

### Mouse Angiopoietin 2 ELISA Kit ab171335

[3 References](#) [3 Images](#)

#### Overview

**Product name** Mouse Angiopoietin 2 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
1	16	210pg/ml	10.5	= 5%
2	16	1256pg/ml	55.06	= 4.4%
3	16	4765pg/ml	274.67	= 5.8%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	230pg/ml	11.5	= 5%
2	24	1185pg/ml	54.51	= 4.6%
3	24	4889pg/ml	303.11	= 6.2%

**Sample type** Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** < 10 pg/ml

**Range** 156 pg/ml - 10000 pg/ml

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** Abcam's mouse Angiopoietin-2 *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of mouse Angiopoietin-2 in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

An Angiopoietin-2 specific rat monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection

polyclonal antibody from goat, specific for Angiopoietin-2 is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the mouse Angiopoietin-2 amount of sample captured in plate.

Get better reproducibility in only 90 minutes with Mouse Angiopoietin 2 ELISA Kit ([ab209883](#)) from our SimpleStep ELISA® range.

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

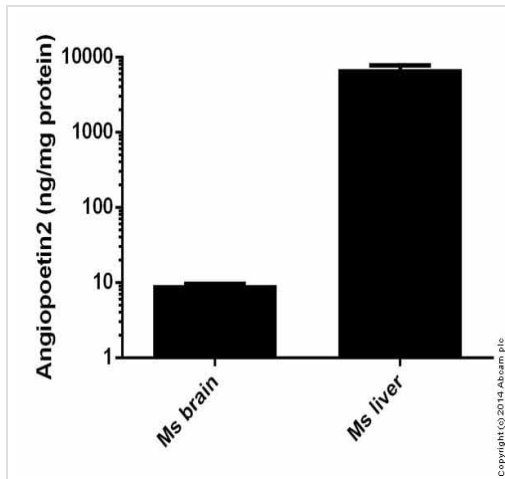
Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-mouse Angiopoietin-2 Antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-mouse Angiopoietin 2 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant mouse Angiopoietin 2 standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

**Function** 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.

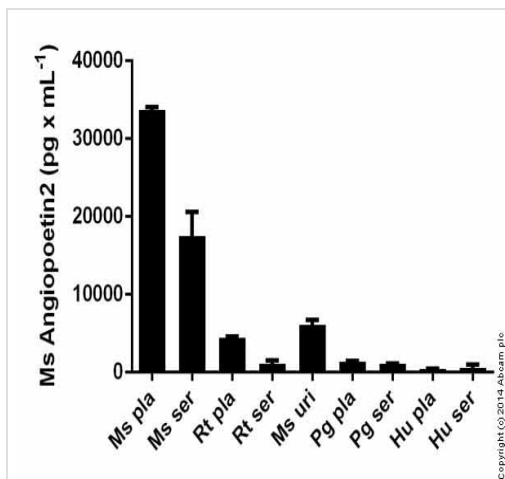
**Sequence similarities** Contains 1 fibrinogen C-terminal domain.

**Domain** The Fibrinogen C-terminal domain mediates interaction with the TEK/TIE2 receptor.

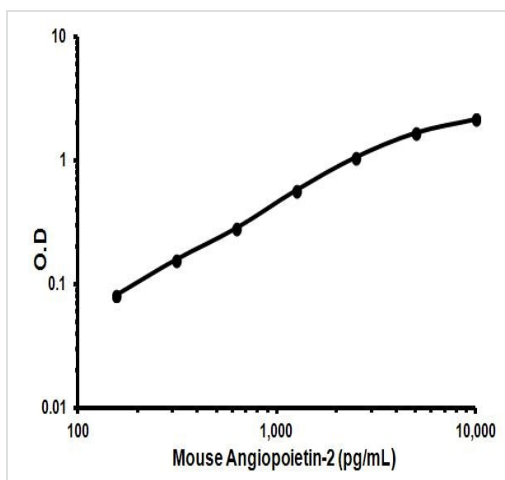
**Cellular localization** Secreted.



ab171335 Angiopoietin 2 Mouse ELISA kit



ab171335 Angiopoietin 2 Mouse ELISA kit



Typical standard curve

Representative standard curve using ab171335

## **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