

Rat Activin A ELISA Kit ab193733

[1 References](#) [2 Images](#)

Overview

Product name Rat Activin A ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 2 pg/ml

Range 2 pg/ml - 400 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	= 143.3	129% - 150%
Plasma	= 138.7	127% - 150%
Cell culture media	= 121.7	105% - 149%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse, Rat, Human

Product overview

Abcam's Rat Activin A ELISA Kit (ab193733) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of rat Activin A in sera, plasma and cell culture supernatants.

This assay employs an antibody specific for rat Activin A coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures Activin A present in the samples. The wells are washed and biotinylated anti-rat Activin A antibody is added. After

washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of Activin A bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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

Properties

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

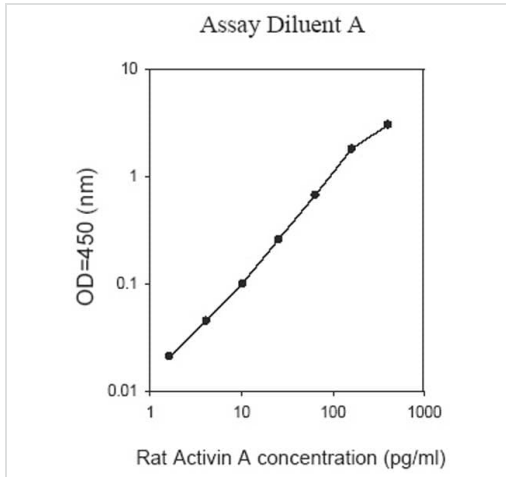
Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated Rat Activin A Detection antibody	2 vials
Pre-coated rat Activin A Microplate (12 strips x 8 wells)	1 unit
Rat Activin A Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

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.¹ 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.² 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.

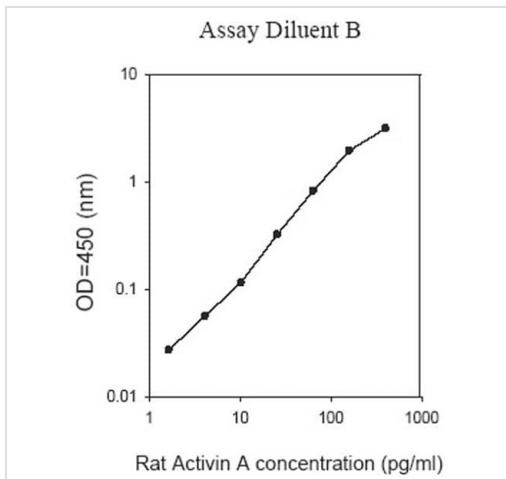
Cellular localization Secreted

Images



A typical standard curve using ab193733 with Assay Diluent A

Typical Standard Curve - Diluent A



A typical standard curve using ab193733 with Assay Diluent B

Typical Standard Curve - Diluent B

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