



# ANCA-P (MPO) ELISA Kit

Catalog Number KA1086

96 assays

Version: 05

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Intended Use .....	3
Background .....	3
Principle of the Assay .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	5
<b>Assay Protocol .....</b>	<b>7</b>
Reagent Preparation .....	7
Sample Preparation .....	7
Assay Procedure .....	8
<b>Data Analysis.....</b>	<b>9</b>
Calculation of Results.....	9
Performance Characteristics .....	9
<b>Resources.....</b>	<b>11</b>
References .....	11
Plate Layout .....	12

## **Introduction**

### **Intended Use**

The ANCA-P (MPO) ELISA Kit is a test system for the quantitative measurement of IgG class autoantibodies against myeloperoxidase (MPO) in human serum or plasma. This product is intended for research use only

### **Background**

Anti-neutrophil cytoplasmic antibodies (ANCA) are diagnostic markers for ANCA-associated vasculitides. Anti-MPO differentiates microscopic polyangiitis (MPA) and eosinophilic granulomatosis with polyangiitis (EGPA) The test supports differential diagnosis of vasculitis, when used in conjunction with other clinical and laboratory findings.

### **Principle of the Assay**

Highly purified myeloperoxidase (MPO) is bound to microwells.

The determination is based on an indirect enzyme linked immune reaction with the following steps:

Specific antibodies in the sample bind to the antigen coated on the surface of the reaction wells. After incubation, a washing step removes unbound and unspecifically bound serum or plasma components. Subsequently added enzyme conjugate binds to the immobilized antibody-antigen-complexes. After incubation, a second washing step removes unbound enzyme conjugate. After addition of substrate solution the bound enzyme conjugate hydrolyses the substrate forming a blue colored product. Addition of an acid stops the reaction generating a yellow end-product. The intensity of the yellow color correlates with the concentration of the antibody-antigen-complex and can be measured photometrically at 450 nm.

## General Information

### Materials Supplied

List of component

Component	Amount
Divisible microplate: Consisting of 12 modules of 8 wells each. Ready to use.	96 (8x12) wells
Calibrators A-F (0, 5, 10, 20, 40, 100 U/mL), containing MPO antibodies in a serum/buffer matrix (PBS, BSA, detergent, NaN <sub>3</sub> 0.09%), yellow. Ready to use.	1.5 mL x 6
Control positive (1) and negative (2), containing MPO antibodies in a serum/buffer matrix (PBS, BSA, detergent, NaN <sub>3</sub> 0.09%), yellow. Ready to use. The concentration is specified on the certificate of analysis.	1.5 mL x 2
Sample buffer, containing PBS, BSA, detergent, preservative NaN <sub>3</sub> 0.09%, yellow, 5x conc.	20 mL
Enzyme Conjugate, containing anti-human IgG antibodies, HRP labelled; PBS, BSA, detergent, preservative ProClin 300 0.05%, light red. Ready to use.	15 mL
TMB Substrate; containing 3,3',5,5'-Tetramethylbenzidin, colorless. Ready to use.	15 mL
Stop Solution: Contains acid. Ready to use.	15 mL
Wash Buffer, containing Tris, detergent, preservative NaN <sub>3</sub> 0.09%; 50x conc.	20 mL

### Storage Instruction

- ✓ Store the kit at 2-8°C in the dark.
- ✓ Do not expose reagents to heat, sun or strong light during storage and usage.
- ✓ Store microplate sealed and desiccated in the clip bag provided.
- ✓ Unopened reagents are stable until expiration of the kit. See labels for individual batch.
- ✓ Diluted Wash Buffer and Sample Buffer are stable for at least 30 days when stored at 2-8°C. We recommend consumption on the same day.

### Materials Required but Not Supplied

- ✓ Microplate reader capable of endpoint measurements at 450 nm; optional: reference filter at 620 nm.
- ✓ Data reduction software
- ✓ Multi-Channel Dispenser or repeatable pipet for 100 µL
- ✓ Vortex mixer
- ✓ Pipets for 10 µL, 100 µL and 1000 µL
- ✓ Laboratory timing device
- ✓ Distilled or deionized water
- ✓ Measuring cylinder for 1000 and 100 mL

- ✓ Plastic container for storage of the wash solution

This ELISA assay is suitable for use on open automated ELISA processors. Each assay has to be validated on the respective automated system. Detailed information is provided upon request.

### **Precautions for Use**

#### ✓ Procedural Notes

- Do not use kit components beyond their expiration dates.
- Do not interchange kit components from different lots and products.
- All materials must be at room temperature (20-28°C) prior to use.
- Prepare all reagents and samples. Once standard, perform the test without interruption.
- Double determinations may be done. By this means pipetting errors may become obvious.
- Perform the assay steps only in the order indicated.
- Always use fresh sample dilutions.
- Pipette all reagents and samples into the bottom of the wells.
- To avoid carryover or contamination, change the tip between samples and different kit controls.
- Wash microwells thoroughly and remove the last droplets of Wash Buffer.
- All incubation steps must be accurately timed.
- Do not re-use microplate wells.

