

Human Prolactin ELISA Kit ab108679

1 Image

Overview

Product name Human Prolactin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Level 1	20	6.32	0.4	6.3%
Level 2	20	19.22	1.19	6.18%
Level 3	20	54.52	2.64	4.84%

Inter-assay

Sample	n	Mean	SD	CV%
Level 1	20	6.58	0.45	6.77%
Level 2	20	21.34	1.3	6.11%
Level 3	20	58.39	2.93	5.01%

Sample type Serum

Assay type Sandwich (quantitative)

Sensitivity 0.25 ng/ml

Range 5 ng/ml - 100 ng/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's Prolactin *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Prolactin in Human serum.

Microtiterstrip wells are precoated with inactivated specific *Corynebacterium diphtheriae* toxin (toxoid) antigens to bind corresponding antibodies of the specimen. After washing the wells to remove all unbound sample material horseradish peroxidase (HRP) labelled anti-human IgG

conjugate is added. This conjugate binds to the captured C. diphtheriae toxin-specific antibodies. The immune complex formed by the bound conjugate is visualized by adding Tetramethylbenzidine (TMB) substrate which gives a blue reaction product. The intensity of this product is proportional to the amount of C. diphtheriae toxin-specific IgG antibodies in the specimen. Sulphuric acid is added to stop the reaction. This produces a yellow endpoint colour. Absorbance at 450 nm is read using an ELISA microwell plate reader.

Platform Microplate

Properties

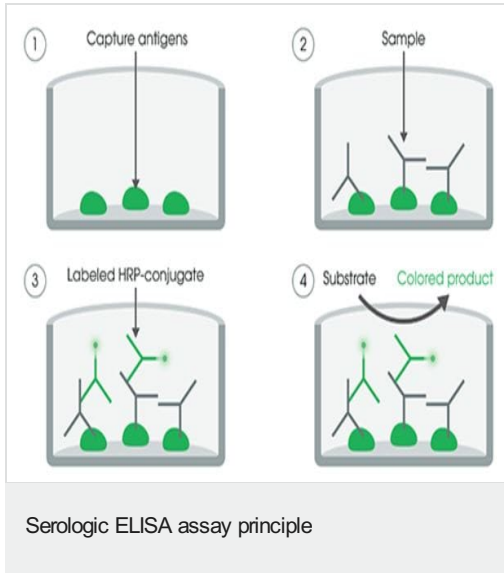
Storage instructions Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
10X Wash Solution		1 x 50ml
Anti-Prolactin-HRP conjugate		1 x 12ml
Cover foil		1 unit
Prolactin control		1 x 1ml
Prolactin Standard 0: 0 ng/ml		1 x 1ml
Prolactin Standard 1: 5 ng/ml		1 x 1ml
Prolactin Standard 2: 10 ng/ml		1 x 1ml
Prolactin Standard 3: 25 ng/ml		1 x 1ml
Prolactin Standard 4: 50 ng/ml		1 x 1ml
Prolactin Standard 5: 100 ng/ml		1 x 1ml
Stop Solution		1 x 15ml
Streptavidin Coated Microplate (12 x 8 wells)	12 Strips of 8 wells	1 unit
Strip holder		1 unit
TMB Substrate Solution		1 x 15ml

Relevance Various hormones are secreted from the anterior pituitary during development and growth. Prolactin is a growth factor secreted by the anterior pituitary that is necessary for the proliferation and differentiation of the mammary glands. Prolactin acts primarily on the mammary gland by promoting lactation.

Cellular localization Secreted

Images



Specific antigens are coated on the 96-well plate, controls or test samples are added to the well and incubated. The wells are washed to remove any unbound Human anti-antigen antibodies (Ig). A horseradish peroxidase (HRP) labelled anti-Human Ig conjugate is added to the wells. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The intensity of yellow coloration is directly proportional to the amount of Human anti-antigen Ig captured on the plate.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors