

# Mouse Albumin ELISA Kit ab108792

★★★★★ [1 Abreviews](#) [78 References](#) [1 Image](#)

### Overview

**Product name** Mouse Albumin ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				5.2%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				9.8%

**Sample type** Cell culture supernatant, Urine, Serum, Plasma, Cell Lysate

**Assay type** Sandwich (quantitative)

**Sensitivity** > 0.44 ng/ml

**Range** 3.125 ng/ml - 50 ng/ml

**Recovery** 97 %

**Assay duration** Multiple steps standard assay

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

**Product overview** Mouse Albumin ELISA kit (ab108792) is a sandwich ELISA kit designed for the quantitative measurement of mouse Albumin levels in urine and cell culture supernatants.

An Albumin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Albumin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase 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 amount of Albumin captured in plate.

**The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid**

**repeated freeze-thaw cycles.**

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

**Platform** Microplate

**Properties**

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

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Mouse Albumin Antibody	1 x 120µl
Albumin Microplate (12 x 8 wells)	1 unit
Albumin Standard (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

**Function** Serum albumin, the main protein of plasma, has a good binding capacity for water, Ca(2+), Na(+), K(+), fatty acids, hormones, bilirubin and drugs. Its main function is the regulation of the colloidal osmotic pressure of blood. Major zinc transporter in plasma, typically binds about 80% of all plasma zinc.

**Tissue specificity** Plasma.

**Involvement in disease** Defects in ALB are a cause of familial dysalbuminemic hyperthyroxinemia (FDH) [MIM:103600]. FDH is a form of euthyroid hyperthyroxinemia that is due to increased affinity of ALB for T(4). It is the most common cause of inherited euthyroid hyperthyroxinemia in Caucasian population.

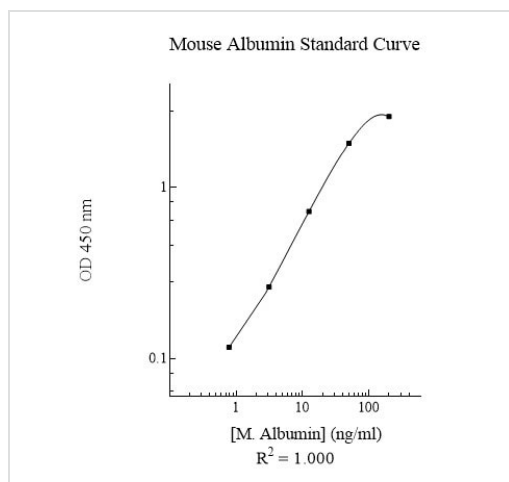
**Sequence similarities** Belongs to the ALB/AFP/VDB family.  
Contains 3 albumin domains.

**Post-translational modifications** Kenitra variant is partially O-glycosylated at Thr-620. It has two new disulfide bonds Cys-600 to Cys-602 and Cys-601 to Cys-606.  
Glycated in diabetic patients.  
Phosphorylation sites are present in the extracellular medium.  
Acetylated on Lys-223 by acetylsalicylic acid.

**Cellular localization** Secreted.

**Images**

Representative Standard Curve Using ab108792.



Typical Standard Curve for ab108792 Mouse Albumin ELISA kit

**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