#### ✓ Warnings and Precautions

- All reagents of this kit are intended for professional research use only.
- Components containing human serum were tested and found negative for HBsAg, HCV, HIV1 and HIV2 by FDA approved methods. No test can guarantee the absence of HBsAg, HCV, HIV1 or HIV2 and so all human serum based reagents in this kit must be handled as though capable of transmitting infection.
- Bovine serum albumin (BSA) used in components has been tested for BSE and found negative.
- Avoid contact with the TMB (3,3',5,5'-Tetramethyl-benzidine).
- Stop Solution contains acid, classification is non-hazardous. Avoid contact with skin.
- Controls, Calibrators, Sample Buffer and Wash Buffer contain sodium azide ( $\text{NaN}_3$ ) 0.09% as preservative. This concentration is classified as non-hazardous.
- Enzyme conjugate contains ProClin 300 0.05% as preservative. This concentration is classified as non-hazardous.
- During handling of all reagents, controls and serum samples observe the existing regulations for laboratory safety regulations and good laboratory practice:
- First aid measures: In case of skin contact, immediately wash thoroughly with water and soap. Remove contaminated clothing and shoes and wash before reuse. If system fluid comes into contact with skin, wash thoroughly with water. After contact with the eyes carefully rinse the opened eye with running water for at least 10 minutes. Get medical attention if necessary.
- Personal precautions, protective equipment and emergency procedures:

- Observe laboratory safety regulations. Avoid contact with skin and eyes. Do not swallow. Do not pipette by mouth. Do not eat, drink, smoke or apply makeup in areas where specimens or kit reagents are handled. When spilled, absorb with an inert material and put the spilled material in an appropriate waste disposal.
- Exposure controls/personal protection: Wear protective gloves of nitrile rubber or natural latex. Wear protective glasses. Used according to intended use no dangerous reactions known.
- Conditions to avoid: Since substrate solution is light-sensitive. Store in the dark.
- For disposal of laboratory waste the national or regional legislation has to be observed.

Observe the guidelines for performing quality control in medical laboratories by assaying control sera.

✓ Limitations of the procedure

The abnormal and normal reference ranges for antibodies in samples should be regarded as recommendations only. Each laboratory should establish its own ranges according to ISO 15189 or other applicable laboratory guidelines.

## Assay Protocol

### Reagent Preparation

✓ Wash Buffer

Dilute the contents of one vial of the buffered wash solution concentrate (50x) with distilled or deionised water to a final volume of 1000 mL prior to use.

✓ Sample buffer

Prior to use dilute the contents (20 mL) of one vial of sample buffer 5x concentrate with distilled or deionized water to a final volume of 100 mL.

### Sample Preparation

✓ Specimen collection, storage and handling

1. Collect whole blood specimens using acceptable medical techniques to avoid hemolysis.
2. Allow blood to clot and separate the serum or plasma by centrifugation.
3. Test serum should be clear and non-hemolyzed. Contamination by hemolysis or lipemia should be avoided, but does not interfere with this assay.
4. Specimens may be refrigerated at 2-8°C for up to five days or stored at -20°C up to six months.
5. Avoid repetitive freezing and thawing of serum or plasma samples. This may result in variable loss of antibody activity.
6. Testing of heat-inactivated sera is not recommended.

✓ Preparation of Samples

Dilute samples 1:100 before the assay: Put 990 µL of prediluted sample buffer in a polystyrene tube and add 10 µL of sample. Mix well.

*Note: Calibrators/Controls are ready to use and need not be diluted.*

### **Assay Procedure**

Prepare enough microplate modules for all calibrators/controls and samples.

1. Pipette 100  $\mu$ L of calibrators, controls and prediluted samples into the wells.
2. Incubate for 30 minutes at room temperature (20-28°C).
3. Discard the contents of the microwells and wash 3 times with 300  $\mu$ L of wash solution.
4. Dispense 100  $\mu$ L of enzyme conjugate into each well.
5. Incubate for 15 minutes at room temperature.
6. Discard the contents of the microwells and wash 3 times with 300  $\mu$ L of wash solution.
7. Dispense 100  $\mu$ L of TMB substrate solution into each well.
8. Incubate for 15 minutes at room temperature.
9. Add 100  $\mu$ L of stop solution to each well of the modules.
10. Incubate for 5 minutes at room temperature.
11. Read the optical density at 450 nm (reference 600-690 nm) and calculate the results. The developed color is stable for at least 30 minutes. Read optical densities during this time.

