

ARBOR ASSAYS™  
Interactive Assay Solutions™



# DetectX<sup>®</sup>

## HIGH SENSITIVITY RBP Enzyme Immunoassay Kit

1 Plate Kit Catalog Number KU04-H1

5 Plate Kit Catalog Number KU04-H5

Species Independent

Formerly Named Urinary RBP EIA Kit

Assay Buffer Change: See pages 5 & 8

Sample Types Validated:

**Dried Blood Spot and Urine**

Please read this insert completely prior to using the product.  
For research use only. Not for use in diagnostic procedures.

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

# TABLE OF CONTENTS

|  |       |
|--|-------|
| Background                                   | 3     |
| Assay Principle                              | 4     |
| Related Products                             | 4     |
| Supplied Components                          | 5     |
| Storage Instructions                         | 5     |
| Other Materials Required                     | 6     |
| Precautions                                  | 6     |
| Sample Types                                 | 7     |
| Sample Preparation                           | 7     |
| Reagent Preparation                          | 8     |
| Assay Protocol                               | 9     |
| Calculation of Results                       | 10    |
| Typical Data                                 | 10-11 |
| Validation Data Sensitivity, Linearity, etc. | 11-13 |
| Sample Values                                | 14    |
| Warranty & Contact Information               | 15    |
| Plate Layout Sheet                           | 16    |

## BACKGROUND

Retinol binding protein (RBP) is from a family of structurally related proteins that bind small hydrophobic molecules such as bile pigments, steroids, odorants, etc<sup>1</sup>. RBP is a 21 kDa highly conserved, single-chain glycoprotein, consisting of 182 amino acids with 3 disulfide bonds, that has a hydrophobic pocket which binds retinol (vitamin A).

RBP binds retinol in a 1:1 stoichiometry, which serves to not only solubilize retinol but also protect it from oxidation. When in serum, the majority of RBP bound with retinol is reversibly complexed with transthyretin (prealbumin)<sup>2,3</sup>. This complex then transports retinol to specific receptors of various tissues in the body. Vitamin A status is reflected by serum concentration as it is hemostatically controlled and does not fall until stores are dramatically reduced<sup>4,5</sup>.

RBP has also been shown to be a useful marker for renal function<sup>6</sup> as it is totally filtered by the glomeruli and reabsorbed by proximal tubules<sup>7</sup>. This has made urinary RBP (uRBP) a tool to study renal function in heart<sup>8</sup> or kidney<sup>9</sup> transplant recipients, type 1 and 2 diabetics<sup>10</sup>, and in people exposed to uranium from mining operations<sup>11</sup>. Measurement of uRBP levels has also been useful in detection and characterization of diseases including hypertension<sup>12</sup> and certain cancers<sup>13,14</sup>.

