



EAGLE
BIOSCIENCES

Prolactin ELISA Kit

Catalog Number:
PRL31-K01 (1 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.

v. 10.0 (effective 23/JAN/2023)

EAGLE BIOSCIENCES, INC.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

WWW.EAGLEBIO.COM

INTENDED USE

The Eagle Biosciences Prolactin ELISA Kit (enzyme-linked immunoassay kit) is intended for the quantitative measurement of prolactin in human serum. The Eagle Biosciences Prolactin ELISA Kit is for research use only and not to be used in diagnostic procedures.

For further information about this kit, its application, or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.

LIMITATIONS RELATED TO INTENDED USE

1. This test is not intended to be used for screening purposes.
2. This test is not intended for home testing or self-testing.
3. The kit is calibrated for the determination of prolactin in human serum. The kit is not calibrated for the determination of prolactin in other specimens of human or animal origin.
4. The results obtained with this kit shall never be used as the basis for a clinical diagnosis or for therapeutic decisions.
5. Although common interfering substances have been evaluated with this test, other substances that have not been evaluated such as drugs and the occurrence of heterophilic antibodies in individuals regularly exposed to animals or animal products have the potential of causing interferences.
6. Some individuals may have antibodies to mouse protein that can possibly interfere in this assay. Therefore, the results from any individuals who have received a preparation of mouse antibodies should be interpreted with caution.
7. The measurement of prolactin may also include the measurement of its other structural forms (big prolactin, macroprolactin, etc.) As a result, individuals exhibiting elevated prolactin levels may require further investigations.

ASSAY BACKGROUND

Prolactin is a polypeptide hormone synthesized by the lactotrophic cells of the anterior pituitary gland. Structurally, it is similar to two other polypeptide hormones namely, growth hormone and placental lactogen. Prolactin is a polypeptide containing 199 amino acids, while growth hormone and placental lactogen each have 191 amino acids. There is approximately 100 µg of prolactin in the human pituitary gland, which is a very small amount when compared to growth hormone, which is present at 8-10 mg.

The target organ of prolactin is the breast (mammary gland). Its main physiological action is not only to initiate but also to sustain lactation. The hypothalamus secretes dopamine, which has a direct effect of inhibition of the secretion of prolactin.

If dopamine is not available or absent the secretion of prolactin is autonomous.

- If the pituitary gland is deficient it leads to failure of lactation.
- In Sheehan's syndrome the pituitary gland is deficient, therefore the prolactin level is reduced.
- A few conditions where increases in prolactin levels are found include hyperprolactinemia, sleep, pregnancy, hypothyroidism and stress.

PRINCIPLE OF THE ASSAY

This Eagle Biosciences Prolactin ELISA Kit is a one-step capture or 'sandwich' type immunoassay. The assay makes use of two highly specific monoclonal antibodies: A monoclonal antibody specific for prolactin is immobilized onto the microplate and another monoclonal antibody specific for a different epitope of prolactin is conjugated to horse radish peroxidase (HRP conjugate).

In the first incubation step, prolactin present in the specimen samples, standards and controls is simultaneously bound by the immobilized antibody and the HRP conjugate antibody, thus forming a

sandwich complex. Excess and unbound materials are removed by a washing step. Next, the TMB substrate (enzyme substrate) is added which reacts with HRP to form a blue coloured product that is directly proportional to the amount of prolactin present. The enzymatic reaction is terminated by the addition of the stopping solution, converting the color from blue to yellow. The absorbance is measured on a microplate reader at 450 nm. A set of standards is used to plot a standard curve from which the amount of prolactin in specimen samples and controls can be directly read.

