

Mouse Relaxin 1 ELISA Kit ab213885

1 References 1 Image

Overview

Product name Mouse Relaxin 1 ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
1	16	104pg/ml	4.16	= 4%
2	16	612pg/ml	47.12	= 7.7%
3	16	2407pg/ml	154.04	= 6.4%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	108pg/ml	7.71	= 5.2%
2	24	616pg/ml	59.13	= 9.6%
3	24	2227pg/ml	144.75	= 6.5%

Sample type Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 78 pg/ml - 5000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Mouse

Product overview The Mouse Relaxin 1 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213885) is designed for the quantitative measurement of Mouse Relaxin 1 in cell culture supernatants, serum and plasma (heparin or EDTA ).

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from rat specific for Relaxin 1 has been pre-coated onto 96-well plates. Standards (Expression system for standard: E.coli; Immunogen sequence: R23-R54 (B) & E161-C185 (A)) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Relaxin 1 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Mouse Relaxin 1 amount of sample captured in plate.

**Notes** Relaxin 1 is a member of relaxin-like peptide family. Relaxin gene maps to mouse chromosome 19 near D19Mit23. Relaxin is a peptide hormone produced by the corpora lutea of ovaries during pregnancy in many mammalian species, including man. Relaxin widens the pubic bone and facilitates labor, it also softens the cervix (cervical ripening), and relaxes the uterine musculature. However, its significance may reach much further. Relaxin affects collagen metabolism, inhibiting collagen synthesis and enhancing its breakdown by increasing matrix metalloproteinases. It also enhances angiogenesis and is a potent renal vasodilator.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

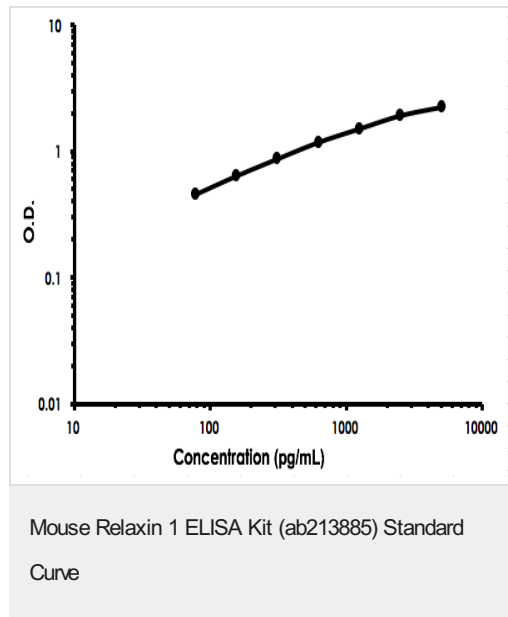
Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Mouse Relaxin 1 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Mouse Relaxin 1 antibody		1 x 100µl
Lyophilized recombinant Mouse Relaxin 1 standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

**Function** Relaxin is an ovarian hormone that acts with estrogen to produce dilatation of the birth canal in many mammals. May be involved in remodeling of connective tissues during pregnancy, promoting growth of pubic ligaments and ripening of the cervix.

**Tissue specificity** Prostate. Not expressed in placenta, decidua or ovary.

<b>Sequence similarities</b>	Belongs to the insulin family.
<b>Cellular localization</b>	Secreted.

## Images



Mouse Relaxin 1 ELISA Kit (ab213885) Standard Curve.

**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