

distributed in the US/Canada by:  
**EAGLE BIOSCIENCES, INC.**  
 20A NW Blvd, Suite 112 Nashua, NH 03063  
 Phone: 617-419-2019 FAX: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



# KRIBIOLISA™ Bovine Serum Albumin ELISA

**REF** : KBBP10

Ver 2.1

**RUO**

Enzyme Immunoassay for the Quantitative Determination of Bovine Serum Albumin in serum, plasma and other biological samples.

|            |                              |            |                                       |
|------------|------------------------------|------------|---------------------------------------|
| <b>RUO</b> | <b>For Research Use Only</b> | <b>REF</b> | <b>Catalog Number</b>                 |
|            | <b>Store At</b>              | <b>LOT</b> | <b>Batch Code</b>                     |
|            | <b>Manufactured By</b>       |            | <b>Biological Risk</b>                |
|            | <b>Expiry Date</b>           |            | <b>Consult Operating Instructions</b> |

*For Research Use Only. Purchase does not include or carry the right to resell or transfer this product either as a stand-alone product or as a component of another product. Any use of this product other than the permitted use without the express written authorization of KRISHGEN BioSystems is strictly prohibited.*

**REF** KBBP10

96 tests

**KRISHGEN BioSystems** For US/Europe Customers: toll free +1(888)-970-0827 | tel +1(562)-568-5005  
 For Asia/India Customers: +91(22)-49198700  
 Email: [sales@krishgen.com](mailto:sales@krishgen.com) | <http://www.krishgen.com>

**Introduction:**

The KRIBIOLISA™ ELISA kits are used for assessing the specific biomarker in samples analytes which may be serum, plasma and cell culture supernatant as validated with the kit. The kit employs a competitive ELISA technique.

**Intended Use:**

The Bovine Serum Albumin KRIBIOLISA™ ELISA kit is used as an analytical tool for quantitative determination of Bovine Serum Albumin in serum, plasma and other biological samples.

**Principle:**

The method employs competitive enzyme-linked immunosorbent assay (ELISA) technique to assay the level of Bovine Serum Albumin in samples. Standards or Samples competes with the biotinylated Bovine Serum Albumin, to form a complex with the Bovine Serum Albumin antibody coated microtiter well. Wells are washed to remove the excess conjugate and Streptavidin:HRP Conjugate is added to the microplate and incubated. After incubation and a washing step TMB Substrate, are added. Blue color develops on incubation and the reaction is stopped with a Stop Solution to form a yellow color. The concentration of the Bovine Serum Albumin in the samples is inversely proportional to the yellow color developed (absorbance) in the wells

**Materials Provided:**

1. Bovine Serum Albumin Antibody Coated Microtiter Plate (8 x 12 wells) - 1 no
2. Bovine Serum Albumin Standard (lyophilized, concentrated, 30000 ng/ml) - 2 vials
3. Biotinylated Bovine Serum Albumin (concentrated) – 120 ul
4. Streptavidin:HRP Conjugate (concentrated) - 120 ul
5. Standard Diluent - 20 ml
6. Biotin Antigen Dilution Buffer - 12 ml
7. HRP Conjugate Dilution Buffer - 12 ml
8. (20X) Wash Buffer - 25 ml
9. TMB Substrate - 12 ml
10. Stop Solution - 12 ml
11. Instruction Manual

**Materials to be provided by the End-User:**

1. Microtiter Plate Reader able to measure absorbance at 450 nm.
2. Adjustable pipettes and multichannel pipettor to measure volumes ranging from 25 ul to 1000 ul
3. Deionized (DI) water
4. Wash bottle or automated microplate washer
5. Clean tubes and Eppendorf tubes
6. Precision single and multi-channel pipette and disposable tips.
7. 37°C incubator
8. Timer.

**Handling/Storage:**

1. All reagents should be stored as indicated on the component label.
2. All the reagents and wash solutions should be used within 12 months from manufacturing date.
3. Before using, bring all components to room temperature (18-25°C). Upon assay completion ensure all components of the kit are returned to appropriate storage conditions.
4. The Substrate is light-sensitive and should be protected from direct sunlight or UV sources.

**Health Hazard Warnings:**

1. Reagents that contain preservatives may be harmful if ingested, inhaled or absorbed through the skin.
2. For Research Use Only.

