

### Mouse Renin 1 ELISA Kit ab193728

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

#### Overview

**Product name** Mouse Renin 1 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** 6 pg/ml

**Range** 6 pg/ml - 50000 pg/ml

**Recovery** Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	= 107.3	87% - 132%
Serum	= 130.5	112% - 145%
Plasma	= 134.4	128% - 142%

**Assay duration** One step assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** Abcam's Renin 1 mouse ELISA kit (ab193728) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of mouse Renin 1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse Renin 1 coated on a 96 well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures Renin 1 present in

the sample. The wells are washed and biotinylated anti-mouse Renin 1 antibody is added. After washing away unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted into the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color develops in proportion to the amount of Renin 1 bound in each well. Addition of the Stop Solution changes 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
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated Mouse Renin 1 detection antibody	2 vials
mouse Renin 1 Standards (lyophilized)	2 vials
Pre-coated mouse Renin1 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

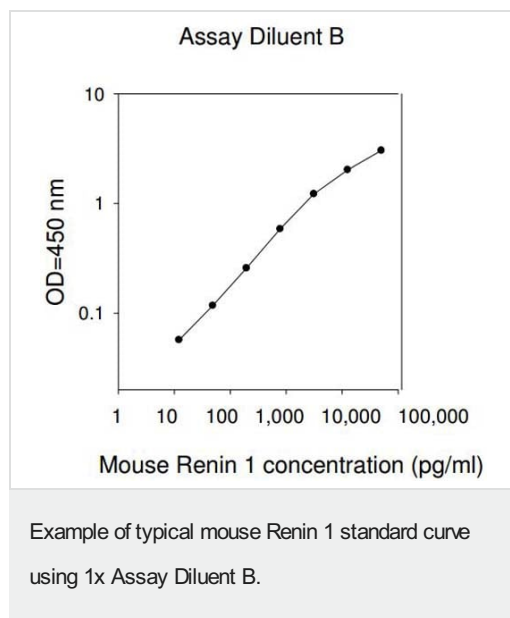
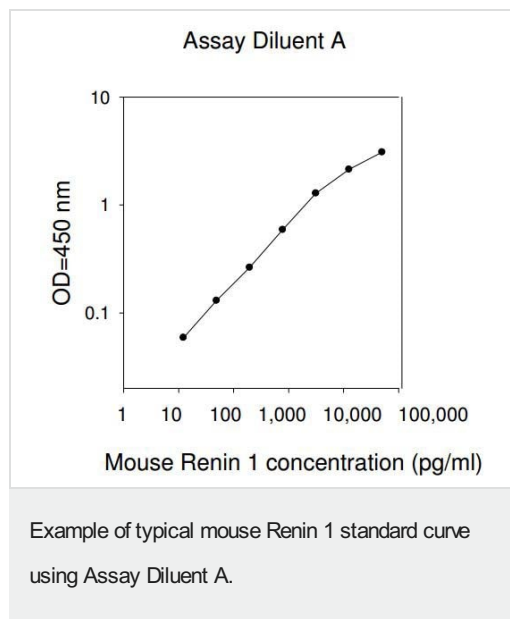
**Function** Functions as a renin and prorenin cellular receptor. May mediate renin-dependent cellular responses by activating ERK1 and ERK2. By increasing the catalytic efficiency of renin in AGT/angiotensinogen conversion to angiotensin I, it may also play a role in the renin-angiotensin system (RAS).

**Tissue specificity** Expressed in brain, heart, placenta, liver, kidney and pancreas. Barely detectable in lung and skeletal muscles. In the kidney cortex it is restricted to the mesangium of glomeruli. In the coronary and kidney artery it is expressed in the subendothelium, associated to smooth muscles where it colocalizes with REN. Expressed in vascular structures and by syncytiotrophoblast cells in the mature fetal placenta.

**Involvement in disease** Defects in ATP6AP2 are a cause of mental retardation X-linked with epilepsy (MRXE) [MIM:300423]. MRXE is a syndromic mental retardation. Patients manifest mild to moderate mental retardation associated with epilepsy, delays in motor milestones and speech acquisition in infancy.

**Post-translational modifications** Phosphorylated.

**Cellular localization** Membrane.



**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