



# CA19-9 (Human) ELISA Kit

Catalog Number KA0207

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>6</b>
Reagent Preparation .....	6
Sample Preparation .....	6
Assay Procedure .....	6
<b>Data Analysis.....</b>	<b>8</b>
Calculation of Results.....	8
Performance Characteristics .....	8
<b>Resources.....</b>	<b>9</b>
References .....	9
Plate Layout .....	10

## **Introduction**

### **Intended Use**

For the quantitative determination of the Cancer Antigen CA19-9 concentration in human serum.

### **Background**

A group of mucin type glycoprotein Sialosyl Lewis Antigens (SLA), such as CA19-9 and CA19-5, has come to be recognized as circulating cancer associated antigens for gastrointestinal cancer. CA19-9 represents the most important and basic carbohydrate tumor marker. The immunohistologic distribution of CA19-9 in tissues is consistent with the quantitative determination of higher CA19-9 concentrations in cancer than in normal or inflamed tissues. Recently reports indicates that the serum CA19-9 level is frequently elevated in the serum of subjects with various gastrointestinal malignancies, such as pancreatic, colorectal, gastric and hepatic carcinomas. Together with CEA, elevated CA19-9 is suggestive of gallbladder neoplasm in the setting of inflammatory gallbladder disease. This tumor-associated antigen may also be elevated in some non-malignant conditions. Research studies demonstrate that serum CA 19-9 values may have utility in monitoring subjects with the above-mentioned diagnosed malignancies. It has been shown that a persistent elevation in serum CA19-9 value following treatment may be indicative of occult metastatic and/or residual disease. A persistently rising serum CA 19-9 value may be associated with progressive malignant disease and poor therapeutic response. A declining CA 19-9 value may be indicative of a favorable prognosis and good response to treatment.

### **Principle of the Assay**

The CA19-9 (Human) ELISA Kit is based on the principle of a solid phase enzyme-linked immunosorbent assay. The assay system utilizes a monoclonal antibody directed against a distinct antigenic determinant on the intact CA19-9 molecule is used for solid phase immobilization (on the microtiter wells). Another CA 19-9 monoclonal antibody conjugated to horseradish peroxidase (HRP) is in the antibody-enzyme conjugate solution. The test sample is allowed to react sequentially with the two antibodies, resulting in the CA19-9 molecules being sandwiched between the solid phase and enzyme-linked antibodies. After two separate incubation steps at 37°C for 90 minutes, the wells are washed with wash buffer to remove unbound labeled antibodies. A solution of TMB Reagent is added and incubated for 20 minutes, resulting in the development of a blue color. The color development is stopped with the addition of Stop Solution changing the color to yellow. The concentration of CA19-9 is directly proportional to the color intensity of the test sample. Absorbance is measured spectrophotometrically at 450 nm.

## General Information

### Materials Supplied

List of component

Component	Amount
Murine monoclonal anti-CA19-9 coated 96 well microtiter plate.	96 wells
CA19-9 reference standards, containing 0, 25, 75, 150, 300, and 600 U/mL CA19-9, liquid, ready to use	0.5 mL x 6 vials
CA 19-9 Assay Buffer	13 mL
Enzyme Conjugate Concentrate (12x)	1.1 mL
CA 19-9 Conjugate Diluent	13 mL
Wash Buffer Concentrate (20x)	50 mL
TMB Reagent (One-Step)	11 mL
Stop Solution (1 N HCl)	11 mL

### Storage Instruction

Unopened test kits should be stored at 2-8°C upon receipt and the microtiter plate should be kept in a sealed bag with desiccants to minimize exposure to damp air. Opened test kits will remain stable until the expiration date shown, provided it is stored as described above.

### Materials Required but Not Supplied

- ✓ Precision pipettes and tips: 10 µL and 100 µL
- ✓ Distilled water
- ✓ Vortex mixer
- ✓ Absorbent paper or paper towel
- ✓ Graph paper
- ✓ A microtiter plate reader with a bandwidth of 10 nm or less and an optical density range of 0-2 OD or greater at a wavelength of 450 nm is acceptable for use in absorbance measurement.

