

### Human Relaxin 1 ELISA Kit ab213820

1 Image

#### Overview

**Product name** Human Relaxin 1 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
1	16	144pg/ml	10.94	= 7.6%
2	16	819pg/ml	37.67	= 4.6%
3	16	2390pg/ml	100.38	= 4.2%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	133pg/ml	10.5	= 7.9%
2	24	796pg/ml	42.98	= 5.4%
3	24	2573pg/ml	118.35	= 4.6%

**Sample type** Cell culture supernatant, Serum, Hep 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:** Human

**Product overview** The Human Relaxin 1 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213820) is designed for the quantitative measurement of Human Relaxin 1 in cell culture supernatants, serum and plasma (heparin).

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

monoclonal antibody from mouse specific for Relaxin 1 has been pre-coated onto 96-well plates. Standards (*E.coli*, R162-Cys185(A) and K26-S53(B)) 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 is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is 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 Human Relaxin 1 amount of sample captured in plate.

**Notes** Relaxin 1 is a member of relaxin-like peptide family. Relaxin gene maps to human 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-Human Relaxin 1 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human Relaxin 1 antibody		1 x 100µl
Lyophilized recombinant Human 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.

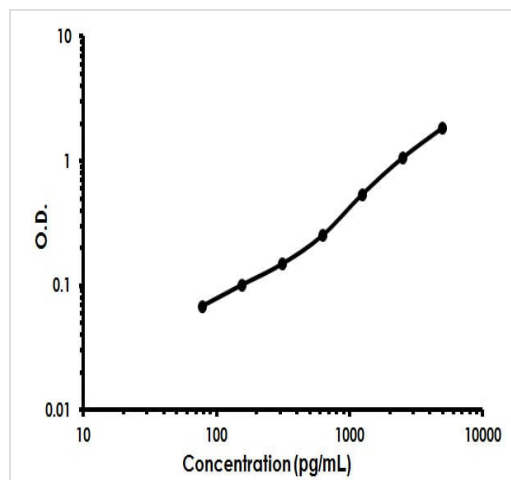
**Tissue specificity** Prostate. Not expressed in placenta, decidua or ovary.

**Sequence similarities** Belongs to the insulin family.

---

## Images

---



Human Relaxin 1 ELISA Kit (ab213820) Standard Curve.

Human Relaxin 1 ELISA Kit (ab213820) 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