

ab324135 – Mouse/Rat ADAMTS1 SimpleStep ELISA® Kit – Extracellular

For the quantitative measurement of ADAMTS1 in mouse/rat serum, plasma (citrate), plasma (EDTA), cell culture supernatant, and CSF.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: www.abcam.com/ab324135

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 1 x 96 tests	Quantity 10 x 96 tests	Storage Condition
Mouse/Rat ADAMTS1 Capture Antibody 10X	600 µL	10 x 600 µL	+4°C
Mouse/Rat ADAMTS1 Detector Antibody 10X	600 µL	10 x 600 µL	+4°C
Mouse/Rat ADAMTS1 Lyophilized Recombinant Protein	2 Vials	10 x 2 Vials	+4°C
Antibody Diluent CPI2	6 mL	10 x 6 mL	+4°C
Sample Diluent NS	12 mL	2 x 50 mL	+4°C
Wash Buffer PT 10X	20 mL	200 mL	+4°C
TMB Development Solution	12 mL	120 mL	+4°C
Stop Solution	12 mL	120 mL	+4°C
SimpleStep Pre-Coated 96-Well Microplate	96 wells	10 x 96 wells	+4°C
Plate Seal	1	10	+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 ADAMTS1 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 120,000 pg/mL **Stock Standard** Solution.
2. Label eight tubes, Standards 1–8.
3. Add 355 µL of Sample Diluent NS into tube number 1 and 150 µL of Sample Diluent NS into numbers 2-8.
4. 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	Stock Standard	20	355	120,000	6,400
2	Standard#1	150	150	6,400	3,200
3	Standard#2	150	150	3,200	1,600
4	Standard#3	150	150	1,600	800
5	Standard#4	150	150	800	400
6	Standard#5	150	150	400	200
7	Standard#6	150	150	200	100
8	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Mouse Serum	≤ 25%
Mouse Plasma – Citrate	≤ 25%
Mouse Plasma – EDTA	≤ 25%
Mouse Cerebrospinal Fluid	≤ 25%
Rat Serum	≤ 25%
Rat Plasma – Citrate	≤ 25%
Rat Plasma – EDTA	≤ 25%
Rat Cerebrospinal Fluid	≤ 25%
Cell Culture Media*	≤ 50%

*Media is DMEM containing 10% fetal bovine 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. Dilute samples to at least 25% 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 or EDTA. Centrifuge samples at 2,000 x g for 10 minutes. Dilute samples to at least 25% in 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. Note: This kit is incompatible with plasma heparin samples.

Cell Culture Supernatants Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants. Dilute samples to at least 50% in Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Cerebrospinal Fluid (CSF) Dilute samples to at least 25% 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 7 – 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

For technical support contact information, visit: www.abcam.com/contactus

ab324135 – Mouse/Rat ADAMTS1 SimpleStep ELISA® Kit – Extracellular

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of mouse/rat ADAMTS1.

The standard protein in this kit is an N-terminal fragment representing approximately 67% of the full-length mass of mature mouse ADAMTS1.

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

Urine, saliva, and milk samples have not been tested with this kit.

This kit is incompatible with plasma (heparin) samples.

For the measurement of ADAMTS1 in cell extract and tissue extract sample types, use Mouse/Rat ADAMTS1 ELISA kit - Intracellular ab324136.

SPECIES REACTIVITY

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.

Δ 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 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	0.099	0.099	0.099
100	0.159	0.156	0.158
200	0.222	0.217	0.219
400	0.321	0.314	0.318
800	0.514	0.505	0.509
1,600	0.886	0.859	0.873
3,200	1.627	1.596	1.612
6,400	2.966	2.905	2.935

Table 1. Example of mouse/rat ADAMTS1 standard curve in Sample Diluent NS. The ADAMTS1 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 54.23 pg/mL. The MDD was determined by calculating the mean of zero standard replicates (n=24) and adding 2 standard deviations then extrapolating the corresponding concentration.

Recovery

Three concentrations of ADAMTS1 were spiked into the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
25% Mouse Serum	85	80 - 91
25% Mouse Plasma – Citrate	89	82 - 94
25% Mouse Plasma – EDTA	83	81 - 86
25% Mouse Cerebrospinal Fluid	96	90 - 103
25% Rat Serum	83	81 - 85
25% Rat Plasma – Citrate	87	83 - 93
25% Rat Plasma – EDTA	94	92 - 96
25% Rat Cerebrospinal Fluid	102	96 - 109
50% Cell Culture Media*	104	100 - 110

*Media is DMEM containing 10% fetal bovine 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 ADAMTS1 was spiked into the following biological samples and then diluted in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	25% Mouse Serum	25% Mouse Plasma (Citrate)	25% Mouse Plasma (EDTA)	25% Mouse CSF	50% Cell Culture Media
Undiluted	pg/mL	2,514.2	2,648.5	2,641.2	3,732.7	2,961.4
	% Expected value	100	100	100	100	100
2	pg/mL	1,260.2	1,355.0	1,324.0	1,715.8	1,560.0
	% Expected value	100	102	100	92	105
4	pg/mL	715.4	758.0	752.2	836.0	850.1
	% Expected value	114	114	114	90	115
8	pg/mL	360.2	377.5	378.9	482.3	444.1
	% Expected value	115	114	115	103	120
16	pg/mL	NL	NL	NL	254.0	NL
	% Expected value	NL	NL	NL	109	NL

NL – Non-Linear

Recombinant ADAMTS1 was spiked into the following biological samples and then diluted in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	25% Rat Serum	25% Rat Plasma (Citrate)	25% Rat Plasma (EDTA)	25% Rat CSF
Undiluted	pg/mL	2,521.7	2,780.0	2,654.7	3,453.9
	% Expected value	100	100	100	100
2	pg/mL	1,403.4	1,451.9	1,486.6	1,653.4
	% Expected value	111	104	112	96
4	pg/mL	737.0	769.1	775.1	822.8
	% Expected value	117	111	117	95
8	pg/mL	375.2	385.5	392.3	447.0
	% Expected value	119	111	118	104
16	pg/mL	190.8	201.0	192.0	218.4
	% Expected value	121	116	116	101

Precision

Mean coefficient of variations of interpolated values of ADAMTS1 from a single concentration of spiked mouse serum within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	3.0	5.6

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

www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus

Copyright © 2025 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.
Version 1 | 2025-05-02