

Rat ADAM12 ELISA Kit ab213919

1 图像

概述

产品名称	大鼠ADAM12 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	1	16	217pg/ml	9.54	= 4.4%
	2	16	1804pg/ml	137.1	= 7.6%
	3	16	6261pg/ml	306.78	= 4.9%
	批次间				
	样品	n	Mean	SD	CV%
	1	24	237pg/ml	10.9	= 4.6%
	2	24	1849pg/ml	144.22	= 7.8%
	3	24	6482pg/ml	388.92	= 6%
样品类型	Cell culture supernatant, Serum, Hep Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 10 pg/ml				
范围	156 pg/ml - 10000 pg/ml				
检测时间	3h 30m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Rat				
产品概述	The Rat ADAM12 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213919) is designed for the quantitative measurement of Rat ADAM12 in cell culture supernatants, serum and plasma (heparin).				

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for ADAM12 has been precoated onto 96-well plates. Standards and test samples are added to the wells; a biotinylated detection polyclonal antibody from goat specific for ADAM12 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Rat ADAM12 amount of sample captured in plate.

说明

ADAM12 (Disintegrin and metalloproteinase domain-containing protein 12), also known as MLTN, is an enzyme that in humans is encoded by the ADAM12 gene. This gene encodes a member of the ADAM (a disintegrin and metalloprotease) protein family. Members of this family are membrane anchored proteins structurally related to snake venom disintegrins, and have been implicated in a variety of biological processes involving cell-cell and cell-matrix interactions, including fertilization, muscle development, and neurogenesis. This gene has two alternatively spliced transcripts: a shorter secreted form and a longer membrane-bound form. The shorter form is found to stimulate myogenesis. By RT-PCR and immunoblot analyses that expression of mouse Adam12 increases during muscle regeneration, while the levels of other ADAMs remain constant. Immunofluorescence analysis revealed staining of small, newly formed muscle fibers in regenerating but not normal adult muscle cells. By using of fluorescence in situ hybridization, the ADAM12 gene is mapped to human chromosome 10q26.3.

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

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

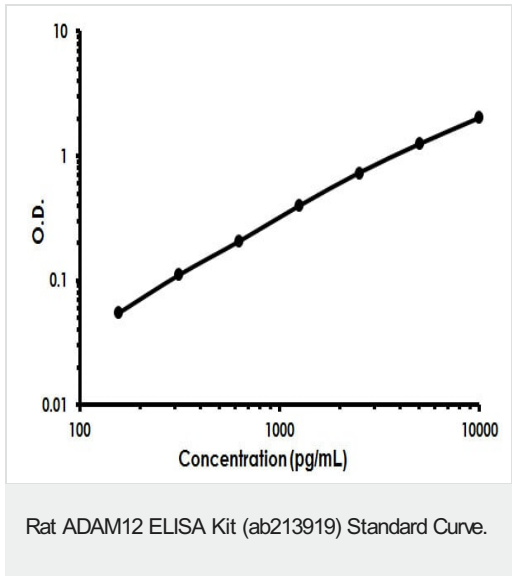
组件	标识符	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-rat ADAM12 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- rat ADAM12 antibody		1 x 100µl
Lyophilized recombinant rat ADAM12 standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

功能

Involved in skeletal muscle regeneration, specifically at the onset of cell fusion. Also involved in

	macrophage-derived giant cells (MGC) and osteoclast formation from mononuclear precursors.
组织特异性	Isoform 1 is expressed in placenta and skeletal, cardiac, and smooth muscle. Isoform 2 seems to be expressed only in placenta or in embryo and fetus. Both forms were expressed in some tumor cells lines. Not detected in brain, lung, liver, kidney or pancreas.
序列相似性	Contains 1 disintegrin domain. Contains 1 EGF-like domain. Contains 1 peptidase M12B domain.
结构域	The cysteine-rich domain supports cell adhesion through syndecans and triggers signaling events that lead to beta-1 integrin-dependent cell spreading. In carcinomas cells the binding of this domain to syndecans does not allow the integrin-mediated cell spreading. The conserved cysteine present in the cysteine-switch motif binds the catalytic zinc ion, thus inhibiting the enzyme. The dissociation of the cysteine from the zinc ion upon the activation-peptide release activates the enzyme.
翻译后修饰	The precursor is cleaved by a furin endopeptidase.
细胞定位	Secreted and Cell membrane.

图片



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