



Canine RBP4 ELISA Kit

Catalog Number NBP2-60464

Novus kits are guaranteed for 6 months from date of receipt

Assay Summary

Step 1. Add 50 μ l of Standard or Sample per well.
Incubate 2 hours.

Step 2. Wash, then add 50 μ l of Biotinylated Antibody per well.
Incubate 1 hour.

Step 3. Wash, then add 50 μ l of SP Conjugate per well.
Incubate 30 minutes.

Step 4. Wash, then add 50 μ l of Chromogen Substrate per well.
Incubate 10 minutes.

Step 5. Add 50 μ l of Stop Solution per well.
Read at 450 nm immediately.

Symbol Key



Consult instructions for use.

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

Assay Template

Canine Retinol-Binding Protein 4 (RBP4) ELISA Kit

Catalog No. NBP2-60464

Sample insert for reference use only

Introduction

Serum retinol-binding protein (RBP4), secreted by the liver and adipocytes, is implicated in systemic insulin resistance. RBP4 transports retinol and circulates in the plasma by binding to the larger transthyretin (TTR) homotetramer, forming a protein complex that reduces renal clearance of RBP4. In insulin-resistant ob/ob mice, urinary fractional excretion of RBP4 was reduced, which is consistent with increased retention, while TTR level is elevated (1). RBP4 is encoded by the *RBP4* gene which maps to chromosome 10q23-q24 (2, 3). Transgenic overexpression of human RBP4 or injection of recombinant RBP4 in normal mice causes insulin resistance. Conversely, genetic deletion of RBP4 enhances insulin sensitivity. Increasing serum RBP4 induces hepatic expression of the gluconeogenic enzyme phosphoenolpyruvate carboxykinase and impairs insulin signaling in muscle tissue (4). Expression of RBP4 is induced in adipose tissue as a consequence of decreased glucose transporter GLUT4 expression.

Principle of the Assay

The Canine RBP4 ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for detection of canine RBP4 in **plasma, serum, urine, and cell culture samples**. This assay employs a quantitative **sandwich enzyme immunoassay** that measures canine RBP4 in less than 4 hours. A polyclonal antibody specific for canine RBP4 has been pre-coated onto a 96-well microplate with removable strips. Canine RBP4 in standards and samples is sandwiched by the immobilized antibody and biotinylated polyclonal antibody specific for canine RBP4, which is recognized by a streptavidin-peroxidase conjugate. All unbound material is then washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

Caution and Warning

- This product is for **Research Use Only** and is **Not For Use In Diagnostic Procedures**.
- Prepare all reagents (working 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.

Reagents

- **Canine RBP4 Microplate:** A 96-well polystyrene microplate (12 strips of 8 wells) coated with a polyclonal antibody against canine RBP4.
- **Sealing Tapes:** Each kit contains 3 precut, pressure sensitive sealing tapes, which can be cut to fit the format of the individual assay.
- **Canine RBP4 Standard:** Canine RBP4 in a buffered protein base (50 ng, lyophilized).
- **Biotinylated Canine RBP4 Antibody (50x):** A 50-fold concentrated biotinylated polyclonal antibody against canine RBP4 (140 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, 2 bottles).
- **Streptavidin-Peroxidase Conjugate (SP Conjugate):** A 100-fold concentrate (80 µl).
- **Chromogen Substrate:** A ready-to-use stabilized peroxidase chromogen substrate tetramethylbenzidine (8 ml).
- **Stop Solution:** A 0.5 N hydrochloric acid to stop the chromogen substrate reaction (12 ml).

Storage Condition

- 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 packs and resealed. May be stored for up to 30 days in a vacuum desiccator.
- Diluent (1x) may be stored for up to 30 days at 2-8°C.
- Store Standard at 2-8°C before reconstituting with Diluent and at -20°C after reconstituting with Diluent.

Other Supplies Required

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

Sample Collection, Preparation, and Storage

- **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. Dilute samples 1:5000 into MIX Diluent. 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. Dilute samples 1:5000 into MIX Diluent and assay. 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. Dilute urine 1:4 into MIX Diluent and assay. Samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.
- **Cell Culture Supernatants:** Centrifuge cell culture media at 3000 x g for 10 minutes to remove debris. Collect supernatants and assay. Store samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Refer to Sample Dilution Guidelines below for further instruction.