PROCEDURAL WARNINGS AND PRECAUTIONS

- This kit is for use by trained laboratory personnel (professional use only). For research use only.
- Practice good laboratory practices when handling kit reagents and specimens. This includes:
- Do not pipette by mouth.
- Do not smoke, drink, or eat in areas where specimens or kit reagents are handled.
- Wear protective clothing and disposable gloves.
- Wash hands thoroughly after performing the test.
- Avoid contact with eyes, use safety glasses; in case of contact with eyes, flush eyes with water immediately and contact a doctor.
- Users should have a thorough understanding of this protocol for the successful use of this kit. Reliable performance will only be attained by strict and careful adherence to the instructions provided.
- Do not use this kit beyond the expiry date stated on the label.
- If the kit reagents are visibly damaged, do not use the test kit.
- Do not use kit components from different kit lots within a test and do not use any component beyond the expiration date printed on the label.
- All kit reagents and specimens must be brought to room temperature and mixed gently but thoroughly before use. Avoid repeated freezing and thawing of specimens.
- When the use of water is specified for dilution or reconstitution, use deionized or distilled water.
- Immediately after use, each individual component of the kit must be returned to the recommended storage temperature stated on the label.
- A standard curve must be established for every run.
- It is recommended to all customers to prepare their own control materials or sample pools which should be included in every run at a high and low level for assessing the reliability of results.
- The controls (if applicable with this kit) must be included in every run and their results must fall within the ranges stated in the quality control certificate; a failed control result might indicate improper reagent storage.
- When dispensing the substrate and stopping solutions, do not use pipettes in which these liquids will come into contact with any metal parts.
- The TMB Substrate is sensitive to light and should remain colorless if properly stored. Instability or contamination may be indicated by the development of a blue color, in which case it should not be used.
- Do not use grossly hemolyzed, grossly lipemic, icteric or improperly stored serum.
- Samples or controls containing azide or thimerosal are not compatible with this kit, they may lead to false results.
- Avoid microbial contamination of reagents.
- To prevent the contamination of reagents, use a new disposable pipette tip for dispensing each reagent, sample, standard, and control.
- To prevent contamination of reagents, do not pour reagents back into the original containers.
- Kit reagents must be regarded as hazardous waste and disposed of according to local and/or national regulations.

- Consumables used with the kit that are potentially biohazardous (e.g., pipette tips, bottles or containers containing human materials) must be handled according to biosafety practices to minimize the risk of infection and disposed of according to local and/or national regulations relating to biohazardous waste.
- This kit contains 1 M sulfuric acid in the stopping solution component. Do not combine acid with waste material containing sodium azide or sodium hypochlorite.
- The use of safety glasses, and disposable plastic, is strongly recommended when manipulating biohazardous or bio-contaminated solutions.
- Proper calibration of the equipment used with the test, such as the pipettes and absorbance microplate reader, is required.
- If a microplate shaker is required for the assay procedure, the type and speed of shaker required is stated in the REAGENTS AND EQUIPMENT NEEDED BUT NOT PROVIDED section. Both the type and speed of shaker used can influence the optical densities and test results. If a different type of shake and/or speed is used, the user is responsible for validating the performance of the kit.
- Do not reuse the microplate wells, they are for SINGLE USE only.
- To avoid condensation within the microplate wells in humid environments, do not open the pouch containing the microplate until it has reached room temperature.
- Any serious incident that has occurred in relation to the device shall be reported to the manufacturer and the competent authority of the Member State in which the user and/or the participant is established.
- When reading the microplate, the presence of bubbles in the wells will affect the optical densities (ODs). Carefully remove any bubbles before performing the reading step.

SAFETY CAUTIONS AND WARNINGS

BIOHAZARDS

The reagents should be considered a potential biohazard and handled with the same precautions applied to human specimens. All human specimens should be considered a potential biohazard and handled as if capable of transmitting infections and in accordance with good laboratory practices.

CHEMICAL HAZARDS

Avoid direct contact with any of the kit reagents. Specifically avoid contact with the TMB Substrate (contains tetramethylbenzidine) and Stopping Solution (contains sulfuric acid). If contacted with any of these reagents, wash with plenty of water and refer to SDS for additional information.

