

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 info@eaglebio.com



KRIBIOLISA™

Pichia pastoris Host Cell Proteins

ELISA

REF: KBBP02

Ver 4.0

RUO

Immunoassay for quantitative determination of *P. pastoris* Host Cell Proteins

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

For Research Purposes Only. Not for use in diagnostic or therapeutic procedures. 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 KBBP02

96 tests



KRISHGEN BioSystems

For US / Europe: toll free +1(888)-970-0827 tel: +1(562)-568-5005

For Asia / India: tel: +91(22)-49198700

Email: sales@krishgen.com

Introduction:

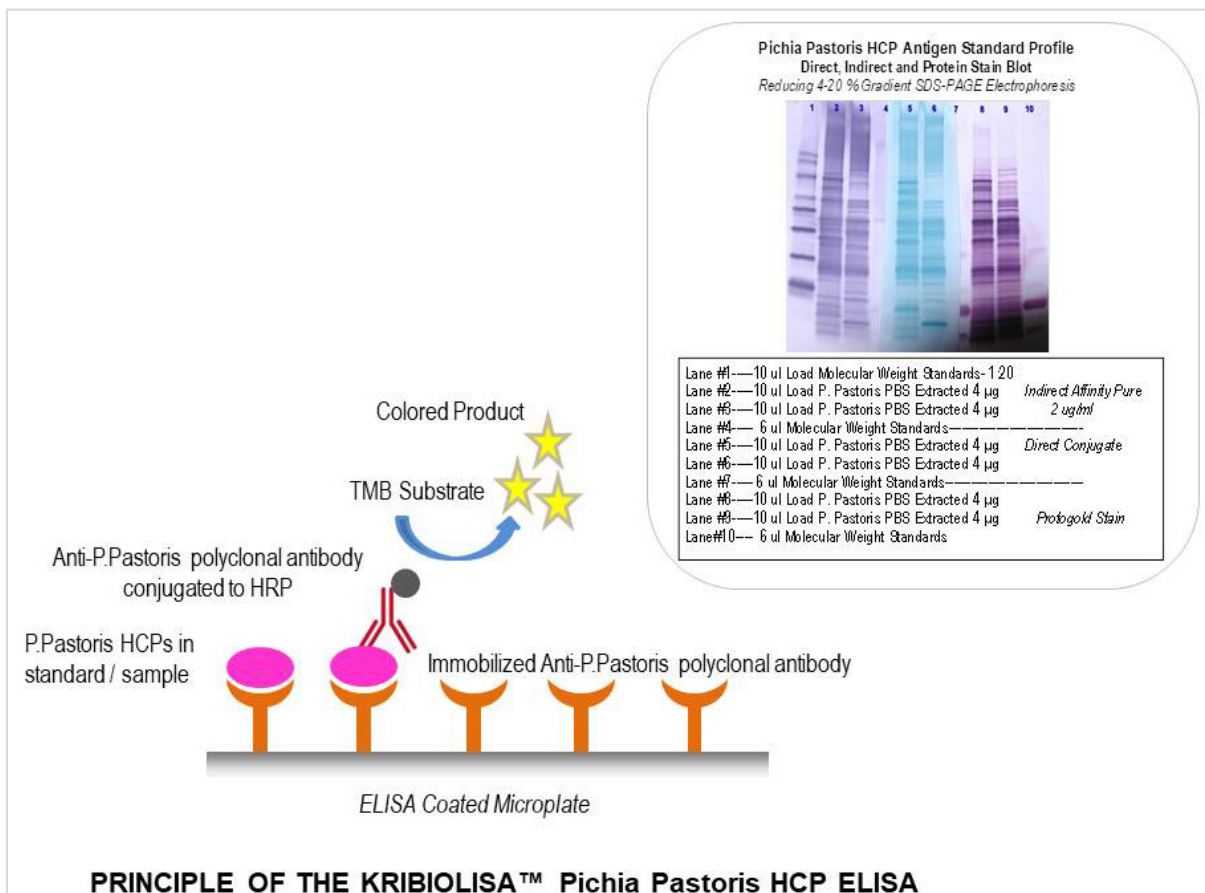
A variety of proteins and pDNA which are used as therapeutic agents in humans and animals are produced through recombinant expression in *P.pastoris*. The manufacturing and purification process of these products tends to leave the potential for contamination by Host Cell Proteins (HCPs) from *P.pastoris* which may result in adverse toxic or immunological reactions that ultimately affect the efficacy of the therapeutic agent. The simple, objective and semi quantitative ELISA is a highly sensitive method that aids in purification process development, process control, quality control and product release testing optimally.

