



# **ab202401 – Survivin (Birc5) Mouse SimpleStep ELISA<sup>®</sup> Kit**

## Instructions for Use

For the quantitative measurement of Survivin (Birc5) in mouse serum, plasma, cell culture supernatant, cell and tissue extracts.

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

# Table of Contents

---

## INTRODUCTION

1. BACKGROUND	2
2. ASSAY SUMMARY	4

## GENERAL INFORMATION

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

## ASSAY PREPARATION

9. REAGENT PREPARATION	8
10. STANDARD PREPARATION	9
11. SAMPLE PREPARATION	12
12. PLATE PREPARATION	15

## ASSAY PROCEDURE

13. ASSAY PROCEDURE	16
---------------------	----

## DATA ANALYSIS

14. CALCULATIONS	18
15. TYPICAL DATA	19
16. TYPICAL SAMPLE VALUES	23
17. ASSAY SPECIFICITY	27
18. SPECIES REACTIVITY	28

## RESOURCES

19. TROUBLESHOOTING	29
20. NOTES	30

## 1. **BACKGROUND**

Survivin *in vitro* SimpleStep ELISA® kit is designed for the quantitative measurement of Survivin protein in mouse serum, plasma, cell culture supernatant, cell and tissue extracts.

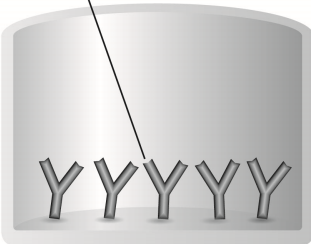
The SimpleStep ELISA® employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating 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.

Survivin is a multitasking protein that has dual roles in promoting cell proliferation and preventing apoptosis. Survivin is a component of a chromosome passage protein complex (CPC) which is essential for chromosome alignment and segregation during mitosis and cytokinesis. Survivin acts as an important regulator of the localization of this complex; directs CPC movement to different locations from the inner centromere during prometaphase to midbody during cytokinesis and it participates in the organization of the center spindle by associating with polymerized microtubules. Survivin complex with RAN plays a role in mitotic spindle formation by serving as a physical scaffold to help deliver the RAN effector molecule TPX2 to microtubules. Survivin may counteract a default induction of apoptosis in G2/M phase. The acetylated form of Survivin represses STAT3 transactivation of target gene promoters. Survivin may play a role in

neoplasia. Survivin may be an inhibitor of CASP3 and CASP7. Isoform 2 and isoform 3 of Survivin do not appear to play vital roles in mitosis. Isoform 3 shows a marked reduction in its anti-apoptotic effects when compared with the displayed wild-type isoform.

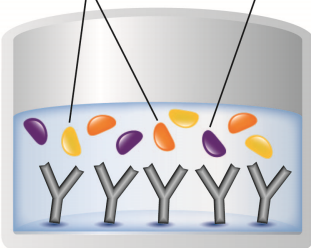
## 2. ASSAY SUMMARY

Immobilization Antibody



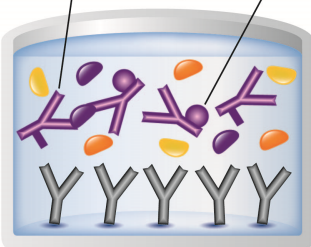
Remove appropriate number of antibody coated well strips. Equilibrate all reagents to room temperature. Prepare all reagents, samples, and standards as instructed.

Matrix Proteins Target Analyte



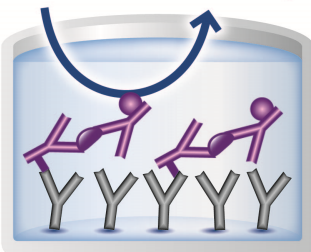
Add standard or sample to appropriate wells.