Guidelines for Dilutions of 1:100 or Greater <i>(for reference only; please follow the insert for specific dilution suggested)</i>	
1:100	1:10000
<p>A) 4 μl sample: 396 μl buffer(100x) = 100 fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 400 μl.</i></p>	<p>A) 4 μl sample : 396 μl buffer (100x) B) 4 μl of A : 396 μl buffer (100x) = 10000 fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 400 μl.</i></p>
1:1000	1:100000
<p>A) 4 μl sample : 396 μl buffer (100x) B) 24 μl of A : 216 μl buffer (10x) = 1000 fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 240 μl.</i></p>	<p>A) 4 μl sample : 396 μl buffer (100x) B) 4 μl of A : 396 μl buffer (100x) C) 24 μl of B : 216 μl buffer (10x) = 100000 fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 240 μl.</i></p>

Reagent Preparation

- Freshly dilute all reagents and bring all reagents to room temperature before use.
- **MIX Diluent Concentrate (10x):** If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the MIX Diluent Concentrate 1:10 with reagent grade water. Store for up to 30 days at 2-8°C.
- **Canine RBP4 Standard:** Reconstitute the 50 ng of Canine RBP4 Standard with 1 ml of MIX Diluent to generate a 50 ng/ml standard stock solution. Allow the standard to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare duplicate or triplicate standard points by serially diluting the standard stock solution (50 ng/ml) 1:2 with MIX Diluent to produce 25, 12.5, 6.25, 3.125, 1.563, and 0.781 ng/ml solutions. MIX Diluent serves as the zero standard (0 ng/ml). Any remaining solution should be frozen at -20°C and used within 30 days.

Standard Point	Dilution	[Canine RBP4] (ng/ml)
P1	1 part Standard (50 ng/ml)	50.00
P2	1 part P1 + 1 part MIX Diluent	25.00
P3	1 part P2 + 1 part MIX Diluent	12.50
P4	1 part P3 + 1 part MIX Diluent	6.250
P5	1 part P4 + 1 part MIX Diluent	3.125
P6	1 part P5 + 1 part MIX Diluent	1.563
P7	1 part P6 + 1 part MIX Diluent	0.781
P8	MIX Diluent	0.000

- **Biotinylated Canine RBP4 Antibody (50x):** Spin down the antibody briefly and dilute the desired amount of the antibody 1:50 with MIX Diluent. Any remaining solution should be frozen at -20°C.
- **Wash Buffer Concentrate (20x):** If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the Wash Buffer Concentrate 1:20 with reagent grade water.
- **SP Conjugate (100x):** Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 1:100 with MIX Diluent. Any remaining solution should be frozen at -20°C.

Assay Procedure

- 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).
- Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccants inside. Reseal the pouch

securely to minimize exposure to water vapor and store in a vacuum desiccator.

- Add 50 μ l of Canine RBP4 Standard or sample per well. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
- Wash five times with 200 μ l of Wash Buffer manually. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a machine, wash six times with 300 μ l of Wash Buffer and then invert the plate, decanting the contents; hit 4-5 times on absorbent material to completely remove the liquid.
- Add 50 μ l of Biotinylated Canine RBP4 Antibody to each well and incubate for 1 hour.
- Wash the microplate as described above.
- Add 50 μ l of Streptavidin-Peroxidase Conjugate per well and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
- Wash the microplate as described above.
- Add 50 μ l of Chromogen Substrate per well and incubate for 10 minutes or till the optimal blue color density develops. Gently tap plate to ensure thorough mixing and break the bubbles in the well with pipette tip.
- Add 50 μ l of Stop Solution to each well. The color will change from blue to yellow.
- 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.

Data Analysis

- 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 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

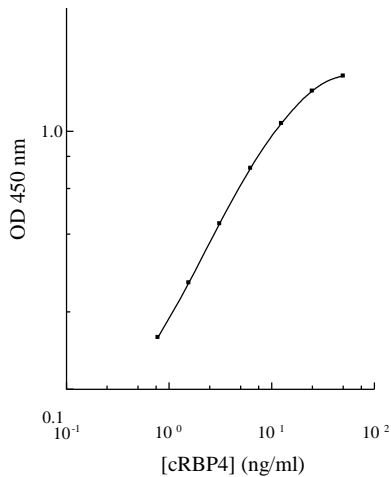
- The typical data is provided for reference only. Individual laboratory means may vary from the values listed. Variations between laboratories may be caused by technique differences.

Standard Point	ng/ml	OD	Average OD
P1	50.00	1.630 1.642	1.636
P2	25.00	1.391 1.350	1.371
P3	12.50	1.000 1.017	1.008
P4	6.250	0.648 0.634	0.641
P5	3.125	0.372 0.350	0.361
P6	1.563	0.215 0.219	0.217
P7	0.781	0.124 0.126	0.125
P8	0.000	0.037 0.033	0.035

Standard Curve

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

Canine RBP4 Standard Curve



Performance Characteristics

- The minimum detectable dose of canine RBP4 as calculated by 2SD from the mean of a zero standard was established to be 0.7 ng/ml.
- Intra-assay precision was determined by testing replicates of three plasma samples in one assay.
- Inter-assay precision was determined by testing three plasma samples in twenty assays.

