



SERPINA1 (Human) ELISA Kit

Catalog Number KA0459

96 assays

Version: 16

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Background	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	5
Assay Protocol	6
Reagent Preparation	6
Sample Preparation	6
Assay Procedure	8
Data Analysis	10
Calculation of Results	10
Performance Characteristics	11
Resources	13
Troubleshooting	13
References	14
Plate Layout	15

Introduction

Background

Alpha-1-antitrypsin (A1AT) is a protein that protects the lungs. The liver usually makes the protein and releases it into the bloodstream. A1AT is a major protease inhibitor that controls tissue degradation. A reduction of A1AT levels can cause a change in collagen metabolism (1). A1AT inhibits neutrophil elastase release into the lungs during inflammatory states (2). A1AT deficiency is an uncommon genetic disease (3) that can lead to emphysema (4), hepatitis, cirrhosis (5), and chronic obstructive pulmonary disease (COPD) [6].

Principle of the Assay

The SERPINA1 (Human) ELISA Kit is designed for detection of A1AT in human plasma, serum, milk, urine, saliva, CSF, cell culture, and cell lysate samples. This assay employs a quantitative sandwich enzyme immunoassay technique that measures human A1AT in approximately 4 hours. A polyclonal antibody specific for human A1AT has been pre-coated onto a 96-well microplate with removable strips. A1AT in standards and samples is sandwiched by the immobilized antibody and a biotinylated polyclonal antibody specific for human A1AT, which is recognized by a streptavidin-peroxidase (SP) conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

General Information

Materials Supplied

List of component

Component	Amount
Human alpha-1-Antitrypsin Microplate: A 96-well polystyrene microplate (12 strips of 8 wells) coated with a polyclonal antibody against human A1AT.	96 (8x12) wells
Sealing Tapes: Pressure sensitive sealing tapes that can be cut to fit the format of the individual assay.	3 slides
Human alpha-1-Antitrypsin Standard: Human A1AT in a buffered protein base, lyophilized.	27.5 ng
Biotinylated Human alpha-1-Antitrypsin Antibody (50x): A 50-fold concentrated biotinylated polyclonal antibody against human A1AT.	120 µL
MIX Diluent Concentrate (10x): A 10-fold concentrated buffered protein base.	30 mL
Wash Buffer Concentrate (20x): A 20-fold concentrated buffered surfactant.	30 mL x 2
SP Conjugate (100x): A 100-fold concentrate.	80 µL
Chromogen Substrate (1x): A stabilized peroxidase chromogen substrate tetramethylbenzidine.	7 mL
Stop Solution (1x): A 0.5 N hydrochloric acid solution to stop the chromogen substrate reaction.	11 mL

Storage Instruction

- ✓ Upon arrival, immediately store components of the kit at recommended temperatures up to the expiration date.
- ✓ Store SP Conjugate and Biotinylated Antibody at -20°C.
- ✓ Store Microplate, Diluent Concentrate (10x), Wash Buffer, Stop Solution, and Chromogen Substrate at 2-8°C.
- ✓ Unused microplate wells may be returned to the foil pouch with the desiccant pack and resealed. May be stored for up to 30 days in a vacuum desiccator.
- ✓ Store Standard at 2-8°C before reconstituting with Diluent and at -20°C after reconstituting with Diluent.

Materials Required but Not Supplied

- ✓ Microplate reader capable of measuring absorbance at 450 nm.
- ✓ Pipettes (1-20 µL, 20-200 µL, 200-1000 µL and multiple channel).
- ✓ Deionized or distilled reagent grade water

Precautions for Use

- ✓ This product is for Research Use Only and is not intended for use in diagnostic procedures.
- ✓ Prepare all reagents (diluent buffer, wash buffer, standard, biotinylated antibody, and SP conjugate) as instructed, prior to running the assay.
- ✓ Prepare all samples prior to running the assay. The dilution factors for the samples are suggested in this insert. However, the user should determine the optimal dilution factor.
- ✓ Spin down the SP conjugate vial and the biotinylated antibody vial before opening and using contents.
- ✓ The Stop Solution is an acidic solution.
- ✓ The kit should not be used beyond the expiration date.