Intended Use:

This KRIBIOLISA *Pichia Pastoris* HCP ELISA is a generic kit intended in determining the presence of *P. pastoris* Host Cell Proteins contamination in various biotechnology products that are manufactured through recombinant expression in *P. pastoris*. The kit has been validated successfully for testing of in process and final product HCPs in variety of products regardless of growth and purification process.

Principle:

This assay is based on the Sandwich ELISA procedure. Samples containing *P.pastoris* HCPs are reacted with already coated affinity purified capture anti-*P.pastoris* antibody and bind to them. Plates are washed with wash buffer to remove unbound reactants. Anti *P.pastoris*-HRP conjugate is added which results in the formation of a sandwich complex of solid phase antibody-HCP-enzyme labeled antibody. The wells are washed to remove any unbound reactants as per the wash procedure. The substrate 3, 3',5, 5' Tetra Methyl Benzidine is then reacted. The amount of hydrolyzed substrate is read on a microtiter plate reader and it is directly proportional to the concentration of *P. pastoris* HCPs present.



Materials Provided:

| Part | Description | Qty |
|-------------------------------------------------|-----------------------------------------------------------------------------------------------------------------------|--------------|
| Anti- <i>P.pastoris</i> Coated Microtiter Plate | 96 well polystyrene microplate (12 strips of 8 wells) coated with Anti- <i>P.pastoris</i> antibody. | 1 x 96 wells |
| <i>P.pastoris</i> HCP Standards (0.5ml/vial) | Recombinant <i>P.pastoris</i> HCP Standard, (0, 15.6, 31.3, 62.5, 125, 250, 500, 1000 ng/ml) | 8 x 0.5ml |
| Anti- <i>P.pastoris</i> HRP Conjugate | Buffered protein base with preservative thiomersol <0.01% | 12 ml |
| (5X) Sample Diluent | Buffered protein base with preservative thiomersol <0.01% | 12 ml |
| (20X) Wash Buffer | 20-fold concentrated solution of buffered surfactant with preservative thiomersol < 0.01%. May turn yellow over time. | 25 ml |
| TMB Substrate | Stabilized Chromogen | 12 ml |
| Stop Solution | 2N Sulfuric Acid | 12 ml |
| Instruction Manual | | 1 no |

Materials to be provided by the End-User:

1. Microplate Reader able to measure absorbance at 450 nm.
2. Adjustable pipettes to measure volumes ranging from 50 ul to 1000 ul.
3. Deionized (DI) water.
4. Wash bottle or automated microplate washer.
5. Graph paper or software for data analysis.
6. Tubes to prepare standard/sample dilutions.
7. Timer.
8. Absorbent paper.

Storage Information:

1. Store main kit components at 2-8°C.
2. All the reagents and wash solution are stable until the expiration date of the kit.
3. Before using, bring all components to room temperature (18-25°C). Upon assay completion return all components to appropriate storage conditions.

Health Hazard Warnings:

Reagents that contain preservatives may be harmful if ingested, inhaled or absorbed through the skin. Refer to the MSDS online for details.

Specimen Collection and Handling:

Specimens should be clear. Samples should be run at a number of dilutions to ensure accurate quantitation. In case additional sample diluent is required, please contact us at +91-22-49198700 | email: sales@krishgen.com to order.

Cell Culture Supernatant: If necessary, centrifuge to remove debris prior to analysis. Samples can be stored at < -20°C. Avoid repeated freeze/thaw cycles.

Procedural Notes:

1. For good assay reproducibility and sensitivity, proper washing of the ELISA plate to remove excess/unbound reagents is essential.

2. If the *P.pastoris* HCP concentration of the undiluted sample is less than the diluted sample, this may be indicative of the Hook Effect. High Dose Hook Effect may be observed in samples with very high concentrations of *P.pastoris* HCP, usually in samples from the initial stages of purifications. To overcome Hook Effect samples to be assayed should be sufficiently diluted with our recommended diluent.
3. Avoid assay of samples containing sodium azide (NaN₃), as it may destroy the HRP activity of the conjugate resulting in the under-estimation of the levels of *P.pastoris* HCP.
4. All Standards and Samples should be assayed at least in duplicates.
5. Maintain a repetitive timing sequence from well to well for all the steps to ensure that the incubation timings are same for each well.
6. If the Substrate has a distinct blue color prior to use it may have been contaminated and use of such substrate can lead to compromisation of the sensitivity of the assay.
7. The plates should be read within 30 minutes after adding the Stop Solution.
8. Make a work list in order to identify the location of Standards and Samples.

Preparation before Use:

Allow samples to reach room temperature prior to assay. Take care to agitate samples gently in order to ensure homogeneity.

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 **(1X) Wash Buffer**; dilute **25 ml of (20X) Wash Buffer** in **475 ml of DI water**.
4. To make **Sample Diluent (1X)**, add **10ml of Sample Diluent (5X)** to **40ml of DI water**. This is the working solution.