1. Blaner WS., "Retinol binding protein: the serum transport protein for vitamin A.", *Endocr Rev.* 1989, 10(3):308–16.
2. Wolf G., "Multiple functions of vitamin A.", *Physiol. Rev.* 1984, 64(3):873-937.
3. Petersen PA., "Characteristics of a Vitamin A-transporting Protein Complex Occuring in Human Serum.", *J. Biol. Chem.* 1971 246:34-43.
4. Goodman DS., Blaner WS., in *The Retinoids*, eds. Sporn MB., Roberts AB., Goodman DS., (Orlando: Academic Press, 1984) vol.2, 1-39.
5. Olson JA., "Vitamin A, retinoids and carotenoids.", *Modern Nutrition in Health and Disease*, eds. Shils ME, Olson JA, Shike M, (Philadelphia: Lea & Febiger, 1994) 8th ed. 287–307.
6. Peterson PA., and Berggård I., "Isolation and Properties of a Human Retinol-transporting Protein.", *J. Biol. Chem.* 1971, 246:25-33.
7. Bernard AM., and Lauwerys RR., "Retinol Bining Protein in Urine: A More Practical Index than Urinary  $\beta$ 2-microglobulin for the Screening of Renal Tubular Function.", *Clin. Chem.* 1981, 27:1781-2.
8. Câmara NO., Matos AC., Rodrigues DA., et al., "Early Detection of Heart Transplant patients with Increased Risk of Cyclosporin Nephrotoxicity.", *The Lancet*, 2001, 357: 856-857.
9. Hosaka B., Park SI., Felipe CR., Garcia RG., Machado PG., Pereira AB., Tedesco-Silva H., and Medina Pestana JO., "Predictive Value of Urinary Retinol Binding Protein for Graft Dysfunction after Kidney Transplantation." *Transplant Proc.*, 2003, 35:1341-1343.
10. Abahusain MA., Wright J., Dickerson JWT., and de Vol EB., "Retinol,  $\alpha$ -Tocopherol and Carotenoids in Diabetes." *Am J Clin Nutr.*, 2004, 79:218-25.
11. Wyatt SA., Reitz LV., Croley TR., Hawkins D., Barrett E., McKeown A., Powell N., West A., Hamner T. and Royster MO., "Biological Monitoring of Uranium Exposure in South Central Virginia." *J Exp Sci Environ Epidemiem.*, 2008, 18:59-75.
12. Bang LE., Holm J., and Svendsen TL. "Retinol-Binding protein and Transferrin in Urine: New Markers of Renal Function in Essential Hypertension and White Coat Hypertension" *Am. J. Hypertens*, 1996, 9:1024-1028.
13. Moise AR, Noy N, Palczewski K, Blaner WS. "Delivery of retinoid-based therapies to target tissues." *Biochemistry*, 2007, 9:1024-1028.
14. Gavrilo V., Yermiahu T., Gorodischer R., "Renal pathology and retinol status in multiple myeloma patients." *Kidney Int.*, 2006, 69(1):173-7.

## ASSAY PRINCIPLE

The DetectX® High Sensitivity Retinol Binding Protein (RBP) kit is designed to quantitatively measure RBP present in dried blood spot (DBS) and urine samples. Please read the complete kit insert before performing this assay. A RBP standard is provided to generate a standard curve for the assay and all samples should be read off the standard curve. Standards or diluted samples are pipetted into a clear microtiter plate coated with an antibody to capture rabbit antibodies. A RBP-peroxidase conjugate is added to the standards and samples in the wells. The binding reaction is initiated by the addition of the RBP polyclonal antibody to each well. After an hour incubation the plate is washed and substrate is added. The substrate reacts with the bound RBP-peroxidase conjugate. After a short incubation, the reaction is stopped and the intensity of the generated color is detected in a microtiter plate reader capable of measuring 450 nm wavelength. The concentration of the RBP in the sample is calculated, after making a suitable correction for the dilution of the sample, using software available with most plate readers.

## RELATED PRODUCTS

| Kits  | Catalog No. |
|---|-------------|
| BCA Protein Dual Range Colorimetric Detection Kit | K041-H1     |
| Human Cystatin C EIA Kit                          | K012-H1     |
| Serum Creatinine Detection Kit                    | KB02-H1     |
| Thiol Detection Kit                               | K005-F1     |
| Urea Nitrogen (BUN) Detection Kits                | K024-H1/H5  |
| Urinary Creatinine Detection Kits                 | K002-H1/H5  |

## SUPPLIED COMPONENTS

### Coated Clear 96 Well Plates

A clear plastic microplate(s) with break-apart strips coated with goat anti-rabbit IgG.

Kit KU04-H1 or -H5 1 or 5 Each

Catalog Number X016-1EA

### RBP Standard

A stock solution of native human RBP at 20 µg/mL.

Kit KU04-H1 or -H5 60 µL or 240 µL

Catalog Number C014-60UL or -240UL

### DetectX® RBP Antibody

A polyclonal antibody specific for RBP.

Kit KU04-H1 or -H5 3 mL or 13 mL

Catalog Number C011-3ML or -13ML

### DetectX® RBP Conjugate

A RBP-peroxidase conjugate.

Kit KU04-H1 or -H5 3 mL or 13 mL

Catalog Number C015-3ML or -13ML

### Assay Buffer Concentrate *now supplied as a concentrate*

A 5X concentrate that must be diluted with deionized or distilled water.