Assay Protocol

Reagent Preparation

- ✓ Freshly dilute all reagents and bring all reagents to room temperature before use.
- ✓ MIX Diluent Concentrate (10x): Dilute the MIX Diluent Concentrate 10-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved. Store for up to 30 days at 2-8°C.
- ✓ Human alpha-1-Antitrypsin Standard: Reconstitute the Human alpha-1-Antitrypsin Standard (27.5 ng) with 1.1 mL of MIX Diluent to generate a 25 ng/mL standard stock solution. Allow the vial to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare duplicate or triplicate standard points by serially diluting from the standard stock solution (25 ng/mL) 2-fold with equal volume of MIX Diluent to produce 12.5, 6.25, 3.125, 1.563, 0.781, and 0.391 ng/mL solutions.. MIX Diluent serves as the zero standard (0 ng/mL). Any remaining stock solution should be stored at -20°C and used within 30 days. Avoid repeated freeze-thaw cycles.

Standard Point	Dilution	[A1AT] (ng/mL)
P1	1 part Standard (25 ng/mL)	25
P2	1 part P1 + 1 part MIX Diluent	12.5
P3	1 part P2 + 1 part MIX Diluent	6.25
P4	1 part P3 + 1 part MIX Diluent	3.125
P5	1 part P4 + 1 part MIX Diluent	1.563
P6	1 part P5 + 1 part MIX Diluent	0.781
P7	1 part P6 + 1 part MIX Diluent	0.391
P8	MIX Diluent	0.0

- ✓ Biotinylated Human alpha-1-Antitrypsin Antibody (50x): Spin down the antibody briefly and dilute the desired amount of the antibody 50-fold with MIX Diluent to produce a 1x solution. The undiluted antibody should be stored at -20°C.
- ✓ Wash Buffer Concentrate (20x): Dilute the Wash Buffer Concentrate 20-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved.
- ✓ SP Conjugate (100x): Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 100-fold with MIX Diluent to produce a 1x solution. The undiluted conjugate should be stored at -20°C.

Sample Preparation

- ✓ Plasma: Collect plasma using one-tenth volume of 0.1 M sodium citrate as an anticoagulant. Centrifuge samples at 3000 x g for 10 minutes and collect plasma. A 200000-fold sample dilution is suggested into

MIX Diluent; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles. (EDTA or Heparin can also be used as an anticoagulant).

- ✓ Serum: Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 3000 x g for 10 minutes and remove serum. A 200000-fold sample dilution is suggested into MIX Diluent; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ Milk: Collect milk using sample tube. Centrifuge samples at 800 x g for 10 minutes. A 2000-fold sample dilution is suggested into MIX Diluent; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ Urine: Collect urine using sample pot. Centrifuge samples at 800 x g for 10 minutes. A 20-fold sample dilution is suggested into MIX Diluent or within the range of 2x - 200x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ Saliva: Collect saliva using sample tube. Centrifuge samples at 800 x g for 10 minutes. A 400-fold sample dilution is suggested into MIX Diluent or within the range of 40x - 4000x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ CSF: Collect cerebrospinal fluid (CSF) using sample pot. Centrifuge samples at 3000 x g for 10 minutes. A 4000-fold sample dilution is suggested into MIX Diluent; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -80°C for up to 3 months. Avoid repeated freeze-thaw cycles.
- ✓ Cell Culture Supernatants: Centrifuge cell culture media at 1500 rpm for 10 minutes at 4°C to remove debris and collect supernatants. Samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.
- ✓ Cell Lysate: Rinse cell with cold PBS and then scrape the cell into a tube with 5 mL of cold PBS and 0.5 M EDTA. Centrifuge suspension at 1500 rpm for 10 minutes at 4°C and aspirate supernatant. Resuspend pellet in ice-cold Lysis Buffer (10 mM Tris, pH 8.0, 130 mM NaCl, 1% Triton X-100, protease inhibitor cocktail). For every 1×10^6 cells, add approximately 100 μ L of ice-cold Lysis Buffer. Incubate on ice for 60 minutes. Centrifuge at 13000 rpm for 30 minutes at 4°C and collect supernatant. Samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.
- ✓ *Note: Applicable samples may also include biofluids, cell culture, and tissue homogenates. If necessary, user should determine optimal dilution factor depending on application needs.*

- ✓ Refer to Dilution Guidelines for further instruction.