**Sample Preparation and Storage:**

Specimens should be clear and non-hemolyzed. Samples should be run at a number of dilutions to ensure accurate quantitation.

1. Extract as soon as possible after specimen collection as per relevant procedure. The samples should be tested as soon as possible after the extraction. Alternately the extracted samples can be kept in -20°C. Avoid repeated freeze-thaw cycles.
2. **Serum-** Coagulate at room temperature for 10-20 minutes; centrifuge for 20-min at 2000-3000 rpm. Remove the supernatant. If precipitation appears, recentrifuge.
3. **Plasma-** Use EDTA or citrate plasma as an anticoagulant, mix for 10-20 minutes; centrifuge for 15-min at 2000-3000 rpm. Remove the supernatant carefully. If precipitation appears, recentrifuge.
4. **Cell Culture Supernatant-** Collect sample in a sterile container. Centrifuge for 20-mins at 2000-3000 rpm. Remove the supernatant carefully. When examining the components within the cell, dilute cell suspension with PBS (pH 7.2-7.4), if cell concentration is greater than 1 million/ml. Damage the cells by repeated freeze-thaw cycles to release intracellular components. Centrifuge for 20-min at 2000-3000 rpm. If precipitation appears, centrifuge again.
5. **Tissue Samples-** Rinse tissues in PBS (pH 7.4) to remove excess blood thoroughly and weigh before homogenization. Mince tissues and homogenize them in PBS (pH7.4) with a glass homogenizer on ice. Thaw at 2-8°C or freeze at -20°C. Centrifuge at 2000-3000 RPM for approximately 20 minutes and collect the supernatant carefully.

**Note:** Grossly hemolyzed samples are not suitable for use in this assay.

**Reagent Preparation (all reagents should be diluted immediately prior to use):**

1. Label any aliquots made with the kit Lot No and Expiration date and store it at appropriate conditions mentioned.
2. Bring all reagents to Room temperature before use.
3. To make **Wash Buffer (1X) 500 ml**; dilute **25 ml of (20X) Wash Buffer in 475 ml of DI water**.
4. **Streptavidin:HRP Conjugate & Biotinylated Bovine Serum Albumin Working Solution** - Briefly spin or centrifuge the streptavidin:HRP Conjugate & Biotinylated Bovine Serum Albumin before use. Dilute them to the working concentration 100-fold with HRP Conjugate Dilution Buffer & Biotin Antigen Dilution Buffer, respectively.
5. **Standards Preparation:** Reconstitute original Bovine Serum Albumin Standard with 1.0 ml of Standard Diluent. Keep the standard for 10 mins with gentle agitation before making further dilutions. Prepare the additional Standards by serially diluting the standard stock solution as per the below table.

| Standard Concentration | Standard Vial | Dilution Particulars                           |
|------------------------|---------------|--|
| 30000 ng/ml            | Standard No.6 | Reconstitute with 1.0 ml Standard Diluent      |
| 10000 ng/ml            | Standard No.5 | 300 ul Standard No.6 + 600 ul Standard Diluent |
| 3333.3 ng/ml           | Standard No.4 | 300 ul Standard No.5 + 600 ul Standard Diluent |
| 1111.1 ng/ml           | Standard No.3 | 300 ul Standard No.4 + 600 ul Standard Diluent |
| 370.4 ng/ml            | Standard No.2 | 300 ul Standard No.3 + 600 ul Standard Diluent |
| 0 ng/ml                | Standard No.1 | 300 ul Standard Diluent only                   |

**Procedural Notes:**

1. In order to achieve good assay reproducibility and sensitivity, proper washing of the plates to remove excess un-reacted reagents is essential.
2. High Dose Hook Effect may be observed in samples with very high concentrations of Bovine Serum Albumin. High Dose Hook Effect is due to excess of antibody for very high concentrations of Bovine Serum Albumin present in the sample.

3. Bovine Serum Albumin concentration of the undiluted sample is less than the diluted sample, this may be indicative of the Hook Effect.
4. Avoid assay of Samples containing sodium azide ( $\text{NaN}_3$ ), as it could destroy the HRP activity resulting in under-estimation of the amount of Bovine Serum Albumin.
5. It is recommended that all Standards and Samples be assayed in duplicates or triplicates.
6. Maintain a repetitive timing sequence from well to well for all the steps to ensure that the incubation timings are same for each well.
7. If the Substrate has a distinct blue color prior to use it may have been contaminated and use of such substrate can lead to compromise of the sensitivity of the assay.
8. The plates should be read within 30 minutes after adding the Stop Solution.
9. Make a work list in order to identify the location of Standards and Samples.