Kit KU04-H1 or -H5 28 mL or 55 mL

Catalog Number X053-28ML or -55ML

### Wash Buffer Concentrate

A 20X concentrate that should be diluted with deionized or distilled water.

Kit KU04-H1 or -H5 30 mL or 125 mL

Catalog Number X007-30ML or -125ML

### TMB Substrate

Kit KU04-H1 or -H5 11 mL or 55 mL

Catalog Number X019-11ML or -55ML

### Stop Solution

A 1M solution of hydrochloric acid. **CAUSTIC.**

Kit KU04-H1 or -H5 5 mL or 25 mL

Catalog Number X020-5ML or -25ML

### Plate Sealer

Kit KU04-H1 or -H5 1 or 5 Each

Catalog Number X002-1EA

## STORAGE INSTRUCTIONS

All components of this kit should be stored at 4°C until the expiration date of the kit.

## OTHER MATERIALS REQUIRED

Distilled or deionized water.

A microplate shaker and a microplate washer.

Repeater pipet with disposable tips capable of dispensing 25  $\mu$ L, 50  $\mu$ L, and 100  $\mu$ L.

Colorimetric 96 well microplate reader capable of reading optical density at 450 nm.

Software for converting raw relative optical density readings from the plate reader and carrying out four parameter logistic curve (4PLC) fitting. Contact your plate reader manufacturer for details.

## PRECAUTIONS

As with all such products, this kit should only be used by qualified personnel who have had laboratory safety instruction. The complete insert should be read and understood before attempting to use the product.

The antibody coated plate needs to be stored desiccated. The silica gel pack included in the foil ziploc bag will keep the plate dry. The silica gel pack will turn from blue to pink if the ziploc has not been closed properly.

The RBP Standard is purified from a human source and as such, should be treated as potentially hazardous. Proper safety procedures must be followed.

This kit utilizes a peroxidase-based readout system. Buffers, including other manufacturers Wash Buffers, containing sodium azide will inhibit color production from the enzyme. Make sure **all** buffers used for samples are **azide free**. Ensure that any plate washing system is rinsed well with deionized water prior to using the supplied Wash Buffer as prepared on Page 8.

The Stop Solution is acid. The solution should not come in contact with skin or eyes. Take appropriate precautions when handling this reagent.

**Laboratory temperature is important.** Please make sure that the kit incubates at a temperature between 22°C and 24°C.



## SAMPLE TYPES

This assay has been fully validated for human urine and dried blood spot samples. Samples containing visible particulate should be centrifuged prior to using.

RBP is a highly conserved protein and we have shown that this kit may measure RBP from sources other than human. Please see page 14 for details of other urine samples tested. The end user should evaluate recoveries of RBP in other urine samples being tested.

## SAMPLE PREPARATION

### Urine Samples

Samples must be diluted 1:2 by adding one part of urine to one part diluted Assay Buffer prior to running in the kit. Any samples with RBP concentrations greater than the standard curve range should be diluted further with diluted Assay Buffer to obtain readings within the standard curve. Samples that are too dilute to be measured should be concentrated prior to measuring in the assay.

### Dried Blood Spots (DBS)

Dried blood spot (DBS) samples should be prepared according to the 2007 Clinical Chemistry paper by Masako Fujita, et al, vol. 53 (11), page 1972-1975. Briefly, whole blood is spotted onto Whatman 309 filter paper and thoroughly dried at room temperature. These can be stored desiccated at  $\leq 4^{\circ}\text{C}$  until use. DBS samples, 1/4" or 1/8", are punched out into clean plastic tubes with caps. For each 1/4" DBS a minimum of 240  $\mu\text{L}$  of diluted Assay Buffer is added. For each 1/8" DBS a minimum of 60  $\mu\text{L}$  of diluted Assay Buffer is added. This is equivalent to a 1:40 dilution of the sample. The tubes are capped and left at  $4^{\circ}\text{C}$  overnight. The following morning, the red solution can be run without centrifugation or further dilution.

For calculation purposes a 1/8 inch DBS is considered to contain 1.5  $\mu\text{L}$  of whole blood sample.

The dilution of any samples that fall outside the standard range should be adjusted to allow samples to read within the standard curve.