SPECIMEN COLLECTION, STORAGE, AND PRE-TREATMENT

Specimen Collection & Storage

Approximately 0.1 mL of serum is required per duplicate determination. Collect 4-5 mL of venous blood into an appropriately labeled tube and allow it to clot. Centrifuge at room temperature and carefully transfer the serum into a new storage tube or container. Serum samples may be stored at 2-8°C for up to 24 hours or at -10°C or lower if the analyses are to be done at a later date.

Specimen Pre-Treatment

Specimen pre-treatment is not required.

REAGENTS AND EQUIPMENT NEEDED BUT NOT PROVIDED

- Calibrated single-channel pipette to dispense 25 µL.
- Calibrated multi-channel pipette to dispense 50 µL, 100 µL, and 150 µL
- Calibrated multi-channel pipettes to dispense 300 µL (if washing manually)

- Automatic microplate washer (recommended)
- Microplate shaker:
 - Orbital shaker (3 mm diameter) set to 600 rpm or
 - Reciprocating shaker (1.5" stroke length) set to 180 oscillations/minute
- Disposable pipette tips
- Distilled or deionized water
- Calibrated absorbance microplate reader with a 450 nm and an upper OD limit of 3.0 or greater

REAGENTS PROVIDED

1. Microplate

Contents: One anti-prolactin monoclonal antibody-coated 96-well (12x8) microplate in a resealable pouch with desiccant.

Format: Ready to Use

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

2. HRP-Conjugate (51x)

Contents: One bottle containing anti-prolactin monoclonal antibody-Horse Radish Peroxidase (HRP) conjugate in a protein-based buffer with a non-mercury preservative.

Format: Concentrated; Requires Preparation

Volume: 0.3 mL/bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

Preparation: **Dilute 1:51** in assay buffer before use (e.g., 40 µL of conjugate concentrate in 2 mL of assay buffer). If the whole plate is to be used dilute 240 µL of conjugate concentrate in 12 mL of assay buffer. Discard any that is left over.

3. Standard A - F

Contents: Six bottles of standard containing specified prolactin concentrations. Protein-based buffer with a non-mercury preservative. Prepared by spiking buffer with defined quantities of prolactin.

Listed below are approximate concentration, please refer to vial labels for exact concentrations.

Concentrations:

0, 20, 100, 400, 800, 3200 µIU/mL

0, 0.77, 3.84, 15.38, 30.75, 123 ng/mL

Format: Ready to Use

Volume: Standard A: 2.0 mL/bottle

Standard B-F: 0.3 mL/bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label.

After Opening: Stable for four weeks.

4. Control 1 - 2

Contents: Two bottles of control containing different prolactin concentrations. Protein-based buffer with a non-mercury preservative. Prepared by spiking buffer with defined quantities of prolactin.
Refer to the QC certificate for the target values and acceptable ranges.

Format: Ready to Use

Volume: 0.3 mL/bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label.
After Opening: Stable for four weeks.

5. **Assay Buffer**

Contents: One bottle containing a protein-based buffer with a non-mercury preservative.

Format: Ready to Use

Volume: 15 mL/bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label.
After Opening: Stable for four weeks.

6. **TMB Substrate**

Contents: One bottle containing a tetramethylbenzidine and hydrogen peroxide in a non-DMF or DMSO containing buffer.

Format: Ready to Use

Volume: 16 mL/bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label.
After Opening: Stable for four weeks.

7. **Stopping Solution**

Contents: One bottle containing 1M sulfuric acid.

