

Product datasheet

Human Progesterone ELISA Kit ab108670

★★★★★ 1 Abreviews 6 References 1 Image

Overview

Product name	Human Progesterone ELISA Kit
Detection method	Colorimetric
Sample type	Serum, Hep Plasma, EDTA Plasma, Cit plasma
Assay type	Competitive
Sensitivity	0.05 ng/ml
Range	0.2 ng/ml - 40 ng/ml
Recovery	106 %
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human

Predicted to work with: Mammals

Product overview

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

A 96-well plate has been precoated with anti-Progesterone. Samples and the Progesterone-HRP conjugate are added to the wells, where Progesterone in the sample competes with the added Progesterone-HRP for antibody binding. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is inversely proportional to the amount of Progesterone in the sample and the intensity is measured at 450 nm.

Platform Microplate

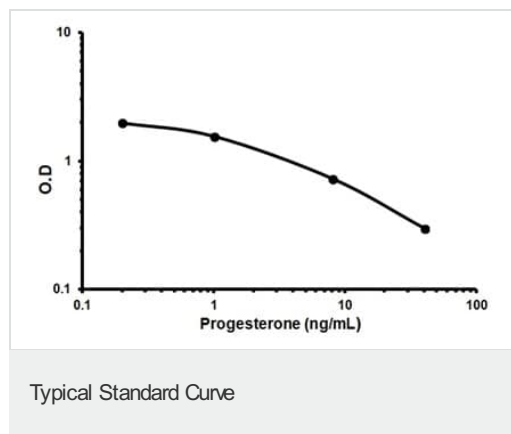
Properties

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

Components	Identifier	1 x 96 tests
10X Wash Solution		1 x 50ml
Anti-Progesterone IgG Coated Wells		1 unit

Components	Identifier	1 x 96 tests
Cover foils		1 unit
Progesterone control		1 x 1ml
Progesterone Standard 0 - 0 ng/mL		1 x 1ml
Progesterone Standard 1 - 0.2 ng/mL		1 x 1ml
Progesterone Standard 2 - 1.0 ng/mL		1 x 1ml
Progesterone Standard 3 - 8.0 ng/mL		1 x 1ml
Progesterone Standard 4 - 40.0 ng/mL		1 x 1ml
Progesterone-HRP Conjugate		1 x 22ml
Stop Solution		1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

Images



Representative Standard Curve using ab108670.

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