Sample	Intra-Assay Precision			Inter-Assay Precision		
	1	2	3	1	2	3
n	20	20	20	20	20	20
CV (%)	3.5%	3.9%	3.6%	9.7%	9.5%	9.8%
Average CV (%)	3.7%			9.7%		

Recovery

Standard Added Value	1.5 – 12.5 ng/ml
Recovery %	83 – 112%
Average Recovery %	96%

Linearity

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

Average Percentage of Expected Value (%)		
Sample Dilution	Plasma	Serum
1:2500	91%	94%
1:5000	97%	99%
1:10000	106%	104%

Cross-Reactivity

Species	Cross Reactivity (%)
Bovine	None
Canine	100%
Monkey	None
Mouse	None
Rat	None
Swine	None
Human	None
Rabbit	None

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

Troubleshooting

Issue	Causes	Course of Action
Low Precision	Use of expired components	<ul style="list-style-type: none"> • Check the expiration date listed before use. • Do not interchange components from different lots.
	Improper wash step	<ul style="list-style-type: none"> • Check that the correct wash buffer is being used. • Check that all wells are dry 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	<ul style="list-style-type: none"> • Pipette properly in a controlled and careful manner.
	Inconsistent volumes loaded into wells	<ul style="list-style-type: none"> • Pipette properly in a controlled and careful manner. • Check pipette calibration. • Check pipette for proper performance.
	Insufficient mixing of reagent dilutions	<ul style="list-style-type: none"> • Thoroughly agitate the lyophilized components after reconstitution. • Thoroughly mix dilutions.
	Improperly sealed microplate	<ul style="list-style-type: none"> • 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	<ul style="list-style-type: none"> • Each step of the procedure should be performed uninterrupted.
	Omission of step	<ul style="list-style-type: none"> • Consult the provided procedure for complete list of steps.
	Steps performed in incorrect order	<ul style="list-style-type: none"> • Consult the provided procedure for the correct order.
	Insufficient amount of reagents added to wells	<ul style="list-style-type: none"> • Check pipette calibration. • Check pipette for proper performance.
	Wash step was skipped	<ul style="list-style-type: none"> • Consult the provided procedure for all wash steps.
	Improper wash buffer	<ul style="list-style-type: none"> • Check that the correct wash buffer is being used.
	Improper reagent preparation	<ul style="list-style-type: none"> • Consult reagent preparation section for the correct dilutions of all reagents.
Insufficient or prolonged incubation periods	<ul style="list-style-type: none"> • Consult the provided procedure for correct incubation time. 	
Deficient Standard Curve Fit	Non-optimal sample dilution	<ul style="list-style-type: none"> • Sandwich ELISA: If samples generate OD values higher than the highest standard point (P1), dilute samples further and repeat the assay. • Competitive ELISA: If samples generate OD values lower than the highest standard point (P1), dilute samples further and repeat the assay. • User should determine the optimal dilution factor for samples.
	Contamination of reagents	<ul style="list-style-type: none"> • A new tip must be used for each addition of different samples or reagents during the assay procedure.
	Contents of wells evaporate	<ul style="list-style-type: none"> • Verify that the sealing film is firmly in place before placing the assay in the incubator or at room temperature.

	Improper pipetting	<ul style="list-style-type: none"> • Pipette properly in a controlled and careful manner. • Check pipette calibration. • Check pipette for proper performance.
	Insufficient mixing of reagent dilutions	<ul style="list-style-type: none"> • Thoroughly agitate the lyophilized components after reconstitution. • Thoroughly mix dilutions.

References

- (1) Mody N *et al.* (2008) *Am. J. Physiol Endocrinol Metab.* 294(4): E785-793
- (2) Meigs JB *et al.* (2002) *Diabetes* 51:833–840
- (3) Duggirala R *et al.* (1998) *Diabetes* 47 (Suppl. 1): A170
- (4) Yang Q *et al.* (2005) *Nature* 436 (7049): 356-362

Related Products

- Human RBP ELISA Kit (Urine, Milk, and Cell Culture samples)
- Human RBP ELISA Kit (Plasma and Serum samples)
- Human RBP4 ELISA Kit (Plasma, Serum, Urine, Saliva, Milk, and Cell Cultures samples)
- Mouse RBP4 ELISA Kit (Plasma, Serum, Urine, and Cell Culture samples)