

Product datasheet

Mouse Albumin ELISA Kit ab108791

1 References 3 图像

概述

产品名称 小鼠Albumin ELISA试剂盒

检测方法 Colorimetric

精确度

批次内				
样品	n	Mean	SD	CV%
Overall				4.4%

批次间				
样品	n	Mean	SD	CV%
Overall				7.1%

样品类型 Serum, Plasma, Tissue Lysate

检测类型 Competitive

灵敏度 = 0.09 µg/ml

范围 0.098 µg/ml - 25 µg/ml

回收率 98 %

检测时间 2h 00m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

预测可用于: Rat

产品概述 Mouse Albumin *in vitro* competitive ELISA (enzyme-linked immunosorbent assay) kit is designed for the quantitative measurement of mouse Albumin levels in plasma and serum.

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 PBS or TBS buffer. Streptavidin-Peroxidase Complex 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.

经测试应用

适用于: Competitive ELISA

平台

Microplate

性能

存放说明

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

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

功能

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.

组织特异性

Plasma.

疾病相关

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.

序列相似性

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

翻译后修饰

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.

细胞定位

Secreted.

应用

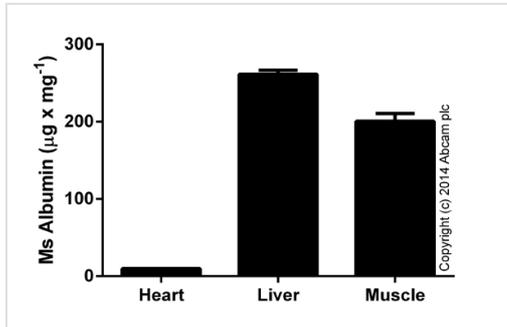
Our [Abpromise guarantee](#) covers the use of **ab108791** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Competitive ELISA

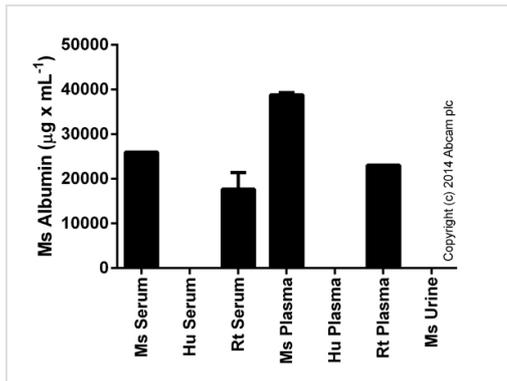
Use at an assay dependent dilution.

图片



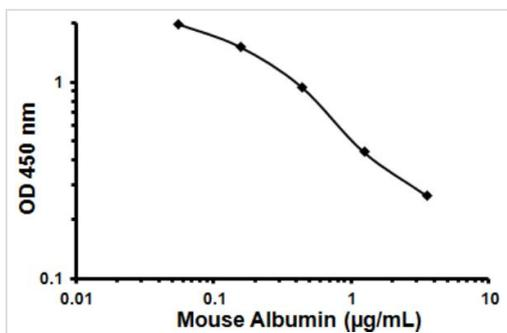
Albumin measured in mouse tissue lysates (expressed as per mg of extracted protein), with background signal subtracted (duplicates; +/- SD).

Competitive ELISA - Albumin Mouse ELISA Kit
(ab108791)



Albumin measured in biological fluids, background signal subtracted (duplicates +/- SD).

Competitive ELISA - Albumin Mouse ELISA Kit
(ab108791)



Representative Standard Curve Using
ab108791.

Typical Standard Curve

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors