

Mouse Angiopoietin 2 ELISA Kit ab209883

重组 SimpleStep ELISA

3 References 4 图像

概述

产品名称 小鼠Angiopoietin 2 ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Mouse plasma	3			8.3%

批次间				
样品	n	Mean	SD	CV%
Mouse plasma	5			10.3%

样品类型 Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

检测类型 Sandwich (quantitative)

灵敏度 17.3 pg/ml

范围 68.75 pg/ml - 4400 pg/ml

特定样本回收率		
样品类型	平均%	范围
Serum	94.5	67.4% - 120.9%
Cell culture media	99.7	94.3% - 103.3%
Hep Plasma	112.4	81.4% - 143.2%
EDTA Plasma	104.9	91.8% - 122.8%
Cit plasma	94.7	87.1% - 103.3%

检测时间 1h 30m

实验步骤 One step assay

种属反应性

与反应: Mouse
不与反应: Cow

产品概述

Mouse Angiopoietin 2 ELISA Kit (ab209883) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Angiopoietin 2 protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse Angiopoietin 2 with 17.3 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

Angiopoietin 2 is a secreted glycosylated protein functioning in angiogenesis, inflammation and vascular development. Angiopoietin 2 binds to receptor tyrosine kinase TEK/TIE2, competing for the ANGPT1 binding site, and modulating ANGPT1 signalling. Angiopoietin 2 can induce tyrosine phosphorylation of TEK/TIE2 in the absence of ANGPT1. In the absence of angiogenic inducers, such as VEGF, Angiopoietin 2-mediated loosening of cell-matrix contacts may induce endothelial cell apoptosis with consequent vascular regression. In concert with VEGF, Angiopoietin 2 may facilitate endothelial cell migration and proliferation, thus serving as a permissive angiogenic signal.

平台

Pre-coated microplate (12 x 8 well strips)

性能

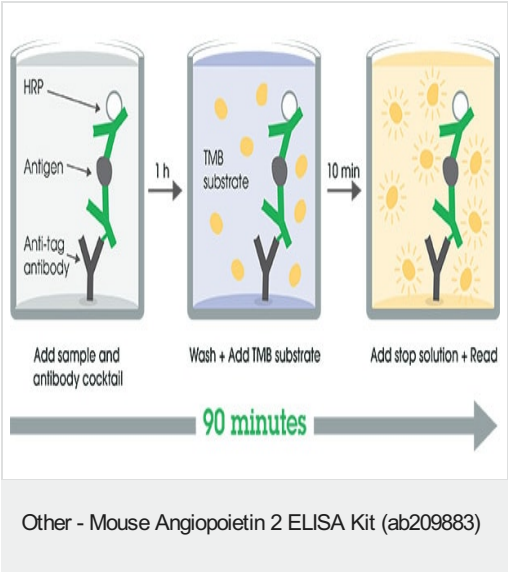
存放说明Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Mouse Angiopoietin 2 Capture Antibody	1 x 600µl
10X Mouse Angiopoietin 2 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse Angiopoietin 2 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit

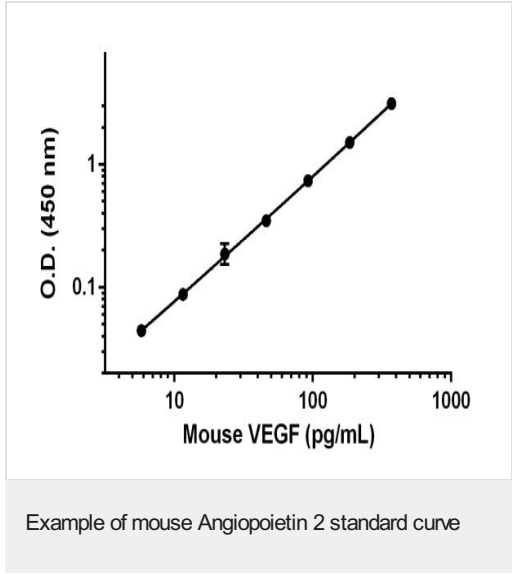
组件	1 x 96 tests
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

功能	Can induce tyrosine phosphorylation of TIE2. Binds to TIE2 receptor and counteracts blood vessel maturation/stability mediated by angiopoietin-1. Its function may be context-dependent. In the absence of angiogenic inducers, such as VEGF, ANG2-mediated loosening of cell-matrix contacts may induce endothelial cell apoptosis with consequent vascular regression. In concert with VEGF, it may facilitate endothelial cell migration and proliferation, thus serving as a permissive angiogenic signal.
序列相似性	Contains 1 fibrinogen C-terminal domain.
结构域	The Fibrinogen C-terminal domain mediates interaction with the TEK/TIE2 receptor.
细胞定位	Secreted.

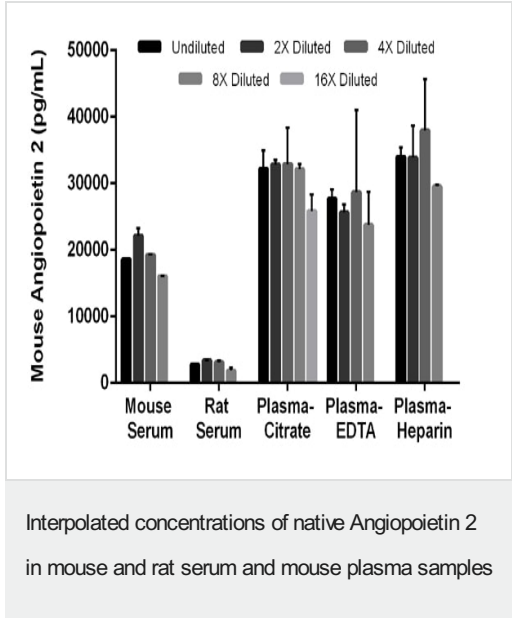
图片



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Background-subtracted data values (mean \pm SD) are graphed.



The concentrations of Angiopoietin 2 were measured in duplicates, interpolated from the Angiopoietin 2 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 25%, plasma (citrate) 6%, plasma (EDTA) 7.5%, plasma (heparin) 7.5%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean Angiopoietin 2 concentration was determined to be 19,017 pg/mL in mouse serum, 2,802 pg/mL in rat serum, 31,192 pg/mL in mouse plasma (citrate), 26,451 pg/mL in mouse plasma (EDTA) and 33,857 pg/mL in mouse plasma (heparin).

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Mouse Angiopoietin 2 ELISA Kit
(ab209883)

To learn more about the advantages of recombinant antibodies see [here](#).

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