**Precautions for Use**

For research use only. Not for use in diagnostic procedures.

- Limitation of procedure
- ✓ Reliable and reproducible results will be obtained when the assay procedure is carried out with a complete understanding of the package insert instructions and with adherence to good laboratory practice.
- ✓ The wash procedure is critical. Insufficient washing will result in poor precision and falsely elevated absorbance readings.
- ✓ Serum samples demonstrating gross lipemia, gross hemolysis, or turbidity should not be used with this test.

## Assay Protocol

### Reagent Preparation

All reagents should be brought to room temperature (18-25°C) before use.

- Wash Buffer (1x):

Add 50 mL of Wash Buffer (20X) to 950 mL of DI water. The diluted Wash Buffer is stable at 2-8°C for 30 days. Mix well before use.

*Note: Any crystals that may be present due to high salt concentration must be redissolved at room temperature before making the dilution.*

- Working CA 19-9 Conjugate Reagent:

For 3.0 mL, which is more than enough for 24 wells: Add 0.25 mL of Conjugate Concentrate (12x) to 2.75 mL of the Enzyme Conjugate Diluent (1:11 dilution) and mix well.

For 6.0 mL, which is more than enough for 48 wells: Add 0.5 mL of Conjugate Concentrate (12x) to 5.5 mL of the Enzyme Conjugate Diluent (1:11 dilution) and mix well.

For 9.0 mL, which is more than enough for 72 wells: Add 0.75 mL of Conjugate Concentrate (12 x) to 8.25 mL of the Enzyme Conjugate Diluent (1:11 dilution) and mix well.

For 12.0 mL, which is more than enough for 96 wells: Add 1.0 mL of Conjugate Concentrate (12x) to 11.0 mL of the Enzyme Conjugate Diluent (1:11 dilution) and mix well.

*The Working CA 19-9 Conjugate Reagent needs to be prepared freshly every time before use.*

*The amount of conjugate diluted depends on your assay size. Discard the excess after use.*

### Sample Preparation

Serum should be prepared from a whole blood specimen obtained by acceptable techniques. This kit is for use with serum samples without additives only.

### Assay Procedure

1. Secure the desired number of coated wells in the holder.
2. Dispense 10 µL of CA19-9 standards, specimens, and controls into appropriate wells.
3. Dispense 100 µL of CA 19-9 Assay Buffer (green-color solution) into each well.
4. Thoroughly mix for 30 seconds. It is very important to mix them completely.
5. Incubate at 37°C for 90 minutes.
6. Remove the incubation mixture by emptying the plate content into a waste container.
7. Rinse and flick the microtiter wells 5 times with diluted Wash Buffer (1x).
8. Strike the microtiter plate sharply onto absorbent paper or paper towels to remove all residual water droplets.
9. Dispense 100 µL of the Working Conjugate Reagent into each well. Mix gently for 30 seconds.

10. Incubate at 37°C for 90 minutes.
11. Remove the incubation mixture by emptying the plate content into a waste container.
12. Rinse and flick the microtiter wells 5 times with diluted Wash Buffer (1x).
13. Strike the microtiter plate sharply onto absorbent paper or paper towels to remove all residual water droplets.
14. Dispense 100  $\mu$ L of the TMB Reagent into each well. Gently mix for 10 seconds.
15. Incubate at room temperature in the dark for 20 minutes without shaking.
16. Stop the reaction by adding 100  $\mu$ L of Stop Solution to each well.
17. Gently mix for 30 seconds. It is important to make sure that all the blue color changes to yellow color completely.
18. Read the optical density at 450 nm with a microtiter plate reader within 15 minutes

## Data Analysis

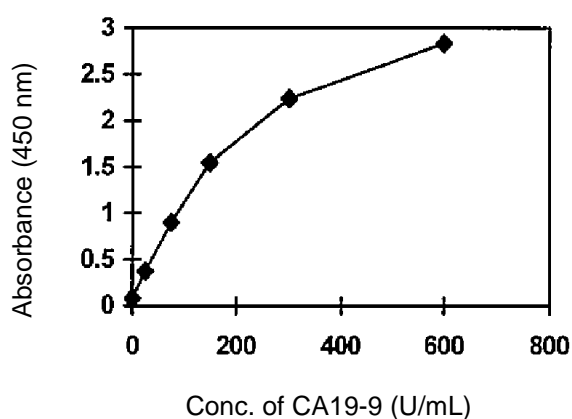
### Calculation of Results

1. Calculate the average absorbance values (A<sub>450</sub>) for each set of reference standards, control, and samples.
2. Construct a standard curve by plotting the mean absorbance obtained for each reference standard against its concentration in U/mL via best fit quadratic on linear graph paper, with absorbance on the vertical (y) axis and concentration on the horizontal (x) axis.
3. Using the mean absorbance value for each sample, determine the corresponding concentration of CA19-9 in U/mL from the standard curve.

✓ Example of standard curve

Results of a typical standard run with optical density readings at 450 nm shown in the Y axis against CA19-9 concentrations shown in the X axis. This standard curve is for the purpose of illustration only, and should not be used to calculate unknowns. Each user should obtain his or her own data and standard curve in each experiment.

CA19-9 (U/mL)	Absorbance (450 nm)
0	0.075
25	0.373
75	0.900
150	1.543
300	2.237
600	2.832



### Performance Characteristics

1. Expected values

Health men and women are expected to have CA19-9 assay value below 35 U/mL. The minimum detectable concentration of CA19-9 in this assay is estimated to be 10 U/mL.



## Resources

### References

1. Glenn, J., Steinberg, W.M., Kurtzman, S.H., et al. Evaluation of the utility of a radioimmunoassay for serum CA 19-9 level in patients before and after treatment of carcinoma of the pancreas. *J. Clin. Oncol.* 1988; 6:462-8.
2. Hayakawa, T., Kondo, T., Shibata, T. et al. Sensitive serum markers for detecting pancreatic cancer. *Cancer* 1988; 61:1827-31.
3. Koprowski, H., Herly, M., Steplewski, Z., et al. Specific antigen in serum of patients with colon carcinoma. *Science* 1981; 212:53-5.
4. Malesci, A., Tommasini, M.A., Bonato, C. et al. Determination of CA19-9 antigen in serum and pancreatic juice for differential diagnosis of pancreatic adenocarcinoma from chronic pancreatitis. *Gastroenterology* 1987; 92:60-7.
5. Safi, F., Roscher, R., Bittner, R., et al. High sensitivity and specificity of CA 19-9 for pancreatic carcinoma in comparison to chronic pancreatitis. Serological and immunohistochemical findings. *Pancreas* 1987; 2:398-403.
6. Steinberg, W. The clinical utility of CA 19-9 tumor associated antigen. *American J. of Gastroenterology* 1990; 85:350-355.
7. Steinberg, W.M., Gelfand, R., Anderson, K.K., et al. Comparison of the sensitivity and specificity of the CA 19-9 and carcinoembryonic antigen assays in detecting cancer of the pancreas. *Gastroenterology* 1986; 90:343-9.
8. Takasaki, H., Uchida, E., Tempero, M.A., et al. Correlative study on expression of CA 19-9 and DU-Pan-2 in tumor tissue and in serum of pancreatic cancer patients. *Cancer Res.* 1988; 48:1435-8.
9. Tatsuta, M., Yamamura, H., Iishi H., et al. Values of CA19-9 in the serum, pure pancreatic juice and aspirated pancreatic material in the diagnosis of malignant pancreatic tumor. *Cancer* 1985; 56:2669-73.
10. Wang, T.H. Lin, J.W., Chen, D.S., et al. Noninvasive diagnosis of advanced pancreatic cancer by real-time ultrasonography, carcinoembryonic antigen, and carbohydrate antigen 19-9. *Pancreas* 1986; 1:219-23.
11. Strom BL, Maislin G, West SL, et al. Serum CEA and CA19-9: potential future diagnostic or screening tests for gallbladder cancer? *Int. J. Cancer* 1990; 45:821.

**Plate Layout**

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