Format: Ready to Use

Volume: 6 mL/bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label.
After Opening: Stable for four weeks

8. **Wash Buffer Concentrate (10x)**

Contents: One bottle containing buffer with a non-ionic detergent and a non-mercury preservative

Format: Concentrated; Requires Preparation

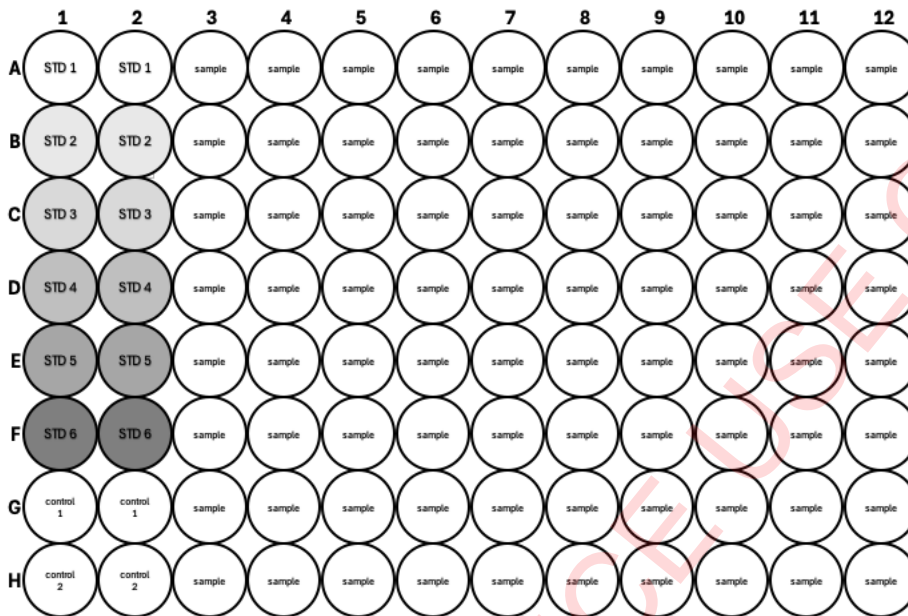
Volume: 50 mL/ bottle

Storage: 2-8°C

Stability: Unopened: Stable until the expiry date printed on the label.
After Opening: Stable for four weeks. Following Preparation: The wash buffer working solution is stable for 2 weeks following preparation, assuming Good Laboratory Practices are adhered to. To prevent microbial growth, prepare the wash buffer working solution in a clean container and store under refrigerated conditions (2-8°C) when not in use.

Preparation: **Dilute 1:10** in distilled or deionized water before use. If the whole microplate is to be used dilute 50 mL of the wash buffer concentrate in 450 mL of distilled or deionized water.

RECOMMENDED ASSAY LAYOUT*



*Layout subject to change based on standard and control quantities

ASSAY PROCEDURE

All kit components, controls, and specimen samples must reach room temperature prior to use. Standards, controls, and specimen samples should be assayed in duplicate. Once the procedure has been started, all steps should be completed without interruption.

1. After all kit components have reached room temperature, mix gently by inversion.
2. Prepare the HRP Conjugate Working Solution and Wash Buffer Working Solution (see Reagents Provided).
3. Plan the microplate wells to be used for standards, controls, and samples. (See Recommended Assay Layout). Remove the strips from the microplate frame that will not be used and place them in the bag with desiccant. Reseal the bag with the unused strips and return it to the refrigerator.
4. **Pipette 25 µL** of each standard, control and specimen sample into assigned wells.
5. **Pipette 100 µL** of the HRP Conjugate Working solution into each well (the use of a multichannel pipette is recommended.)
6. **Incubate** the microplate on a microplate shaker** for **60 minutes** at room temperature.
7. Wash the microplate wells with an automatic microplate washer (preferred) or manually as stated below.
 - a. Automatic: Using an automatic microplate washer, perform a 3-cycle wash using **300 µL/well** of Wash Buffer Working Solution (3 x 300 µL). One cycle consists of aspirating all wells then filling each well with 300 µL of Wash Buffer Working Solution. After the final wash cycle, aspirate all wells and then tap the microplate firmly against absorbent paper to remove any residual liquid.
 - b. Manually: Using an automatic microplate washer, perform a 3-cycle wash using **300 µL/well** of Wash Buffer Working Solution (3 x 300 µL). One cycle consists of aspirating all

wells by briskly emptying the contents of the wells over a waste container, then pipetting 300 μ L of Wash Buffer Working Solution into each well using a multi-channel pipette. After the final wash cycle, aspirate all wells by briskly emptying the contents over a waste container and then tap the microplate firmly against absorbent paper to remove any residual liquid.