Assay Procedure:

1. Bring all reagents to room temperature prior to use. It is strongly recommended that all standards and samples be run in duplicate or triplicate. A standard curve is required for each assay.
2. Pipette **100 ul** of **Standards** or **Samples** into the respective wells.
3. Cover the plate and incubate at 37°C for 2 hours 30 minutes.
4. Aspirate and wash plate 4 times with **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.
5. Add **100 ul** of **Anti-*P.pastoris*:HRP Conjugate** into each well.
6. Cover the plate and incubate at 37°C for 1 hour.
7. Aspirate and wash plate 4 times with **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.
8. Add **100 ul** of **TMB Substrate** in each well.
9. Incubate the plate at 37°C for 15 minutes in dark. DO NOT SHAKE or else it may result in higher backgrounds and worse precision. Positive wells should turn bluish in color.
10. Pipette out **100 ul** of **Stop Solution**. Wells should turn from blue to yellow in color.

11. Read the absorbance at 450 nm with a microplate reader within 30 minutes.

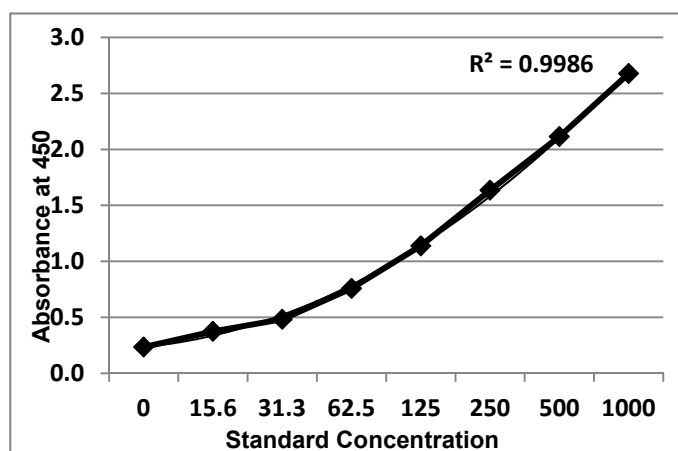
Calculation of Results:

Determine the mean absorbance for each set of duplicate or triplicate standards and samples. Subtract the mean absorbance of the zero standards (background) from each well. Using semi-log graph paper or computer programs, plot the optical densities of each standard on the Y-axis versus the corresponding concentration of the standards on the X-axis. Draw the best fit straight line through the standard points. To determine the unknown *P.pastoris* concentrations, find the unknowns 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 *P.pastoris* concentration. If samples were diluted, multiply by the appropriate dilution factor. Computer based curve-fitting software may be preferred.

Typical Data

| Standard Concentration (ng/ml) | Abs A | Abs B | Mean Abs | Interpolated Concentration | % Interpolated Concentration against Actual Concentration |
|--------------------------------|-------|-------|----------|----------------------------|-----------------------------------------------------------|
| 0 | 0.242 | 0.228 | 0.235 | 1.4 | -- |
| 15.6 | 0.371 | 0.377 | 0.374 | 15.0 | 96.5 |
| 31.3 | 0.490 | 0.475 | 0.482 | 27.1 | 86.6 |
| 62.5 | 0.756 | 0.761 | 0.759 | 63.5 | 101.6 |
| 125 | 1.154 | 1.123 | 1.139 | 129.2 | 103.3 |
| 250 | 1.626 | 1.645 | 1.635 | 257.7 | 103.1 |
| 500 | 2.107 | 2.122 | 2.114 | 470.9 | 94.2 |
| 1000 | 2.743 | 2.612 | 2.678 | 1031.4 | 103.1 |

Typical Graph



Abs = absorbance at 450nm

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 as per EMA/FDA guidelines in line with ICH Code for Harmonization of Biological Assays.

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 15 ng/ml.

Limit Of Detection: It is defined as the lowest detectable concentration corresponding to a signal of Mean of '0' standard plus 2* SD. 10 replicates of '0' standards were evaluated and the LOD was found to be 12 ng/ml

Precision:

Precision is defined as the percent coefficient of variation (%CV) i.e. standard deviation divided by the mean and multiplied by 100. Assay precision was determined by both intra (n=5 assays) and inter assay (n=5 assays) reproducibility on two pools with low (15.6 ng/ml), medium (125 ng/ml) and high (1000 ng/ml) concentrations. While actual precision may vary from laboratory to laboratory and technician to technician, it is recommended that all operators achieve precision below these design goals before reporting results.

| Pool | Intra Assay %CV | Inter Assay %CV |
|--------|-----------------|-----------------|
| Low | <12% | <15% |
| Medium | <8% | <12% |
| High | <8% | <12% |

Recovery/ Interference Studies:

Known amount of *P.pastoris* was spiked into sample diluent (1X). The resulting concentration or “recovery” of the spiked material, demonstrates if the expected value can be measured accurately. If the recovered value differs significantly from the amount expected this may be sign that some factor in the sample matrix may be causing a falsely elevated or falsely depressed value.