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 Survivin Capture Antibody	600 µL	+2-8°C
10X Mouse Survivin Detector Antibody	600 µL	+2-8°C
Mouse Survivin Lyophilized Recombinant Protein	2 Vials	+2-8°C
Antibody Diluent 5BC	6 mL	+2-8°C
10X Wash Buffer PT	20 mL	+2-8°C
5X Cell Extraction Buffer PTR	10 mL	+2-8°C
50X Cell Extraction Enhancer Solution	1 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
Sample Diluent NBP	20 mL	+2-8°C
Sample Diluent 25BP	20 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 KH<sub>2</sub>PO<sub>4</sub>, 8 mM Na<sub>2</sub>HPO<sub>4</sub>, 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.
- 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.
- The provided 50X Cell Extraction Enhancer Solution may precipitate when stored at + 4°C. To dissolve, warm briefly at + 37°C and mix gently. The 50X Cell Extraction Enhancer Solution can be stored at room temperature to avoid precipitation.
- **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 Cell Extraction Buffer PTR (For cell and tissue extracts only)

If required, prepare 1X Cell Extraction Buffer PTR by diluting 5X Cell Extraction Buffer PTR to 1X with deionized water. To make 10 mL 1X Cell Extraction Buffer PTR combine 8 mL deionized water and 2 mL 5X Cell Extraction Buffer PTR. Mix thoroughly and gently. If required protease inhibitors can be added.

### 9.2 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.

### 9.3 Antibody Cocktail

Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 5BC. 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 5BC. Mix thoroughly and gently.

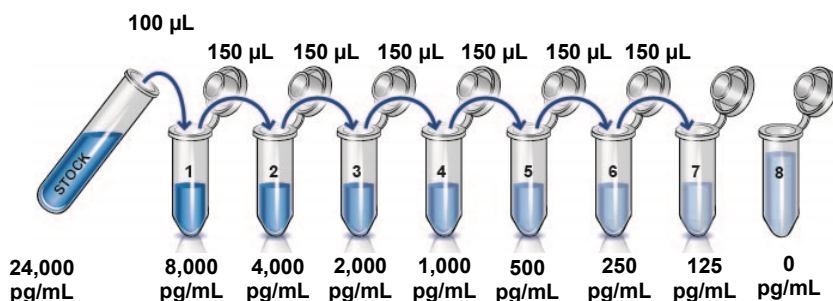
## 10. STANDARD PREPARATION

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).

**IMPORTANT:** If the protein standard vial has a volume identified on the label, reconstitute the Survivin standard by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the Survivin standard by adding 200  $\mu$ L Diluent. Hold at room temperature for 10 minutes and mix gently. This is the 24,000 pg/mL **Stock Standard** Solution.

- 10.1 For **cell and tissue extract samples** measurements follow this section to prepare the standard.
  - 10.1.1 Reconstitute the Survivin standard by adding 1X Cell Extraction Buffer PTR.
  - 10.1.2 Label eight tubes, Standards 1–8.
  - 10.1.3 Add 200  $\mu$ L 1X Cell Extraction Buffer PTR into tube number 1 and 150  $\mu$ L of 1X Cell Extraction Buffer PTR into numbers 2-8.
  - 10.1.4 Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



## ASSAY PREPARATION

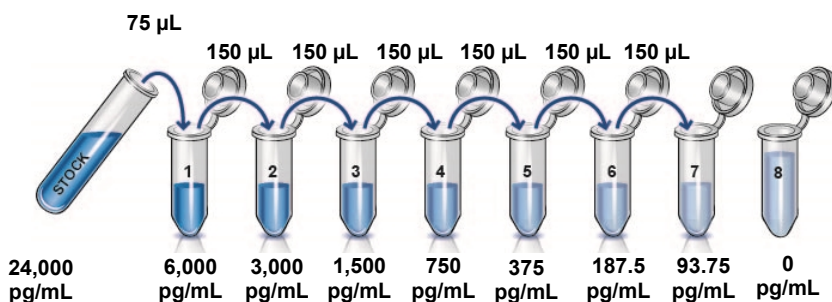
10.2 For **cell culture supernatant samples** measurements follow this section to prepare the standard

10.2.1 Reconstitute the Survivin standard by adding water.

10.2.2 Label eight tubes, Standards 1– 8.

10.2.3 Add 225  $\mu\text{L}$  Sample Diluent NS into tube number 1 and 150  $\mu\text{L}$  of Sample Diluent NS into numbers 2-8.