**Use all samples within 2 hours of dilution.**

## REAGENT PREPARATION

Allow the kit reagents to come to room temperature for 30 minutes. Ensure that all samples have reached room temperature and have been diluted as appropriate prior to running them in the kit.

### Assay Buffer *now supplied as a concentrate*

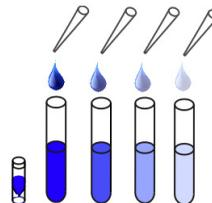
Dilute Assay Buffer Concentrate 1:5 by adding one part of the concentrate to four parts of deionized water. Once diluted this is stable for 3 months at 4°C.

### Wash Buffer

Dilute Wash Buffer Concentrate 1:20 by adding one part of the concentrate to nineteen parts of deionized water. Once diluted this is stable at room temperature for 3 months.

### Standard Preparation

Label tubes #1 through #5. Briefly spin vial of standard in a microcentrifuge to ensure contents are at bottom of vial. Pipet 475  $\mu\text{L}$  of Assay Buffer into tube #1 and 300  $\mu\text{L}$  into tubes #2 to #5. Carefully add 25  $\mu\text{L}$  of the RBP stock solution to tube #1 and vortex completely. Take 100  $\mu\text{L}$  of the RBP solution in tube #1 and add it to tube #2 and vortex completely. Repeat these serial dilutions for tubes #3 through #5. The concentration of RBP in tubes 1 through 5 will be 1,000, 250, 62.5, 15.625, and 3.906 ng/mL.



**Use all Standards within 2 hours of preparation.**

|   | Std 1 | Std 2 | Std 3 | Std 4  | Std 5 |
|---|-------|-------|-------|--------|-------|
| <b>Assay Buffer Volume (<math>\mu\text{L}</math>)</b> | 475   | 300   | 300   | 300    | 300   |
| <b>Addition</b>                                       | Stock | Std 1 | Std 2 | Std 3  | Std 4 |
| <b>Volume of Addition (<math>\mu\text{L}</math>)</b>  | 25    | 100   | 100   | 100    | 100   |
| <b>Final Conc (ng/mL)</b>                             | 1,000 | 250   | 62.5  | 15.625 | 3.906 |



## ASSAY PROTOCOL

**We recommend that all standards and samples be run in duplicate to allow the end user to accurately determine RBP concentrations.**

1. Use the plate layout sheet on the back page of the kit insert to aid in proper sample and standard identification. Determine the number of wells to be used and return unused wells to foil pouch with desiccant. Seal the ziploc plate bag and store at 4°C.
2. Pipet 50  $\mu$ L of samples or standards into wells in the plate. Pipet 75  $\mu$ L of Assay Buffer into the non-specific binding (NSB) wells. Pipet 50  $\mu$ L of Assay Buffer into the maximum binding (B0 or Zero standard) wells.
3. Add 25  $\mu$ L of the DetectX<sup>®</sup> RBP Conjugate to each well, using a repeater pipet.
4. Add 25  $\mu$ L of the DetectX<sup>®</sup> RBP Antibody to each well, **except the NSB wells**, using a repeater pipet.
5. Gently tap the sides of the plate to ensure adequate mixing of the reagents. Cover the plate with the plate sealer and shake at room temperature for 1 hour.
6. Aspirate the plate and wash each well 4 times with 300  $\mu$ L wash buffer. Tap the plate dry on clean absorbent towels.
7. Add 100  $\mu$ L of the TMB Substrate to each well, using a repeater pipet.
8. Incubate the plate at room temperature for 30 minutes without shaking.
9. Add 50  $\mu$ L of the Stop Solution to each well, using a repeater pipet.
10. Read the optical density generated from each well in a plate reader capable of reading at 450 nm.
11. Use the plate reader's built-in 4PLC software capabilities to calculate RBP concentration for each sample.

*NOTE: If you are using only part of a strip well plate, at the end of the assay throw away the used wells and retain the plate frame for use with the remaining unused wells.*

## CALCULATION OF RESULTS

Average the duplicate OD readings for each standard and sample. Create a standard curve by reducing the data using the 4PLC fitting routine on the plate reader, after subtracting the mean OD's for the NSB. The sample concentrations obtained, calculated from the %B/B0 curve, should be multiplied by the dilution factor to obtain neat sample values. Urine RBP values should be normalized to creatinine levels by running the same urines in the DetectX<sup>®</sup> Urinary Creatinine Detection Kit, K002-H1.