## Data Analysis

### Calculation of Results

✓ Calculation of results

For quantitative results plot the optical density of each calibrator versus the calibrator concentration to create a calibration curve. The concentration of samples may then be estimated from the calibration curve by interpolation.

Using data reduction software a 4-Parameter-Fit with lin-log coordinates for optical density and concentration is the data reduction method of choice.

✓ Validation

Test results are valid if the optical densities at 450 nm for calibrators / controls and the results for controls comply with the reference ranges indicated on the Certificate of Analysis enclosed in each test kit. If these quality control criteria are not met the assay run is invalid and should be repeated.

### Performance Characteristics

✓ Calibration

This assay system is calibrated in relative arbitrary units, since no international reference preparation is available for this assay.

✓ Measuring range

The calculation range of this ELISA assay is 0-100 U/mL.

✓ Expected values

In a normal range study with samples from healthy blood donors the following ranges have been established with this ELISA assay: Cut-off 5 U/mL.

✓ Interpretation of results

Negative: < 5 U/mL

Positive: ≥ 5 U/mL

✓ Linearity

Samples containing high levels of specific antibody were serially diluted in sample buffer to demonstrate the dynamic range of the assay and the upper/lower end of linearity. Activity for each dilution was calculated from the calibration curve using a 4-Parameter-Fit with lin-log coordinates.

Sample No	Dilution	Observed [U/mL]	Expected [U/mL]	Observed/Expected (%)
1	1:100	87.3	87.3	100
	1:200	44.1	43.7	101
	1:400	21.5	21.8	99
	1:800	9.7	10.9	89
	1:1600	5.0	5.5	91

2	1:100	79.9	79.9	100
	1:200	39.3	40.0	98
	1:400	19.0	20.0	95
	1:800	8.5	10.0	85
	1:1600	4.3	5.0	86

✓ Limit of detection

Functional sensitivity was determined to be: 0.5 U/mL.

✓ Reproducibility

Intra-assay precision: Coefficient of variation (CV) was calculated for each of three samples from the results of 24 determinations in a single run. Results for precision-within-assay are shown in the table below.

Inter-assay precision: Coefficient of variation (CV) was calculated for each of three samples from the results of 6 determinations in 5 different runs. Results for run-to-run precision are shown in the table below.

Intra-Assay		
Sample	Mean U/mL	CV%
1	7.5	6.4
2	30.2	4.1
3	59.9	3.1

Inter-Assay		
Sample	Mean U/mL	CV%
1	7.0	5.0
2	33.8	4.9
3	78.3	6.3

✓ Interfering substances

No interference has been observed with haemolytic (up to 1000 mg/dL) or lipemic (up to 3 g/dL triglycerides) sera or plasma, or bilirubin (up to 40 mg/dL) containing sera or plasma. Nor have any interfering effects been observed with the use of anticoagulants (Citrate, EDTA, Heparin). However for practical reasons it is recommended that grossly hemolyzed or lipemic samples should be avoided.

✓ Study results

Study population	n	n Pos	%
Crescendic glomerulonephritis	55	53	96.4
Morbus Wegener (cANCA pos)	20	1	5.0
Non-ANCA kidney disease	10	1	10.0
Normal human sera	120	3	2.5

	Pos	Neg	
Pos	54	5	
Neg	1	145	
	55	150	205

Sensitivity: 98.2%

Specificity: 96.7%

Overall agreement: 97.1%

## Resources

### References

1. Jennette, J.C. and Falk, R.J. Antineutrophil Cytoplasmic Autoantibodies and Associated Diseases: a Review. Am. J. Kidney Dis. 1990, Vol. XV, No. 6: 517 - 529.
2. Gross, W.L. et al. Antineutrophil Cytoplasmic Autoantibody-Associated Diseases: A Rheumatologist's Perspective. Am. J. Kidney Dis. 1991, Vol. XVIII, No. 2: 175 - 179.
3. Wieslander, J. How are Antineutrophil Cytoplasmic Autoantibodies Detected ? Am. J. Kidney Dis. 1991, Vol. XVIII, No. 2: 154 - 158.
4. Lesavre, P. Antineutrophil cytoplasmic antibodies antigen specificity. Am. J. Kidney Dis. 1991, Vol. XVIII, No. 2: 159 - 163.
5. Hagen, E.C. et al. Antineutrophil cytoplasmic autoantibodies: a review of the antigens involved, the assays, and the clinical and possible pathogenic consequences. Blood 1993, Vol.81: 1996 - 2000.
6. Gross, W.L. et al. Immunodiagnostische und immunopathogenetische Bedeutung von Anti-Neutrophilen-Cytoplasma-Antikörpern. Deutsche Medizinische Wochenschrift 1993, Vol. 118: 191 - 199.

**Plate Layout**

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