10.2.4 Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



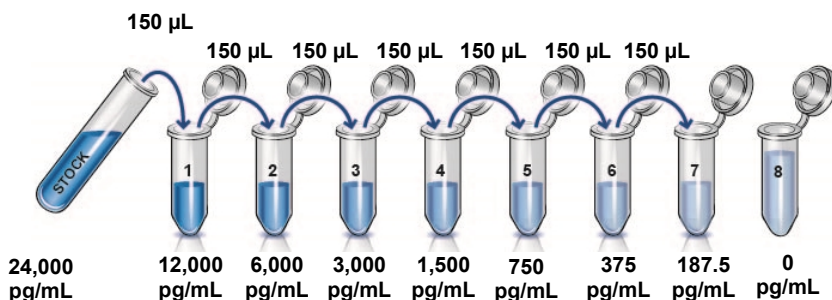
10.3 For **serum and plasma samples** measurements follow this section to prepare the standard.

10.3.1 For **serum samples** measurements, reconstitute the Survivin standard by adding Sample Diluent 25BP.

For **plasma samples** measurements, reconstitute the Survivin standard by adding Sample Diluent NBP.

10.3.2 Label eight tubes, Standards 1– 8 and add 150  $\mu$ L of appropriate diluent (see Step 10.3) into each tube.

10.3.3 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
NIH/3T3 cell extract	16 – 1000 µg/mL
Mouse Serum	Neat
Mouse Plasma - Citrate	Neat
Mouse Plasma - Heparin	Neat

### 11.1 Plasma

Collect plasma using citrate or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Plasma samples can be assayed without dilution. If needed, dilute plasma samples in normal mouse plasma (not provided) 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. Serum samples can be assayed without dilution. If needed, dilute serum samples in normal mouse serum (not provided) 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. Dilute supernatant samples at least 2 X into Sample Diluent NS and assay. Store samples at -20°C or below. Avoid repeated freeze-thaw cycles.

### 11.4 Preparation of extracts from cell pellets

- 11.4.1 Collect non-adherent cells by centrifugation or scrape to collect adherent cells from the culture flask. Typical centrifugation conditions for cells are 500 x g for 5 minutes at 4°C.
- 11.4.2 Rinse cells twice with PBS.
- 11.4.3 Solubilize pellet at  $2 \times 10^7$  cell/mL in chilled 1X Cell Extraction Buffer PTR.
- 11.4.4 Incubate on ice for 20 minutes.
- 11.4.5 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.4.6 Transfer the supernatants into clean tubes and discard the pellets.
- 11.4.7 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.4.8 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

### 11.5 Preparation of extracts from adherent cells by direct lysis (alternative protocol)

- 11.5.1 Remove growth media and rinse adherent cells 2 times in PBS.
- 11.5.2 Solubilize the cells by addition of chilled 1X Cell Extraction Buffer PTR directly to the plate (use 750  $\mu$ L - 1.5 mL 1X Cell Extraction Buffer PTR per confluent 15 cm diameter plate).
- 11.5.3 Scrape the cells into a microfuge tube and incubate the lysate on ice for 15 minutes.
- 11.5.4 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.5.5 Transfer the supernatants into clean tubes and discard the pellets.
- 11.5.6 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.