8. **Pipette 150 μ L** of TMB substrate into each well (the use of a multi-channel pipette is recommended).
9. **Incubate** the microplate on a microplate shaker** for **10-15 minutes** at room temperature. Do not cover the microplate.
10. **Pipette 50 μ L** of Stopping Solution into each well (the use of a multi-channel pipette is recommended) in the same order and speed as was used for addition of the TMB Substrate. Gently tap the microplate frame to mix the contents of the wells.
11. Measure the optical density (absorbance) in the microplate wells using an absorbance microplate reader set to 450 nm, within 20 minutes after addition of the Stopping Solution.

** See Reagents and Equipment Needed But Not Provided for microplate shaker options.

CALCULATIONS

1. Calculate the mean optical density of each standard, control, and specimen sample duplicate.
2. Use a 4-parameter or 5-parameter curve fit with immunoassay software to generate a standard curve.
3. The immunoassay software will calculate the concentrations of the controls and specimen samples using the mean optical density values and the standard curve.
4. If a sample reads more than 3200 μ IU/mL and needs to be diluted and retested, then dilute with standard A at a dilution of no more than 1:8. The result obtained should be multiplied by the dilution factor.

QUALITY CONTROL

When assessing the validity of the test results, the following criteria should be evaluated:

1. The standard A mean optical density meets the acceptable range as stated in the QC Certificate.
2. The standard with the highest concentration meets the optical density acceptable range as stated in the QC Certificate.
3. The values obtained for the kit controls are within the acceptable ranges as stated in the QC certificate.
4. The results of any external controls that were used meet the acceptable ranges.

TYPICAL DATA

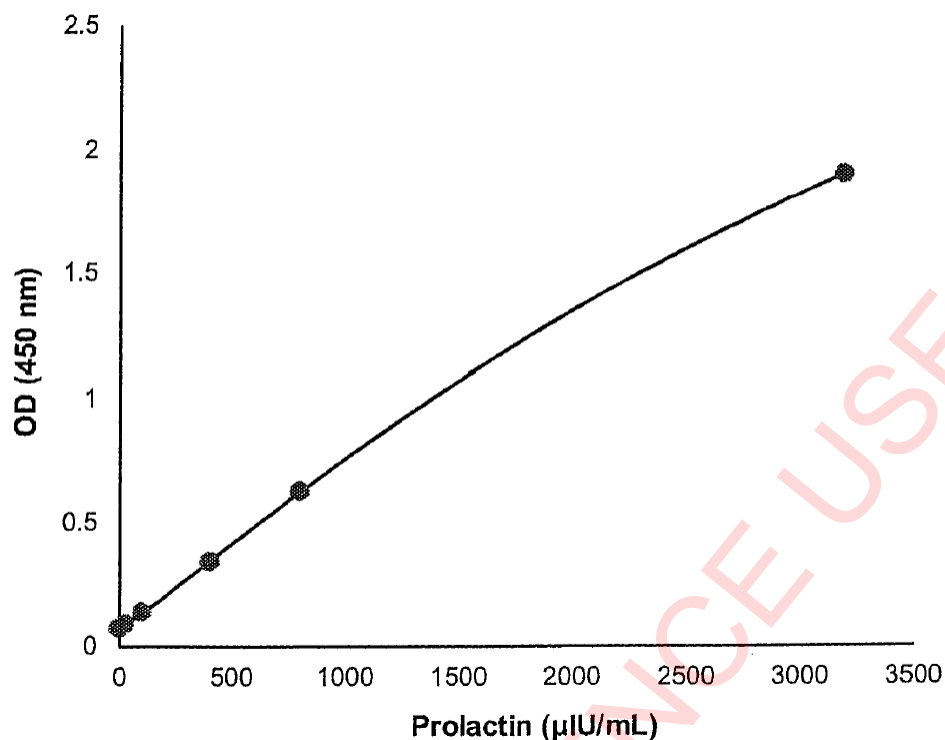
Sample data only. Do not use to calculate results.