Guidelines for Dilutions of 100-fold or Greater (for reference only; please follow the insert for specific dilution suggested)	
100x	10000x
A. 4 μ L sample: 396 μ L buffer (100x) = 100-fold dilution Assuming the needed volume is less than or equal to 400 μ L.	A. 4 μ L sample: 396 μ L buffer (100x) B. 4 μ L of A : 396 μ L buffer (100x) = 10000-fold dilution Assuming the needed volume is less than or equal to 400 μ L.
1000x	100000x
A. 4 μ L sample: 396 μ L buffer (100x) B. 24 μ L of A : 216 μ L buffer (10x) = 1000- fold Assuming the needed volume is less than or equal to 240 μ L.	A. 4 μ L sample: 396 μ L buffer (100x) B. 4 μ L of A : 396 μ L buffer (100x) C. 24 μ L of B : 216 μ L buffer (10x) = 100000-fold Assuming the needed volume is less than or equal to 240 μ L.

Assay Procedure

1. Prepare all reagents, standard solutions, and samples as instructed. Bring all reagents to room temperature before use. The assay is performed at room temperature (20-25°C).
2. Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccant inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
3. Add 50 μ L of Human alpha-1-Antitrypsin Standard or sample to each well. Gently tap plate to thoroughly coat the well. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
4. Wash the microplate manually or automatically using a microplate washer. Invert the plate and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If washing manually, wash five times with 200 μ L of Wash Buffer per well. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a microplate washer, wash six times with 300 μ L of Wash Buffer per well; invert the plate and hit 4-5 times on absorbent material to completely remove the liquid.
5. Add 50 μ L of Biotinylated Human alpha-1-Antitrypsin Antibody to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 1 hour.
6. Wash the microplate as described above.
7. Add 50 μ L of SP Conjugate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles

that may have formed. Cover wells with a sealing tape and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.

8. Wash the microplate as described above.
9. Add 50 μ L of Chromogen Substrate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate in ambient light for 25 minutes or until the optimal blue color density develops.
10. Add 50 μ L of Stop Solution to each well. The color will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.
11. Read the absorbance on a microplate reader at a wavelength of 450 nm immediately. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. Please note that some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

✓ Assay summary

1. Add 50 μ L of Standard or Sample per well. Incubate 2 hours.
2. Wash, then add 50 μ L of Biotinylated Antibody per well. Incubate 1 hour.
3. Wash, then add 50 μ L of SP Conjugate per well. Incubate 30 minutes.
4. Wash, then add 50 μ L of Chromogen Substrate per well. Incubate 25 minutes.
5. Add 50 μ L of Stop Solution per well. Read at 450 nm immediately.

Data Analysis

Calculation of Results

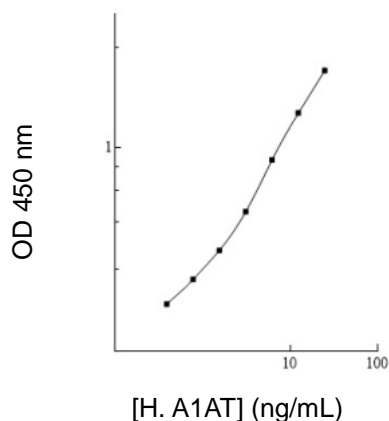
- ✓ Calculate the mean value of the duplicate or triplicate readings for each standard and sample.
- ✓ To generate a standard curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance (OD) on the y-axis. The best fit line can be determined by regression analysis using log-log or four-parameter logistic curve fit.
- ✓ Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.
- ✓ Typical Data

The typical data is provided for reference only. Individual laboratory means many vary from the values listed. Variations between laboratories may be caused by technique differences.

Standard Point	ng/mL	OD	Average OD
P1	25	2.243 2.199	2.221
P2	12.5	1.456 1.394	1.425
P3	6.25	0.877 0.875	0.876
P4	3.125	0.518 0.506	0.512
P5	1.563	0.343 0.341	0.342
P6	0.781	0.258 0.248	0.253
P7	0.391	0.200 0.192	0.196
P8	0.0	0.153 0.149	0.151
Sample: Pooled Normal Sodium Citrate Plasma (200000x)		0.910 0.882	0.896

- ✓ Standard Curve

The curve is provided for illustration only. A standard curve should be generated each time the assay is performed.



✓ **Reference Value**

Normal human A1AT plasma levels range from 700-1900 µg/mL.

Plasma and serum samples from healthy adults were tested (n=40). On average, human A1AT level was 1338 µg/mL.

Sample	n	Average Value (µg/mL)
Pooled Normal Plasma	10	1307
Normal Plasma	20	1284
Pooled Normal Serun	10	1423

Performance Characteristics

- ✓ Kit standard has been calibrated against WHO international Standard.
- ✓ The minimum detectable dose of human A1AT as calculated by 2SD from the mean of a zero standard was established to be 0.31 ng/mL.
- ✓ Intra-assay precision was determined by testing three plasma samples twenty times in one assay.
- ✓ Inter-assay precision was determined by testing three plasma samples in twenty assays.

