

Human ADAM9 ELISA Kit ab193690

2 Images

Overview

Product name	Human ADAM9 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	45 pg/ml				
Range	45 pg/ml - 12000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		96.06		88% - 104%
	Plasma		91.5		81% - 101%
	Cell culture media		122.2		114% - 130%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's ADAM9 Human ELISA Kit (ab193690) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Human ADAM9 in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for Human ADAM9 coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures ADAM9 present in

the samples. The wells are washed and biotinylated anti-Human ADAM9 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 ADAM9 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
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Detection Antibody ADAM9 (biotinylated anti-Human ADAM9)	2 vials
Human ADAM9 Standards (lyophilized)	2 vials
Pre-coated Human ADAM9 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Probable zinc protease. May mediate cell-cell or cell-matrix interactions. Isoform 2 displays alpha-secretase activity for APP.

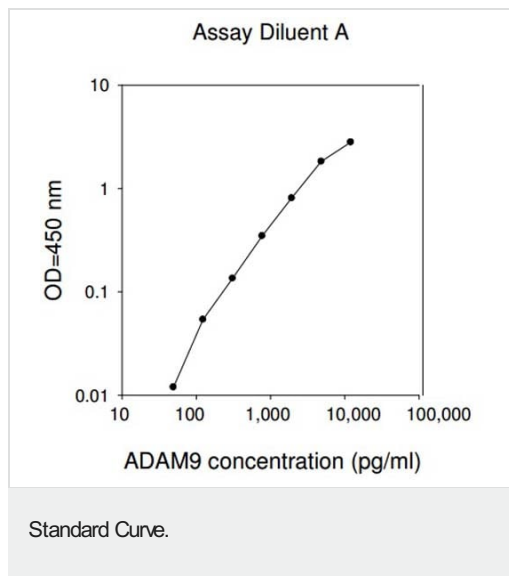
**Tissue specificity** Widely expressed. Expressed in chondrocytes. Isoform 2 is highly expressed in liver and heart.

**Involvement in disease** Cone-rod dystrophy 9

**Sequence similarities** Contains 1 disintegrin domain.  
Contains 1 EGF-like domain.  
Contains 1 peptidase M12B domain.

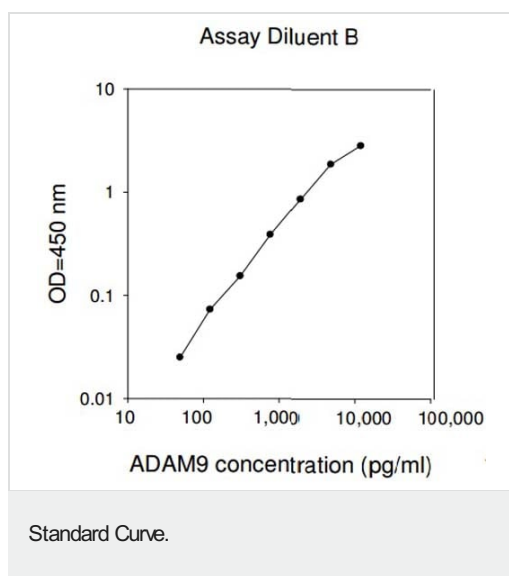
**Cellular localization** Secreted and Cell membrane.

## Images



Standard Curve in Assay Diluent A using ab193690 ADAM9 Human ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Standard Curve in Assay Diluent B using ab193690 ADAM9 Human ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

**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