

ab315311 – Human D-Dimer CatchPoint SimpleStep ELISA® Kit

For the quantitative measurement of D-Dimer in Human serum, plasma (citrate), plasma (EDTA), plasma (heparin), cell culture supernatant, and urine.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: www.abcam.com/ab315311

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.

Limitations: All data, except typical standard curve and sensitivity were collected using the colorimetric version of this kit (ab315310).

Materials Supplied

Item	Quantity	Storage Condition
Human D-Dimer Capture Antibody 10X	600 µL	+4°C
Human D-Dimer Detector Antibody 10X	600 µL	+4°C
Human D-Dimer Lyophilized Native Protein	2 Vials	+4°C
Antibody Diluent 4BI	6 mL	+4°C
Sample Diluent NS	2 x 50 mL	+4°C
Wash Buffer PT 10X	20 mL	+4°C
Stoplight Red Substrate Buffer	12 mL	+4°C
100X Stoplight Red Substrate	120 µL	+4°C
500X Hydrogen Peroxide (H ₂ O ₂ , 3%)	50 µL	+4°C
SimpleStep Pre-Coated Black 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:

Fluorescence microplate reader capable of Ex/Cutoff/Em 530/570/590 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 4BI. 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 4BI. Mix thoroughly and gently.

CatchPoint HRP Development Solution: Just prior to use prepare CatchPoint HRP Development Solution by diluting the 100X Stoplight Red Substrate and the 500X Hydrogen Peroxide in Stoplight Red Substrate Buffer. For example, to make 6 mL of the CatchPoint HRP Development Solution combine 60 µL 100X Stoplight Red Substrate and 12 µL of 500X Hydrogen Peroxide with 5.928 mL Stoplight Red Substrate Buffer. 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 D-Dimer 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 800 ng/mL **Stock Standard** Solution.
2. Label eleven tubes, Standards 1–11.
3. Add 350 µL of Sample Diluent NS into tube number 1 and 150 µL of Sample Diluent NS into numbers 2-11.
4. Use the **Stock Standard** to prepare the following dilution series. Standard #11 contains no protein and is the Blank control:

Standard #	Dilution Sample	Volume to Dilute (µL)	Volume of Diluent (µL)	Starting Conc. (ng/mL)	Final Conc. (ng/mL)
1	Stock Standard	50	350	800	100
2	Standard#1	150	150	100	50
3	Standard#2	150	150	50	25
4	Standard#3	150	150	25	12.5
5	Standard#4	150	150	12.5	6.25
6	Standard#5	150	150	6.25	3.13
7	Standard#6	150	150	3.13	1.56
8	Standard#7	150	150	1.56	0.78
9	Standard#8	150	150	0.78	0.39
10	Standard#9	150	150	0.39	0.20
11	Blank Control	0	150	0	0

Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Serum	1:3,200 - 1:200
Plasma – EDTA	1:200 - 1:25
Plasma – Citrate	1:20,000 - 1:1,250
Plasma – Heparin	1:100 - 1:25
Urine	≤25%
Cell Culture Supernatant	≤50%

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 at least 1:200 into 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. Dilute EDTA and heparin samples at least 1:25 and citrate samples at least 1:1,250 into Sample Diluent NS. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

Urine Centrifuge urine at 2,000 x g for 10 minutes to remove debris. Dilute samples at least 1:4 into Sample Diluent NS and assay. Store un-diluted urine samples at -20°C or below. 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. Dilute samples at least 1:2 into 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, reseal 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 prepared CatchPoint HRP Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm. Further optimization of incubation time vs signal strength can be performed if needed.
8. Record the fluorescence at Ex/Cutoff/Em 530/570/590 nm. If using a Molecular Devices' plate reader supported by SoftMax® Pro software, a preconfigured protocol for these CatchPoint SimpleStep ELISA Kits is available with all the protocol and analysis settings at www.softmaxpro.org.
9. Analyze the data as described below.

Mode	Fluorescence
Instrument settings:	Endpoint
Excitation:	530 nm
Cutoff:	570 nm
Emission:	590 nm
Sensitivity:	6 flashes/read or 200ms
PMT:	Auto
Auto calibrate:	On
Read:	Top
Read Height:	1*

Note For microplate readers with Pre-Read Optimization option, the Read Height as well as Microplate Optimization is recommended before the first read.