Standard	Mean OD (450 nm)	% Binding	Value (μ IU/mL)
A	0.078	3	0
B	0.090	4	20
C	0.137	6	100
D	0.344	15	400
E	0.624	26	800
F	2.361	100	3200
Unknown	0.345	-	402.9



TYPICAL STANDARD CURVE

Sample curve only. Do not use to calculate results.



PERFORMANCE CHARACTERISTICS

Sensitivity

The lower detection limit is calculated from the standard curve by determining the resulting concentration of the mean OD of standard A (based on 10 replicate analyses) plus 2 SD. Therefore, the sensitivity of the Prolactin ELISA Kit is 10 μIU/mL.

Specificity (Cross-Reactivity)

The specificity of the Prolactin ELISA Kit was determined by measuring the apparent prolactin values of the following compounds:

Substance	Concentration Range	Apparent PRL Value (μIU/mL)
hCG (WHO 75/537)	100 – 2500 IU/L	Not Detected
FSH (WHO 1 st 83/575)	25 – 4000 IU/L	Not Detected
hGH (WHO 80/505)	10 – 1000 mg/L	Not Detected
PL	0.1 – 50 mg/L	Not Detected
TSH (WHO 80/558)	25 – 1000 mIU/L	Not Detected

Intra-Assay

Three serum samples were assayed ten time each on the same standard curve. The results (in μIU/mL) are tabulated below:

Sample	Mean	SD	CV%
1	202	14	6.9
2	586	68	11.6

3	1320	136	10.3
---	------	-----	------

Inter-Assay Precision

Three serum samples were assayed ten times over a period of four weeks. The results (in $\mu\text{IU/mL}$) are tabulated below:

Sample	Mean	SD	CV%
1	237	18	7.6
2	589	85	14.4
3	1725	277	13.2

Linearity

Three serum samples were diluted with standard A. The results (in $\mu\text{IU/mL}$) are tabulated below:

Sample	Observed Result	Expected Result	Recovery %
1	292	-	-
1:2	124	146	84.9
1:4	87	73	119.2
1:8	41	37	110.8
2	444	-	-
1:2	202	222	91.0
1:4	116	111	104.5
1:8	69	56	123.2
3	1965	-	-
1:2	1014	983	103.2
1:4	427	491	87.0
1:8	209	246	85.0

Recovery

Spiked samples were prepared by adding defined amounts of prolactin to three serum samples. The results (in $\mu\text{IU/mL}$) are tabulated below:

Sample	Observed Result	Expected Result	Recovery %
1 Unspiked	55	-	-
+ 62	95	117	81.2
+ 130	172	185	93.0
+542	508	597	85.1
2 Unspiked	59	-	-
+ 49	127	108	117.6
+ 145	258	204	126.5
+775	792	834	95.0
3 Unspiked	707	-	-
+ 145	807	852	94.7
+ 385	1356	1092	124.2
+ 775	1868	1482	126.0

REFERENCE RANGES


As for all assays each laboratory should collect data and establish their own range of expected normal values.