- 11.5.7 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

### 11.6 Preparation of extracts from tissue homogenates

- 11.6.1 Tissue lysates are typically prepared by homogenization of tissue that is first minced and thoroughly rinsed in PBS to remove blood (dounce homogenizer recommended).
- 11.6.2 Homogenize 100 to 200 mg of wet tissue in 500  $\mu$ L – 1 mL of chilled 1X Cell Extraction Buffer PTR. For lower amounts of tissue adjust volumes accordingly.
- 11.6.3 Incubate on ice for 20 minutes.
- 11.6.4 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.6.5 Transfer the supernatants into clean tubes and discard the pellets.
- 11.6.6 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.6.7 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

## 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.*

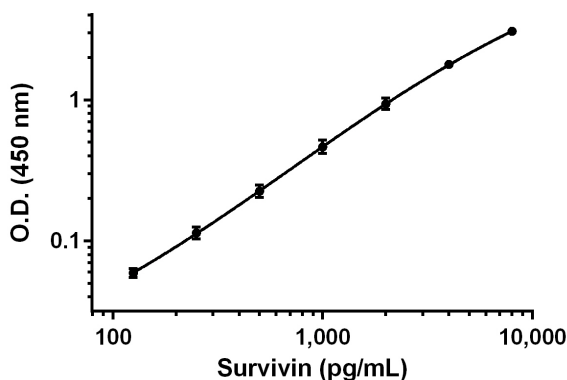
13.9 Analyze the data as described below.

### 14. CALCULATIONS

Subtract average zero standard absorbance measurement from all readings. Average absorbance measurements of the duplicate wells of the standards 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 determine 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.

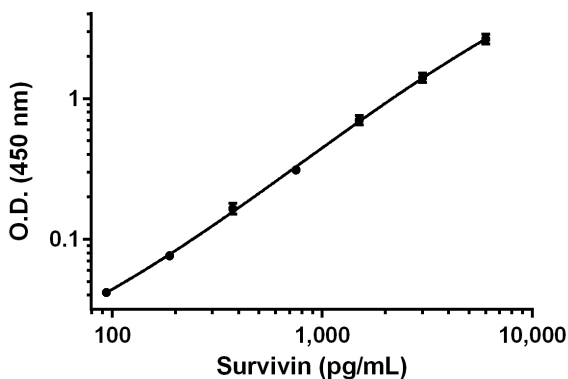
## 15. TYPICAL DATA

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



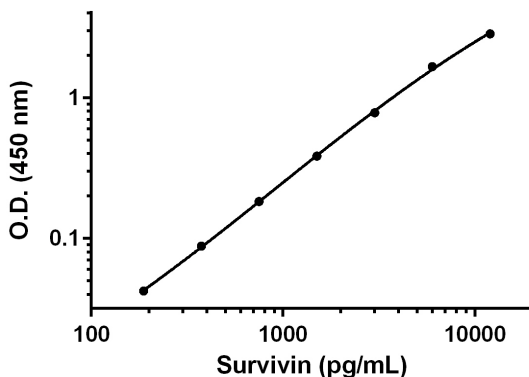
Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.050	0.049	0.050
125	0.112	0.106	0.110
250	0.172	0.156	0.164
500	0.292	0.259	0.276
1,000	0.551	0.479	0.515
2,000	1.049	0.925	0.988
4,000	1.925	1.745	1.836
8,000	3.268	2.958	3.113

**Figure 1.** Example of Survivin standard curve prepared in 1X Cell Extraction Buffer PTR. The Survivin 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.



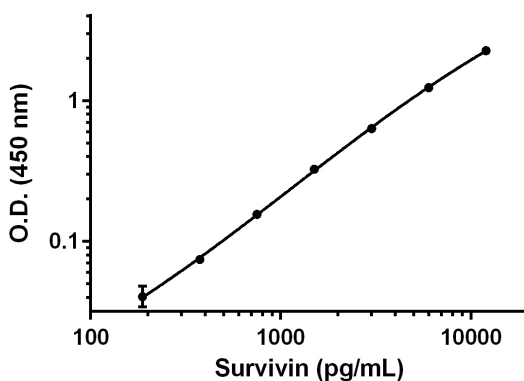
Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.051	0.053	0.052
93.75	0.093	0.095	0.094
187.5	0.127	0.130	0.129
375	0.207	0.228	0.218
750	0.369	0.358	0.364
1,500	0.717	0.794	0.756
3,000	1.375	1.532	1.454
6,000	2.550	2.860	2.706

**Figure 2.** Example of Survivin standard curve prepared in Sample Diluent NS. The Survivin 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.



Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.126	0.140	0.134
187.5	0.175	0.176	0.176
375	0.221	0.221	0.222
750	0.317	0.314	0.316
1,500	0.528	0.508	0.518
3,000	0.892	0.938	0.916
6,000	1.780	1.816	1.799
12,000	2.977	2.985	2.982

**Figure 3.** Example of Survivin standard curve prepared in Sample Diluent 25BP. The Survivin standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.085	0.077	0.082
187.5	0.127	0.117	0.123
375	0.155	0.156	0.156
750	0.238	0.236	0.237
1,500	0.413	0.402	0.408
3,000	0.717	0.715	0.716
6,000	1.336	1.305	1.321
12,000	2.367	2.352	2.360

**Figure 4.** Example of Survivin standard curve for prepared in Sample Diluent NBP. The Survivin standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

## 16. TYPICAL SAMPLE VALUES

### SENSITIVITY –

The calculated minimal detectable dose (MDD) is determined by calculating the mean of zero standard replicates and adding 2 standard deviations then extrapolating the corresponding concentrations. The MDD is dependent on the Sample Diluent buffer used:

Sample Diluent Buffer	n=	Minimal Detectable Dose
Sample Diluent NS	15	9 pg/mL
1X Cell Extraction Buffer PTR	27	25 pg/mL
Sample Diluent 25BP	11	47 pg/mL
Sample Diluent NBP	19	9 pg/mL

### RECOVERY –

For **cell and tissue extract samples** measurements, three concentrations of mouse recombinant Survivin 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 (%)
10% Human Serum	177	172.5 – 181.7
50% Cell Culture Media	133.3	122.5 – 149.8

For **cell culture supernatant, serum and plasma samples** measurements, three concentrations of mouse recombinant Survivin 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 (%)
50% Cell Culture Media	138.6	134.9 – 142.4
100% Mouse Serum	112.3	109.8 – 113.5
100% Mouse Plasma - Citrate	90	85.4 – 98.3
100% Mouse Plasma - Heparin	112.3	105.9 – 121.4

## 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 Survivin was measured in the following biological samples in a 2-fold dilution series. Sample dilutions were made in 1X Cell Extraction Buffer PTR.

<b>Dilution Factor</b>	<b>Interpolated value</b>	<b>1000 µg/mL NIH/3T3 Extract</b>
Undiluted	pg/mL	11,310.36
	<b>% Expected value</b>	<b>100</b>
2	pg/mL	4,674.96
	<b>% Expected value</b>	<b>82.7</b>
4	pg/mL	2,307.37
	<b>% Expected value</b>	<b>81.6</b>
8	pg/mL	1,239.27
	<b>% Expected value</b>	<b>87.7</b>
16	pg/mL	704.51
	<b>% Expected value</b>	<b>99.7</b>
32	pg/mL	351.41
	<b>% Expected value</b>	<b>99.4</b>

Recombinant mouse Survivin was spiked into the following biological samples and diluted in a 2-fold dilution series.

For cell culture supernatant samples measurements, Sample Diluent NS was used as diluent.

For serum samples measurements, normal mouse serum was used as diluent. For plasma (citrate) samples measurements, normal mouse plasma (citrate) was used as diluent.

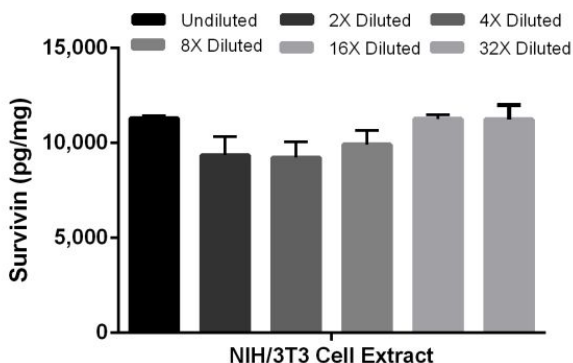
For plasma (heparin) samples measurements, normal mouse plasma (heparin) was used as diluent.