Or use the online tool from MyAssays to calculate the data:

[www.myassays.com/arbor-assays-retinol-binding-protein-urinary-eia-kit.assay](http://www.myassays.com/arbor-assays-retinol-binding-protein-urinary-eia-kit.assay)



### TYPICAL DATA

| Sample     | Mean OD | Net OD | % B/B0 | RBP Conc. (ng/mL) |
|------------|---------|--------|--------|-------------------|
| NSB        | 0.098   | 0      | -      | -                 |
| Standard 1 | 0.194   | 0.096  | 10.7   | 1,000             |
| Standard 2 | 0.319   | 0.221  | 24.5   | 250               |
| Standard 3 | 0.516   | 0.418  | 46.4   | 62.5              |
| Standard 4 | 0.740   | 0.642  | 71.3   | 15.625            |
| Standard 5 | 0.907   | 0.809  | 89.8   | 3.906             |
| B0         | 0.999   | 0.901  | 100    | 0                 |
| Sample 1   | 0.318   | 0.220  | 24.4   | 251.1             |
| Sample 2   | 0.598   | 0.500  | 55.4   | 38.25             |

**Always run your own standard curve for calculation of results. Do not use this data.**

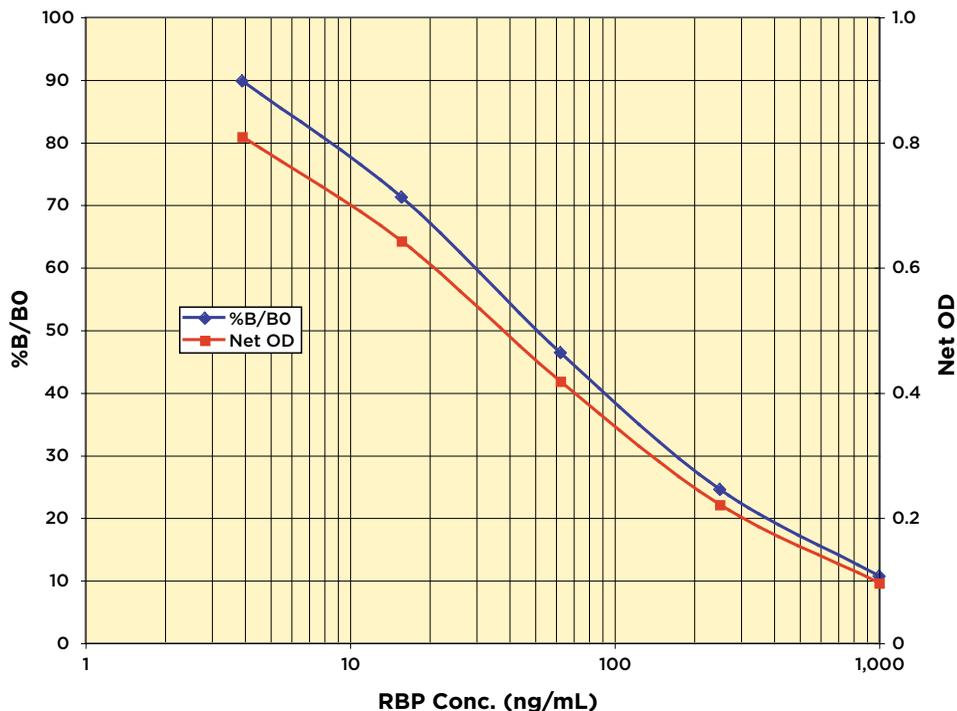
**Conversion Factor: 1 ng/mL of human RBP is equivalent to 47.62 pM RBP.**



\*The MyAssays logo is a registered trademark of MyAssays Ltd.

ARBOR  
ASSAYS

## Typical Standard Curve



Always run your own standard curves for calculation of results. Do not use this data.