	Intra-Assay Precision			Inter-Assay Precision		
Sample	1	2	3	1	2	3
n	20	20	20	20	20	20
CV (%)	4.8%	5.0%	4.1%	9.8%	10.2%	9.1%
Average CV (%)	4.6%			9.7%		

✓ Spiking Recovery

Recovery was determined by spiking two plasma samples with different A1AT concentrations.

Sample	Unspiked Sample (ng/mL)	Spiked Value (ng/mL)	Expected	Observed	Recovery (%)
1	3.4	1.5	4.9	5.4	110%
		5.0	8.4	9.2	110%
		10.0	13.4	13.3	99%
2	6.2	1.5	7.7	8.0	104%
		5.0	11.2	11.1	99%
		10.0	16.2	15.9	98%
Average Recovery (%)					103%

✓ Linearity

Plasma and serum samples were serially diluted to test for linearity.

Average Percentage of Expected Value (%)		
Sample Dilution	Plasma	Serum
100000x	102%	110%
200000x	99%	100%
400000x	99%	93%

✓ Cross-Reactivity

Species	Cross Reactivity (%)
Canine	None
Bovine	None
Monkey	< 5%
Mouse	None
Rat	None
Swine	None
Rabbit	None

10% FBS in culture media will not affect the assay.

Resources

Troubleshooting

Issue	Causes	Course of Action
Low Precision	Use of improper components	Check the expiration date listed before use. Do not interchange components from different lots.
	Improper wash step	Check that the correct wash buffer is being used. Check that all wells are empty after aspiration. Check that the microplate washer is dispensing properly. If washing by pipette, check for proper pipetting technique.
	Splashing of reagents while loading wells	Pipette properly in a controlled and careful manner.
	Inconsistent volumes loaded into wells	Pipette properly in a controlled and careful manner. Check pipette calibration. Check pipette for proper performance.
	Insufficient mixing of reagent dilutions	Thoroughly agitate the lyophilized components after reconstitution. Thoroughly mix dilutions.
	Improperly sealed microplate	Check the microplate pouch for proper sealing. Check that the microplate pouch has no punctures. Check that three desiccants are inside the microplate pouch prior to sealing.
Unexpectedly Low or High Signal Intensity	Microplate was left unattended between steps	Each step of the procedure should be performed uninterrupted.
	Omission of step	Consult the provided procedure for complete list of steps.
	Step performed in incorrect order	Consult the provided procedure for the correct order.
	Insufficient amount of reagents added to wells	Check pipette calibration. Check pipette for proper performance.
	Wash step was skipped	Consult the provided procedure for all wash steps.
	Improper wash buffer	Check that the correct wash buffer is being used.
	Improper reagent preparation	Consult reagent preparation section for the correct dilutions of all reagents.
	Insufficient or prolonged incubation periods	Consult the provided procedure for correct incubation time.

Deficient Standard Curve Fit	Non-optimal sample dilution	<p>Sandwich ELISA: If samples generate OD values higher than the highest standard point (P1), dilute samples further and repeat the assay.</p> <p>Competitive ELISA: If samples generate OD values lower than the highest standard point (P1), dilute samples further and repeat the assay.</p> <p>User should determine the optimal dilution factor for samples.</p>
	Contamination of reagents	A new tip must be used for each addition of different samples or reagents during the assay procedure.
	Contents of wells evaporate	Verify that the sealing film is firmly in place before placing the assay in the incubator or at room temperature.
	Improper pipetting	<p>Pipette properly in a controlled and careful manner.</p> <p>Check pipette calibration.</p> <p>Check pipette for proper performance.</p>
	Insufficient mixing of reagent dilutions	<p>Thoroughly agitate the lyophilized components after reconstitution.</p> <p>Thoroughly mix dilutions.</p>

References

1. Hauck EW *et al.* (2004) *Eur Urol.* 46(5):623-8; discussion 628.
2. Chappell *et al.* (2004) *Hum Mutat.* 24(6):535-6.
3. Strange C *et al.* (2006) *Respiration.* 73(2):185-90.
4. Abboud RT *et al.* (2005) *Treat Respir Med.* 4(1):1-8.
5. Kok KF *et al.* (2005) *Ned Tijdschr Geneesk.* 149(37):2057-61.
6. Teramoto S. (2007) *Intern Med.* 46(2):77-9.

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H