Dilution Factor	Interpolated value	50% Cell Culture Media	100% Mouse Serum	100% Mouse Plasma (Citrate)	100% Mouse Plasma (Heparin)
Undiluted	pg/mL	2,705.8	8,274.2	13,228.3	14,210
	% Expected value	100	100	100	100
2	pg/mL	1,226.5	3,594.7	6,225.4	6,861
	% Expected value	90.7	86.9	94.1	96.6
4	pg/mL	543.7	1,841.3	2,836.9	3,412.5
	% Expected value	80.4	89.0	85.8	96.1
8	pg/mL	273.4	964.8	1,532.5	1,828.2
	% Expected value	80.8	93.3	92.7	102.9
16	pg/mL	144.9	501	696	863.4
	% Expected value	85.7	96.9	84.2	97.2
32	pg/mL	74.1	269.1	471.2	553.4
	% Expected value	87.6	104.1	114	124.6

## PRECISION –

Mean coefficient of variations of interpolated values from 3 concentrations of NIH/3T3 cell extract samples within the working range of the assay.

	Intra-Assay	Inter-Assay
n=	5	3
CV (%)	3.1	13.1



**Figure 5.** Interpolated concentrations of Survivin in NIH/3T3 cell extract samples. The concentrations of Survivin were measured in duplicates, interpolated from the Survivin standard curve and corrected for sample dilution. Note that 1X Diluted extract samples were at 1 mg/mL. The interpolated, dilution factor-corrected values are plotted (mean  $\pm$  SD,  $n=2$ ).

Serum and Plasma. Neat pooled female mouse serum and plasma samples were measured in duplicates for Survivin concentrations using this kit. Means of interpolated sample values are shown in pg of Survivin per mL of serum/plasma in the table below.

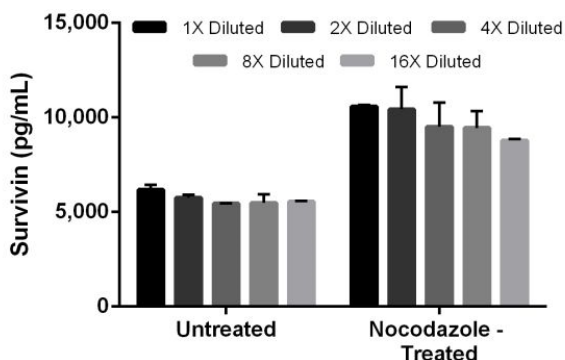
Sample Type	Mean (pg/mL)	Range (pg/mL)
Mouse Serum	190	190 - 190
Mouse Plasma - Citrate	124	112 - 135
Mouse Plasma - Heparin	163	139 - 187

## 17. ASSAY SPECIFICITY

This kit recognizes both native and recombinant mouse Survivin protein in serum, plasma (citrate and heparin), cell culture supernatant, and cell and tissue extract samples only.

Plasma (EDTA) and urine are not compatible with this assay.

Milk and saliva samples have not been tested with this kit.



**Figure 6.** Nocodazole treatment of NIH/3T3 cells stimulates expression of Survivin. NIH/3T3 cells were cultured in the absence or presence of 200 ng/mL nocodazole for 17 hours. The extracts of these cells (adjusted to the same protein concentration) were measured in dilution series (as indicated) with this kit. Interpolated concentrations of Survivin adjusted for sample dilution are graphed in pg of Survivin per mL of 1X Diluted extract (mean  $\pm$  SD,  $n = 2$ ).

### 18. SPECIES REACTIVITY

This kit recognizes mouse Survivin protein.

Other species reactivity was not determined.

Please contact our Technical Support team for more information

## 19. 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.
Precipitate in Diluent	Precipitation and/or coagulation of components within the Diluent.	Precipitate can be removed by gently warming the Diluent to 37°C.

### 20. NOTES

## Technical Support

Copyright © 2021 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)