## VALIDATION DATA

### Sensitivity and Limit of Detection

Sensitivity was calculated by comparing the OD's for twenty wells run for each of the B0 and standard #5. The detection limit was determined at two (2) standard deviations from the B0 along the standard curve.

**Sensitivity was determined as 2.90 ng/mL.**

The Limit of Detection for the assay was determined in a similar manner by comparing the OD's for twenty replicates for each of the zero standard and a low concentration human urine sample.

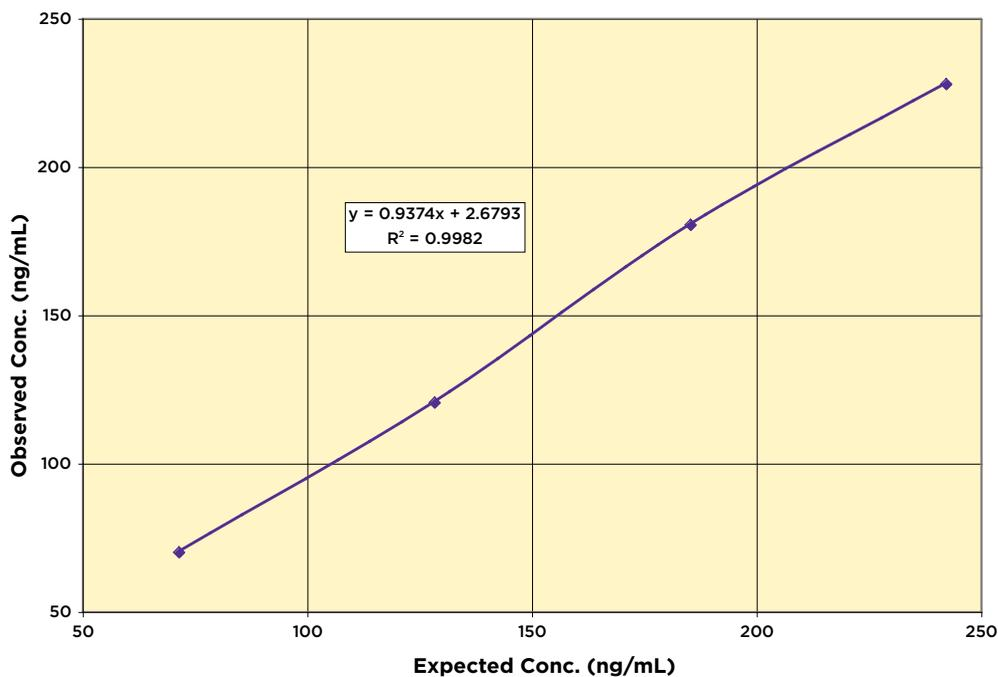
**Limit of Detection was determined as 4.09 ng/mL\*.**

\* Note: Due to the dilute nature of this sample it was run neat instead of being diluted 1:2.

## Linearity

Linearity was determined by taking two human urine samples diluted 1:2, one with a low diluted RBP level of 14.5 ng/mL and one with a higher diluted level of 299.2 ng/mL, and mixing them in the ratios given below. The measured concentrations were compared to the expected values based on the ratios used.

| High Urine           | Low Urine | Expected Conc. (ng/mL) | Observed Conc. (ng/mL) | % Recovery   |
|----------------------|-----------|------------------------|------------------------|--------------|
| 80%                  | 20%       | 242.2                  | 227.9                  | 94.1         |
| 60%                  | 40%       | 185.3                  | 180.4                  | 97.3         |
| 40%                  | 60%       | 128.4                  | 120.5                  | 93.9         |
| 20%                  | 80%       | 71.5                   | 70.0                   | 98.0         |
| <b>Mean Recovery</b> |           |                        |                        | <b>95.8%</b> |



### Intra Assay Precision

Four human urine samples were diluted 1:2 with Assay Buffer and run in replicates of 8 in an assay. The mean and precision of the calculated RBP concentrations were:

| Sample | RBP Conc. (ng/mL) | %CV |
|--------|-------------------|-----|
| 1      | 14.8              | 4.5 |
| 2      | 28.0              | 7.5 |
| 3      | 53.5              | 7.3 |
| 4      | 323.7             | 2.1 |