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 © 2024 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.

Version 1 | 2024-1-12

ab315311 – Human D-Dimer CatchPoint SimpleStep ELISA® Kit

Additional information

ASSAY SPECIFICITY

This kit is designed for the quantification of Human D-Dimer and other high molecular weight fibrin degradation products. This kit does not detect Human Fibrinogen.

The standard protein in this kit is natively purified Human D-Dimer.

Native signal was detected in serum, plasma (citrate), plasma (EDTA), and plasma (heparin).

Spiked protein experiments were used to validate cell culture supernatant and urine sample types.

Saliva, milk, CSF, cell extract, and tissue extract samples have not been tested with this kit.

CROSS REACTIVITY

The following proteins were prepared at 200 ng/mL and tested for cross reactivity. No cross reactivity was observed.

- native human Fibrinogen
- native human Plasminogen
- recombinant human t-Plasminogen Activator (tPA)
- recombinant human Angiostatin
- recombinant Human Serpin E1 (PAI-1)

INTERFERENCE

The following proteins were prepared at 200 ng/mL and tested for interference with 25 ng/mL of native Human D-Dimer. No interference was observed.

- native human Fibrinogen
- native human Plasminogen
- recombinant human t-Plasminogen Activator (tPA)
- recombinant human Angiostatin
- recombinant Human Serpin E1 (PAI-1)

SPECIES REACTIVITY

Other species reactivity was determined by measuring 10% serum samples of various species.

No signal was observed for the following species: Monkey, Mouse, Rat, Cow

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 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 (ng/mL)	RFU		Mean RFU
	1	2	
0	275582	289704	282643
0.20	593626	606575	600101
0.39	958330	855097	906714
0.78	1744709	1722107	1733408
1.56	3269324	3514436	3391880
3.13	10094286	8139713	9117000
6.25	21321396	17065378	19193387
12.5	37105080	39473936	38289508
25	89834048	76497960	83166004
50	137844128	145585664	141714896
100	256881488	252881440	254881464

Table 1. Example of Human D-Dimer standard curve in Sample Diluent NS. The D-Dimer 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 0.065 ng/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 D-Dimer 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 (%)
1:400 Serum	95	92 - 98
1:200 Plasma – EDTA	105	101 - 113
1:5,000 Plasma – Citrate	106	99 - 113
1:80 Plasma – Heparin	101	97 - 104
25% Urine	90	85 - 93
50% Cell Culture Media*	92	86 - 99

*Media is RPMI 1640 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.

Native D-Dimer was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	1:200 Human Serum	1:1,250 Human Plasma (Citrate)	1:25 Human Plasma (EDTA)	1:25 Human Plasma (Heparin)
Undiluted	ng/mL	14.4	29.9	42.3	41.8
	% Expected value	100	100	100	100
2	ng/mL	6.7	16.9	21.6	19.7
	% Expected value	93	113	102	94
4	ng/mL	3.3	7.5	8.7	9.1
	% Expected value	91	100	82	87
8	ng/mL	1.9	3.9	4.3	NL
	% Expected value	103	103	81	NL
16	ng/mL	0.8	2.1	NL	NL
	% Expected value	90	113	NL	NL

NL – Non-Linear

Native D-Dimer was spiked into the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Dilution Factor	Interpolated value	25% Human Urine	50% Cell Culture Media
Undiluted	ng/mL	23.2	20.4
	% Expected value	100	100
2	ng/mL	12.1	11.0
	% Expected value	104	108
4	ng/mL	6.3	6.1
	% Expected value	109	120
8	ng/mL	2.6	3.2
	% Expected value	91	124
16	ng/mL	1.4	1.5
	% Expected value	99	114

Precision

Mean coefficient of variations of interpolated values of D-Dimer from a single concentration of human serum within the working range of the assay.

	Intra-assay	Inter-assay
N=	8	3
CV (%)	7.4	3.8

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 © 2024 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.
Version 1 | 2024-1-12