low Signal                                   | Incubation times too brief                      | Ensure sufficient incubation times; increase to 2 or 3 hour standard/sample incubation   |
|  | Inadequate reagent volumes or improper dilution | Check pipettes and ensure correct preparation  |
|  | Incubation times with TMB too brief             | Ensure sufficient incubation time until blue color develops prior addition of Stop solution                                      |
| Large CV                                     | Plate is insufficiently washed                  | Review manual for proper wash technique. If using a plate washer, check all ports for obstructions.                              |
|  | Contaminated wash buffer                        | Prepare fresh wash buffer  |
| Low sensitivity                              | Improper storage of the ELISA kit               | Store your reconstituted standards at -80°C, all other assay components 4°C. Keep TMB Development Solution protected from light. |

### 19. NOTES



## Technical Support

Copyright © 2022 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)