It was observed that sample diluent obtained the best recoveries (+/- 20%):

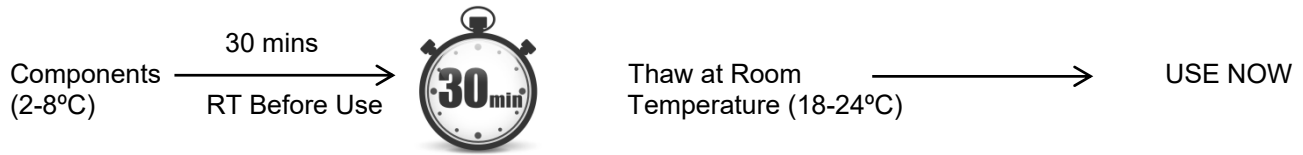
Safety Precautions:

- **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 human 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.

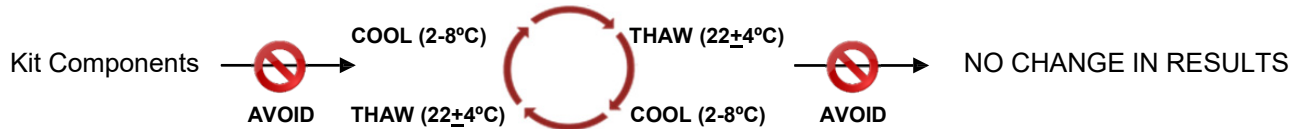


SCHEMATIC ASSAY PROCEDURE

1. Remove all components, 30 minutes before adding into the assay plate.



2. Avoid repeated cool-thaw of the components as there will be a loss of activity and this can affect the results.



3. Pipette **100 ul Standards/Samples** into the respective wells.

4. Cover plate and **incubate** for at 37°C.

5. Aspirate and wash wells 4 times with **(1X) Wash Buffer**.

6. Pipette **100 ul Anti P. Pastoris:HRP Conjugate** into each well.

7. Cover plate and **incubate** for at 37°C.

8. Aspirate and wash wells 4 times with **(1X) Wash Buffer**.

9. Pipette **100 ul TMB Substrate** into each well.

10. Cover plate and **incubate** for at 37°C.

11. Pipette **100 ul Stop Solution** into each well.

12. Read absorbance at 450nm with a microplate reader within of stopping reaction.

Typical Example of a Work List

| Well # | Contents | Absorbance at 450nm | Mean Absorbance | Results |
|--------|------------|---------------------|-----------------|---------|
| 1A | Zero Std | | | |
| 2A | Zero Std | | | |
| 1B | 15.6 ng/ml | | | |
| 2B | 15.6 ng/ml | | | |
| 1C | 31.3 ng/ml | | | |
| 2C | 31.3 ng/ml | | | |
| 1D | 62.5 ng/ml | | | |
| 2D | 62.5 ng/ml | | | |
| 1E | 125 ng/ml | | | |
| 2E | 125 ng/ml | | | |
| 1F | 250 ng/ml | | | |
| 2F | 250 ng/ml | | | |
| 1G | 500 ng/ml | | | |
| 2G | 500 ng/ml | | | |
| 1H | 1000 ng/ml | | | |
| 2H | 1000 ng/ml | | | |
| 1A | Sample | | | |
| 2A | Sample | | | |
| 1B | Sample | | | |
| 2B | Sample | | | |
| 1C | Sample | | | |
| 2C | Sample | | | |
| 1D | Sample | | | |
| 2D | 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 product; 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 product 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 product 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, 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 product. 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. 2021

THANK YOU FOR USING KRISHGEN PRODUCT!

KRISHGEN BIOSYSTEMS®, GENLISA®, DHARMAPLEX™, GENBULK™, GENLISA™, 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

| | |
|-------------------------------------------------------------------------------------|-------------------------------------------------------|
|  | Anti- <i>P.pastoris</i> Microtiter Plate (12x8 wells) |
|  | <i>P.pastoris</i> Standard |
|  | <i>P.pastoris</i> :HRP Conjugate |
|  | (5X) Sample Diluent |
|  | (20X) Wash Buffer |
|  | TMB Substrate |
|  | Stop Solution |
|  | Consult Instructions for Use |
|  | Catalogue Number |
|  | Expiration Date |
|  | Storage Temperature |