**Assay Procedure:**

1. It is strongly recommended that all Standards and Samples be run in duplicates or triplicates. A standard curve is required for each assay.
2. Add **50 ul prepared Standards and Samples** to respective wells.
3. Pipette **50 ul Biotinylated Bovine Serum Albumin Antigen Working Solution** to all wells.
4. Cover the plate with a sealer and incubate for 60 minutes at 37°C.
5. Aspirate and wash plate 4 times with diluted Wash Buffer (1X) and blot residual buffer by firmly tapping plate upside down on absorbent paper. Wipe of any liquid from the bottom outside of the microtiter wells as any residue can interfere in the reading step.
6. Pipette **100 ul Streptavidin:HRP Conjugate Working Solution** to all wells. Mix well.
7. Cover the plate with a sealer and incubate for 30 minutes at 37°C.
8. Aspirate and wash as per Step (5) above.
9. Pipette **100 ul TMB Substrate** in all the wells.
10. Incubate the plate at **37°C for 10 minutes**. DO NOT SHAKE or else it may result in higher backgrounds and worse precision.
11. Pipette **100 ul of Stop Solution** to all wells. The wells should turn from blue to yellow in color.
12. Read the absorbance at 450 nm with a microplate within 10-15 minutes after addition of Stop solution.

**Calculation of Results:**

Determine the Mean Absorbance for each set of duplicate or triplicate Standards and Samples. Using Graph Paper, plot the average value (absorbance 450nm) of each standard on the Y-axis versus the corresponding concentration of the standards on the X-axis. Draw the best fit curve through the standard points. To determine the unknown Bovine Serum Albumin concentrations, find the unknown's Mean Absorbance value on the Y-axis and draw a horizontal line to the standard curve. At the point of intersection, draw a vertical line to the X-axis and read the Bovine Serum Albumin Concentration.

If samples were diluted, multiply by the appropriate dilution factor. Software which is able to generate a cubic spline curve-fit is best recommended for automated results.

**Note:**

It is recommended to repeat the assay at a different dilution factor in the following cases:

- If the sample absorbance value is below the first standard.

**Quality Control:**

It is recommended that for each laboratory assay appropriate quality control samples in each run to be used to ensure that all reagents and procedures are correct.

**Performance Characteristics of the Kit:**

This kit has been validated. Please view the details herein below.

**Standard Calibration Range:**

370.4 ng/ml - 30000 ng/ml

**Sensitivity:****Limit Of Quantification:**

It is defined as the lowest detectable concentration that can be determined with an acceptable repeatability and the LOQ was found to be 161.2 ng/ml.

**Specificity:**

This assay has high sensitivity and excellent specificity for detection of Bovine Serum Albumin. No significant cross-reactivity or interference between Bovine Serum Albumin and analogues was observed.

**Recovery**

Matrices listed below were spiked with certain level of Bovine Serum Albumin and the recovery rates were calculated by comparing the measured value to the expected amount of Bovine Serum Albumin in samples.

| Matrix              | Recovery Range (%) | Average (%) |
|---------------------|--------------------|-------------|
| Serum(n=5)          | 85-96              | 90          |
| EDTA Plasma(n=5)    | 86-98              | 92          |
| Heparin Plasma(n=5) | 96-105             | 101         |

**Precision:**

Intra-Assay: CV<10%

Inter-Assay: CV<12%

**Linearity**

The linearity of the kit was assayed by testing samples spiked with appropriate concentration of Bovine Serum Albumin and their serial dilutions. The results were demonstrated by percentage of calculated concentration to the expectation.

| Sample              | 1:2    | 1:4     | 1:8     | 1:16    |
|---------------------|--------|---------|---------|---------|
| Serum(n=5)          | 81-93% | 97-105% | 82-97%  | 84-103% |
| EDTA Plasma(n=5)    | 90-99% | 87-94%  | 95-102% | 87-98%  |
| Heparin Plasma(n=5) | 83-97% | 82-92%  | 80-104% | 85-101% |