Group	Absolute Range ($\mu\text{IU/mL}$)
Males	67 – 360
Females	55 – 2500
Postmenopausal	< 400

LITERATURE

1. Forsyth IA, Myres RP. Human Prolactin. Evidence Obtained by the Bioassay of Human Plasma. *J Endocrinol.* 1971;51(1):157-68.
2. Friesen HG, et al. Biosynthesis of Human Growth Hormone and Prolactin. *J Clin Endocrinol.* 1970;31(6):611-624.
3. Friesen HG, et al. Prolactin Synthesis in Primates. *Nature.* 1971;232:19-20.
4. Hwang P, et al. A Radioimmunoassay for Human Prolactin. *Proc Natl Acad Sci USA.* 1971 ;68(8) 1902-6.
5. Hwang P, et al. Purification of Human Prolactin. *J Biol Chem.* 1972;247(7):1955-8.
6. Laurence S, et al. A Mixed Heterologous Radioimmunoassay for Human Prolactin. *J Clin Endocrinol.* 1972;34:484-490.
7. Kleinberg DL, Frantz AG. Human Prolactin: Measurement in Plasma by In Vitro Bioassay. *J Clin Invest.* 1971;50(8):1557-68.
8. Kelly PA, et al. The Interaction of Prolactin with its Receptors in Target Tissues and its Mechanism of Action. *Recent Prog Horm Res.* 1984;40:379-439.
9. Kelly PA, et al. Studies of Insulin, Growth Hormone and Prolactin Binding: Ontogenesis, Effects of Sex and Pregnancy. *Endocrinology.* 1974;95(2):532-9.
10. Lewis UJ, et al. Human Prolactin: Isolation and Some Properties. *Biochem Biophys Res Commun.* 1971;44(5):1169-76.
11. Linzer DI, et al. Nucleotide Sequence of a Growth-related mRNA Encoding a Member of the Prolactin-growth Hormone Family. *Proc Natl Acad Sci USA.* 1984;81(14):4255-9.
12. McLeod RM. Regulation of Prolactin Secretion. In: Martin L, Ganong WF, eds: *Frontiers in Neuroendocrinology.* New York: Raven Press; 1976:169-194.
13. Nikolics K, et al. A Prolactin-inhibiting Factor Within the Precursor for Human Gonadotropin-releasing Hormone. *Nature.* 1985; 316:511-7.
14. Rasmussen C, et al. CV 205-502: A New Long-acting Drug for Inhibition of Prolactin Hypersecretion. *Clin Endocrinol (OxQ).* 1987;26:321-6.
15. Reyes FI, et al. Pituitary-ovarian Relationships Preceding the Menopause. A Cross-sectional Study of Serum Follicle-stimulating Hormone, Luteinizing Hormone, Prolactin, Estradiol, and Progesterone Levels. *Am J Obstet Gynecol.* 1977;129:557-64.
16. Riddle O, et al. The Preparation, Identification and Assay of Prolactin-a Hormone of the Anterior Pituitary. *Am J Physiol.* 1933;105:191-216.
17. Shiu RP, Iwasiow BM. Prolactin-inducible Proteins in Human Breast Cancer Cells. *J Biol Chem.* 1985;260(20):11307-13.
18. Allolio B, et al. Size Heterogeneity of Immunoreactive Prolactin in Patients with Prolactinoma. *Acta Endocrinologica (Copenh).* 1987;114(4):475-82.
19. Cavaco B, et al. Some Forms of Big Big Prolactin Behave as a Complex of Monomeric Prolactin with an Immunoglobulin G in Patients with Macroprolactinemia or Prolactinoma. *J Clin Endocrinol Metab.* 1995;80(8):2342-46.
20. Bonhoff A, et al. Identification of Macroprolactin in a Patient with Asymptomatic Hyperprolactinemia as a Stable PRL-IgG Complex. *Exp Clin Endocrinol Diabetes.* 1995;103(4):252-5.
21. Fahie-Wilson MN. Detection of Macroprolactin Hyperprolactinemia in Commercial Assays for Prolactin. *Clin Chem.* 2000;46(12):2022-3.
22. Schneider W, et al. Reactivity of Macroprolactin in Common Automated Immunoassays. *Clin Biochem.* 2001 ;34(6):469-73.

WARRANTY INFORMATION



Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.