

## ab309282 – Human Tau (4R Isoforms) SimpleStep ELISA® Kit

For the quantitative measurement of Tau (4R Isoforms) in human serum, plasma (heparin), plasma (EDTA), plasma (citrate), cell culture supernatant, CSF.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [www.abcam.com/ab309282](http://www.abcam.com/ab309282)

**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 Standard Preparation and Reagent preparation sections.

### Materials Supplied

Item	Quantity	Storage Condition
Human Tau (4R Isoforms) Capture Antibody 10X	600 µL	+4°C
Human Tau (4R Isoforms) Detector Antibody 10X	600 µL	+4°C
Human Tau (4R Isoforms) Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent CPI2	6 mL	+4°C
Sample Diluent NS	12 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
SimpleStep Pre-Coated 96-Well Microplate	96 wells	+4°C
Plate Seal	1	+4°C

### 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.  
Deionized water.  
Multi- and single-channel pipettes.  
Tubes for standard dilution.  
Plate shaker for all incubation steps.  
Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

### 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

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

**Antibody Cocktail:** Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent CPI2. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent CPI2. Mix thoroughly and gently.

### Standard Preparation

Always prepare a fresh set of standards for every use. Discard working standard dilutions after use as they do not store well. The following section describes the preparation of a standard curve for duplicate measurements (recommended).

1. Reconstitute the Tau (4R Isoforms) standard sample by adding the volume of Sample Diluent NS indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 16,000 pg/mL **Stock Standard** Solution.
2. Add 370 µL of Sample Diluent NS into tube number 1 and 150 µL of Sample Diluent NS into numbers 2-8.
3. Use the **Stock Standard** to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:

Standard #	Dilution Sample	Volume to Dilute (µL)	Volume of Diluent (µL)	Starting Conc. (pg/mL)	Final Conc. (pg/mL)
1	<b>Stock Standard</b>	30	370	16,000	1,200
2	Standard#1	150	150	1,200	600
3	Standard#2	150	150	600	300
4	Standard#3	150	150	300	150
5	Standard#4	150	150	150	75
6	Standard#5	150	150	75	37.5
7	Standard#6	150	150	37.5	18.75
8	Blank Control	0	150	N/A	N/A

## Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	≤ 100%
Plasma - Heparin	≤ 50%
Plasma - EDTA	≤ 50%
Plasma - Citrate	≤ 25%
Cell Culture Media*	≤ 100%
Cerebrospinal Fluid (CSF)	≤ 100%

\*Media is RPMI 1640 containing 10% fetal calf serum.

**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. Assay or dilute serum samples as desired in Sample Diluent NS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

**Plasma** Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. plasma (heparin and EDTA) at least 1:2 into Sample Diluent NS and assay. Dilute plasma (citrate) at least 1:4 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.

**Cell Culture Supernatants** Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants and assay. Or dilute samples into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

**Cerebrospinal Fluid (CSF)** Directly assay cerebrospinal fluid or dilute as desired in Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

## 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.

## Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use.

We recommend that you assay all standards, controls and samples in duplicate

1. Prepare all reagents, working standards, and samples as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, resealed and return to 4°C storage.
3. Add 50 µL of all sample or standard to appropriate wells.
4. Add 50 µL of the Antibody Cocktail to each well.
5. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
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. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
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.

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.
9. Alternative to step 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 µL Stop Solution to each well and recording the OD at 450 nm.

**Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:**

[www.abcam.com/protocols/the-complete-elisa-guide](http://www.abcam.com/protocols/the-complete-elisa-guide)

**For technical support contact information, visit:** [www.abcam.com/contactus](http://www.abcam.com/contactus)

# ab309282 – Human Tau (4R Isoforms) SimpleStep ELISA® Kit

## Additional information

### ASSAY SPECIFICITY

This kit is designed for the quantification of human Tau (4R Isoforms).

The standard protein in this kit is full length, recombinant human Tau (4R Isoforms) isoform E (1N4R).

This kit reacts with all CNS Tau 4R isoforms; it does not react with any Tau 3R isoforms.

Spiked protein experiments were used to validate serum, plasma (heparin), plasma (EDTA), plasma (citrate), cell culture supernatant, and CSF sample types.

Pooled 100% serum, 50% plasma (EDTA and heparin), 25% plasma (citrate), and 100% CSF samples from healthy donors were measured in duplicate. All values were below the detectable range of the assay.