**Safety Precautions:**

- **This kit is For Research Use only.** Follow the working instructions carefully.
- The expiration dates stated on the kit are to be observed. The same relates to the stability stated for reagents
- Do not use or mix reagents from different lots.
- Do not use reagents from other manufacturers.
- Avoid time shift during pipetting of reagents.
- All reagents should be kept in the original shipping container.
- Some of the reagents contain small amount of sodium azide (< 0.1 % w/w) as preservative. They must not be swallowed or allowed to come into contact with skin or mucosa.
- Source materials maybe derived from Bovine body fluids or organs used in the preparation of this kit were tested and found negative for HBsAg and HIV as well as for HCV antibodies. However, no known test guarantees the absence of such viral agents. Therefore, handle all components and all patient samples as if potentially hazardous.
- Since the kit contains potentially hazardous materials, the following precautions should be observed
  - Do not smoke, eat or drink while handling kit material
  - Always use protective gloves
  - Never pipette material by mouth



- Wipe up spills promptly, washing the affected surface thoroughly with a decontaminant.
- In any case GLP should be applied with all general and individual regulations to the use of this kit.

## Typical Example of a Work List

| Well #   | Contents      | Absorbance at 450nm | Mean Absorbance | Interpolated Concentration |
|----------|---------------|---------------------|-----------------|----------------------------|
| 1A<br>2A | Standard No.1 |                     |                 |                            |
| 1B<br>2B | Standard No.2 |                     |                 |                            |
| 1C<br>2C | Standard No.3 |                     |                 |                            |
| 1D<br>2D | Standard No.4 |                     |                 |                            |
| 1E<br>2E | Standard No.5 |                     |                 |                            |
| 1F<br>2F | Standard No.6 |                     |                 |                            |
| 1G<br>2G | Sample        |                     |                 |                            |
| 1H<br>2H | Sample        |                     |                 |                            |
| 3A<br>4A | Sample        |                     |                 |                            |
| 3B<br>4B | Sample        |                     |                 |                            |

**LIMITED WARRANTY**

Krishgen Biosystems does not warrant against damages or defects arising in shipping or handling, or out of accident or improper or abnormal use of the Products; against defects in products or components not manufactured by Krishgen Biosystems, or against damages resulting from such non-Krishgen Biosystems made products or components. Krishgen Biosystems passes on to customer the warranty it received (if any) from the maker thereof of such non Krishgen made products or components. This warranty also does not apply to Products to which changes or modifications have been made or attempted by persons other than pursuant to written authorization by Krishgen Biosystems.

THIS WARRANTY IS EXCLUSIVE. The sole and exclusive obligation of Krishgen Biosystems shall be to repair or replace the defective Products in the manner and for the period provided above. Krishgen Biosystems shall not have any other obligation with respect to the Products or any part thereof, whether based on contract, tort, and strict liability or otherwise. Under no circumstances, whether based on this Limited Warranty or otherwise, shall Krishgen Biosystems be liable for incidental, special, or consequential damages.

This Limited Warranty states the entire obligation of Krishgen Biosystems with respect to the Products. If any part of this Limited Warranty is determined to be void or illegal, the remainder shall remain in full force and effect.

Krishgen Biosystems. 2022

**THANK YOU FOR USING KRISHGEN PRODUCT!**

KRISHGEN BIOSYSTEMS®, GENLISA®, DHARMAPLEX™, GENBULK™, KRIBIOLISA™, KRISHZYME®, KRISHGEN®, KRIBIOLISA®, KRISHPLEX®, TITANIUM®, QUALICHEK® are registered trademarks of KRISHGEN BIOSYSTEMS. ©KRISHGEN BIOSYSTEMS. ALL RIGHTS RESERVED.

KRISHGEN BIOSYSTEMS | OUR REAGENTS | YOUR RESEARCH |

## SYMBOLS KEY

|   |                                      |
|---|--------------------------------------|
|    | Coated Microtiter Plate (8x12 wells) |
|    | Standard                             |
|    | Biotinylated Antibody                |
|    | Conjugate Horseradish Peroxidase     |
|    | Biotin Antigen Dilution Buffer       |
|    | HRP Conjugate Dilution Buffer        |
|    | Standard Diluent                     |
|  | (20X) Wash Buffer                    |
|  | TMB Substrate                        |
|  | Stop Solution                        |
|  | Consult Instructions for Use         |
|  | Catalog Number                       |
|  | Expiration Date                      |
|  | Storage Temperature                  |

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) info@eaglebio.com



**EAGLE**  
BIOSCIENCES