### Inter Assay Precision

Four human urine samples were diluted 1:2 with Assay Buffer and run in duplicates in twenty-one assays run over multiple days by three operators. The mean and precision of the calculated RBP concentrations were:

| Sample | RBP Conc. (ng/mL) | %CV |
|--------|-------------------|-----|
| 1      | 14.4              | 9.1 |
| 2      | 28.0              | 9.2 |
| 3      | 50.8              | 8.1 |
| 4      | 309.2             | 14  |

## SAMPLE VALUES

Fourteen random human urine samples were tested in the assay. Values ranged from 6.8 to 788.5 ng/mL with a mean of 114.7 ng/mL. These samples were also run in the DetectX® Urinary Creatinine Detection Kit, K002-H1, and the RBP levels normalized to creatinine levels. Normalized values ranged from 35.9 to 573.9 µg RBP/g creatinine.

Normal ranges for urinary RBP are < 130 µg RBP/g creatinine for individuals under 50 years of age and < 172 µg RBP/g creatinine for those equal to or older than 50<sup>15</sup>.

Five random human whole blood DBS samples were punched out as 1/4" or 1/8" and tested in the assay. Values ranged from 7.54 to 19.7 µg/mL.

### Other Species

We have tested a range of urines from other species for RBP. These include rat, dog and rhesus monkey urine. Because of the difficulty in obtaining urine from animals with known medical history, the values we obtained may not be representative of normal or diseased states. Rat urine diluted 1:2 with Assay Buffer had a neat sample value of 43.37 ng/mL and when normalized to urinary creatine gave a reading of 176.1 µg/g creatinine.

Samples of urine from healthy dog and rhesus monkey read below the 3.906 ng/mL standard. We therefore concentrated these samples by freeze drying them and reconstituting them in one-tenth their original volume with Assay Buffer. At that concentration, neat dog urine read at 29.95 ng/mL and when normalized to urinary creatine gave a reading of 32.27 µg/g creatinine. The neat monkey urine read at 10.50 ng/mL and when normalized to urinary creatine gave a reading of 395.5 µg/g creatinine.

15. 84447 Overview: Retinol-Binding Protein, Random, Urine." Mayo Medical Laboratories: Reference Laboratory services for hospitals worldwide. Web. 02 Sept. 2009.  
[www.mayomedicallaboratories.com/test-catalog/Overview/84447](http://www.mayomedicallaboratories.com/test-catalog/Overview/84447)



## LIMITED WARRANTY

Arbor Assays warrants that at the time of shipment this product is free from defects in materials and workmanship. This warranty is in lieu of any other warranty expressed or implied, including but not limited to, any implied warranty of merchantability or fitness for a particular purpose.

We must be notified of any breach of this warranty within 48 hours of receipt of the product. No claim shall be honored if we are not notified within this time period, or if the product has been stored in any way other than outlined in this publication. The sole and exclusive remedy of the customer for any liability based upon this warranty is limited to the replacement of the product, or refund of the invoice price of the goods.

## CONTACT INFORMATION

For details concerning this kit or to order any of our products please contact us:

### Arbor Assays

1514 Eisenhower Place  
Ann Arbor, Michigan 48108 USA

Phone: 734-677-1774

Fax: 734-677-6860

Web: [www.ArborAssays.com](http://www.ArborAssays.com)

### E Mail Addresses:

[Info@ArborAssays.com](mailto:Info@ArborAssays.com)

[Orders@ArborAssays.com](mailto:Orders@ArborAssays.com)

[Technical@ArborAssays.com](mailto:Technical@ArborAssays.com)

[Contracts@ArborAssays.com](mailto:Contracts@ArborAssays.com)



## OFFICIAL SUPPLIER TO ISWE

Arbor Assays and the International Society of Wildlife Endocrinology (ISWE) signed an exclusive agreement for Arbor Assays to supply ISWE members with EIA kits for wildlife conservation research.

*DetectX<sup>®</sup>, ThioStar<sup>®</sup> and the Arbor Assays logo are all registered trademarks.*



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



Printed on Forest Stewardship Council certified paper