

### Human Activin A ELISA Kit ab119568

#### 画像数 1

#### 製品の概要

製品名	Human Activin A ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Other biological fluids, Tissue Extracts
アッセイタイプ	Sandwich (quantitative)
検出感度	< 12 pg/ml
検出範囲	62.5 pg/ml - 4000 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's Human Activin A <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human Activin A in cell culture supernatants, serum, plasma (heparin, EDTA) and saliva.</p> <p>A Activin A specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat, specific for Activin A is then added 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. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human Activin A amount of sample captured in plate.</p>
試験プラットフォーム	Microplate

#### 製品の特性

**保存方法** Store at -20°C. Please refer to protocols.

内容	ラベル	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Activin A Antibody Microplate (12 x 8 wells)		1 unit	1 unit

内容	ラベル	1 x 96 tests	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Human Activin A antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Activin A standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

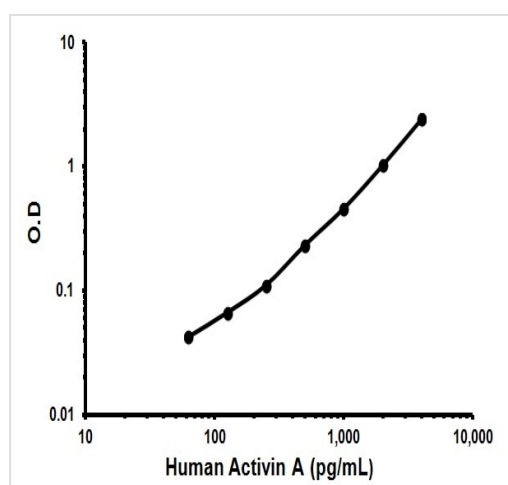
## 関連性

Activin A is a homodimer of 14kDa beta-A. Activin A, a cytokine member of the transforming growth factor-beta superfamily, is expressed locally by the mesenchymal component of the hemopoietic microenvironment. Its expression is regulated on the mRNA level by different cytokines, and the biological activity of the protein is tightly controlled by several inhibitory molecules.<sup>1</sup> Inhibins and activins are members of the transforming growth factor beta superfamily and are known to modulate the growth and differentiation of several cell types.<sup>2</sup> Inhibins and activins inhibit and activate, respectively, the secretion of follitropin by the pituitary gland. Inhibins/activins are involved in regulating a number of diverse functions such as hypothalamic and pituitary hormone secretion, gonadal hormone secretion, germ cell development and maturation, erythroid differentiation, insulin secretion, nerve cell survival, embryonic axial development or bone growth, depending on their subunit composition. Inhibins appear to oppose the functions of activins. The standard product used in this kit is recombinant Activin A, which is composed of two single chains of 116 amino acids with the molecular mass of 26kDa.

## 細胞内局在

Secreted

## 画像



Representative Standard Curve using ab119568.

Typical 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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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