

Product datasheet

Albumin Rabbit ELISA Kit ab108793

1 Image

Overview

Product name Albumin Rabbit ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall 4.8%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall 9.8%

Sample type Cell culture supernatant, Urine, Serum, Plasma

Assay type Competitive

Sensitivity > 0.19 µg/ml

Range 0.4 µg/ml - 40 µg/ml

Recovery 98 %

Assay time 3h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rabbit

Product overview

Abcam's Albumin rabbit *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of rabbit Albumin in plasma, serum, urine, and cell culture supernatant.

An Albumin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Albumin 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 inversely 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
2X Biotinylated Albumin	1 unit
20X Wash Buffer Concentrate	1 x 30ml
Albumin Microplate (12 x 8 well strips)	1 unit
Albumin Standard	1 unit
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml

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.

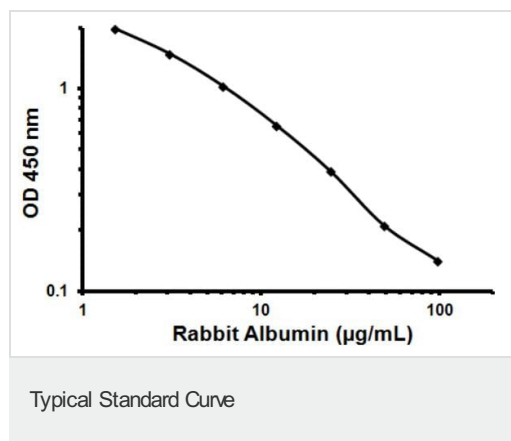
Glycated in diabetic patients.

Phosphorylation sites are present in the extracellular medium.

Acetylated on Lys-223 by acetylsalicylic acid.

Cellular localization Secreted.

Images



Representative standard curve using ab108793

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