

Mouse ACE ELISA Kit (CD143) ab155452

2 References 2 图像

概述

产品名称 小鼠ACE ELISA试剂盒(CD143)

检测方法 Colorimetric

精确度					批次内
	样品	n	Mean	SD	CV%
	overall				< 10%

					批次间
	样品	n	Mean	SD	CV%
	overall				< 12%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 0.12 ng/ml

范围 0.123 ng/ml - 30 ng/ml

回收率			特定样本回收率
	样品类型	平均%	范围
	Serum	104.6	96% - 112%
	Plasma	101.4	94% - 109%
	Cell culture media	99.78	90% - 109%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's ACE (CD143) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse ACE (CD143) in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse ACE coated on a 96-well plate. Standards and samples are pipetted into the wells and ACE present in a sample is bound to the wells by the

immobilized antibody. The wells are washed and biotinylated anti-mouse ACE antibody is added. After washing away unbound biotinylated antibody,

HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of ACE bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
ACE Microplate (12 x 8 wells)	1 unit
Assay Diluent A	1 x 30ml
Biotinylated anti-mouse ACE	2 vials
Recombinant Mouse ACE Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 8ml

功能 Converts angiotensin I to angiotensin II by release of the terminal His-Leu, this results in an increase of the vasoconstrictor activity of angiotensin. Also able to inactivate bradykinin, a potent vasodilator. Has also a glycosidase activity which releases GPI-anchored proteins from the membrane by cleaving the mannose linkage in the GPI moiety.

组织特异性 Ubiquitously expressed, with highest levels in lung, kidney, heart, gastrointestinal system and prostate. Isoform Testis-specific is expressed in spermatocytes and adult testis.

疾病相关 Ischemic stroke (ISCHSTR) [MIM:601367]: A stroke is an acute neurologic event leading to death of neural tissue of the brain and resulting in loss of motor, sensory and/or cognitive function. Ischemic strokes, resulting from vascular occlusion, is considered to be a highly complex disease consisting of a group of heterogeneous disorders with multiple genetic and environmental risk factors. Note=Disease susceptibility is associated with variations affecting the gene represented in this entry.

Renal tubular dysgenesis (RTD) [MIM:267430]: Autosomal recessive severe disorder of renal tubular development characterized by persistent fetal anuria and perinatal death, probably due to pulmonary hypoplasia from early-onset oligohydramnios (the Potter phenotype). Note=The disease is caused by mutations affecting the gene represented in this entry.

Microvascular complications of diabetes 3 (MVCD3) [MIM:612624]: Pathological conditions that develop in numerous tissues and organs as a consequence of diabetes mellitus. They include

diabetic retinopathy, diabetic nephropathy leading to end-stage renal disease, and diabetic neuropathy. Diabetic retinopathy remains the major cause of new-onset blindness among diabetic adults. It is characterized by vascular permeability and increased tissue ischemia and angiogenesis. Note=Disease susceptibility is associated with variations affecting the gene represented in this entry.

Intracerebral hemorrhage (ICH) [MIM:614519]: A pathological condition characterized by bleeding into one or both cerebral hemispheres including the basal ganglia and the cerebral cortex. It is often associated with hypertension and craniocerebral trauma. Intracerebral bleeding is a common cause of stroke. Note=Disease susceptibility is associated with variations affecting the gene represented in this entry.

序列相似性

Belongs to the peptidase M2 family.

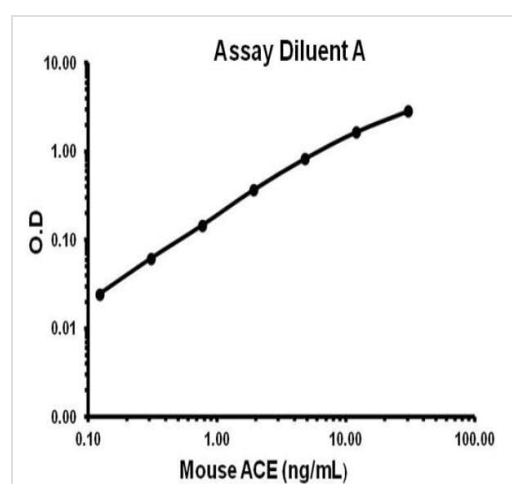
翻译后修饰

Phosphorylated by CK2 on Ser-1299; which allows membrane retention.

细胞定位

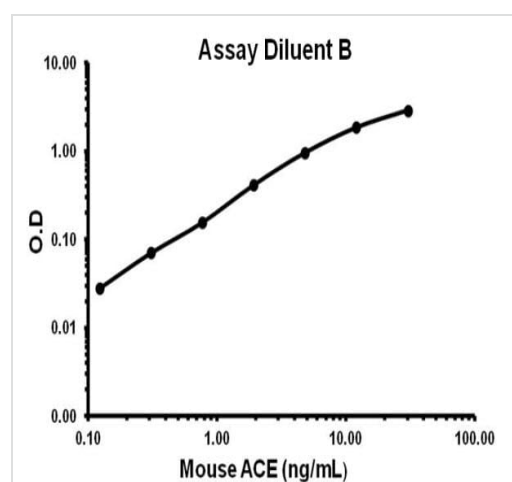
Secreted and Cell membrane.

图片



Representative standard curve using ab155452 - Assay Diluent A.

Typical Standard Curve



Representative standard curve using ab155452 - Assay Diluent B.

Typical Standard Curve

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.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