Saliva, urine, and milk samples have not been tested with this kit.

For the measurement of Tau (4R Isoforms) in cell extract and tissue extract sample types, use Human Tau (4R Isoforms) ELISA kit ab303750.

### CROSS REACTIVITY

1250 pg/mL of recombinant human Tau isoform A and 1250 pg/mL of recombinant human Tau isoform B (1N3R) were tested for cross reactivity. No cross reactivity was observed.

### INTERFERENCE

1250 pg/mL of recombinant human Tau isoform A was tested for interference with 1250 pg/mL of recombinant human Tau Isoform E (1N4R). No interference was observed.

1250 pg/mL and 12,250 pg/mL of recombinant human isoform B (1N3R) was tested for interference with 1250 pg/mL of recombinant human Tau isoform E (1N4R). No interference was observed.

### SPECIES REACTIVITY

Other species reactivity was determined by measuring 100% serum samples of various species, interpolating the protein concentrations from the human standard curve. Interpolated protein concentrations are as follows: rat (148.2 pg/mL), mouse (152.7 pg/mL), and cow (108.9 pg/mL). No signal was observed in serum samples from the following species: Monkey.

Other species reactivity not determined.

## CALCULATION

- Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
- Create a standard curve by plotting the average blank control subtracted absorbance value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.

△ Note: Most microplate reader software or graphing software will plot these values and fit a curve to the data. A four parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4 parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.

- Determine the concentration of the target protein in the sample by interpolating the blank control subtracted absorbance values against the standard curve. Multiply the resulting value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.
- Samples generating absorbance values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at an absorbance values less than that of the lowest standard should be retested in a less dilute form.

## 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			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.046	0.046	0.046
18.75	0.090	0.088	0.089
37.5	0.138	0.138	0.138
75	0.221	0.246	0.233
150	0.402	0.420	0.411
300	0.739	0.760	0.749
600	1.422	1.406	1.414
1,200	2.991	2.947	2.969

Table 1. Example of human Tau (4R Isoforms) standard curve in Sample Diluent NS. The Tau (4R Isoforms) standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

## TYPICAL SAMPLE VALUES

### Sensitivity:

The calculated minimal detectable dose (MDD) is 5.26 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=18) and adding 2 standard deviations then extrapolating the corresponding concentration.

### Recovery

3 concentrations of Tau (4R Isoforms) 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% Serum	85	85 - 86
50% Plasma - Heparin	98	94 - 100
50% Plasma - EDTA	86	85 - 87
25% Plasma - Citrate	95	93 - 97
100% Cell Culture Media*	105	103 - 108
50% CSF	105	103 - 106

\*Media is RPMI 1640 containing 10% fetal calf serum.

### 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.

Recombinant Tau (4R Isoforms) was spiked in in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	100% Human Serum	25% Human Plasma (Citrate)	50% Human Plasma (EDTA)	50% Human Plasma (Heparin)	100% Cell Culture Media*	100% CSF
Undiluted	pg/mL	646	459	395	400	961	831
	<b>% Expected value</b>	100	100	100	100	100	100
2	pg/mL	355	274	205	229	420	355
	<b>% Expected value</b>	110	119	104	115	87	110
4	pg/mL	181.2	137.2	97.3	103.6	217.7	181
	<b>% Expected value</b>	112	119	99	104	91	112
8	pg/mL	92.6	65.8	47.5	48.8	117.1	93
	<b>% Expected value</b>	115	115	96	98	97	115
16	pg/mL	45.5	30.5	22.6	24.2	56.3	45
	<b>% Expected value</b>	113	106	92	97	94	113

\*Media is RPMI 1640 containing 10% fetal calf serum.

## Precision

Mean coefficient of variations of interpolated values of Tau (4R Isoforms) from two concentrations of human recombinant Tau 4R protein spiked into 25% human serum within the working range of the assay.

	Intra-assay	Inter-assay
<b>N=</b>	8	3
<b>CV (%)</b>	8.2	7.1

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

[www.abcam.com/protocols/the-complete-elisa-guide](http://www.abcam.com/protocols/the-complete-elisa-guide)

For technical support contact information, visit: [www.abcam.com/contactus](http://www.abcam.com/contactus)

Copyright © 2023 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.  
Version 1a | 2023-04-14