

Mouse Arginase 1 ELISA Kit ab269541

Recombinant SimpleStep ELISA

3 References 5 Images

Overview

Product name	Mouse Arginase 1 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Serum	8			4.1%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Serum	3			4.7%
Sample type	Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Cit plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	2.02 pg/ml				
Range	7.81 pg/ml - 500 pg/ml				
Recovery	Sample specific recovery				
	Sample type	Average %		Range	
	Serum	99		83% - 111%	
	Cell Lysate	111		106% - 118%	
	Hep Plasma	95		82% - 102%	
	EDTA Plasma	101		97% - 104%	
	Cit plasma	106		101% - 113%	
Assay time	1h 30m				
Assay duration	One step assay				

Species reactivity**Reacts with:** Mouse**Product overview**

Mouse Arginase 1 ELISA Kit (ab269541) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Arginase 1 protein in edta plasma, hep plasma, serum, cit plasma, and cell lysate. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse Arginase 1 with 2.02 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.

Notes

Arginase is a key element of the urea cycle involving the conversion of L-arginine to urea and L-ornithine, which is further metabolized into metabolites that drive collagen synthesis and bioenergetic pathways critical for cell proliferation. Arginine metabolism is a critical regulator of innate and adaptive immune responses, involved in an antimicrobial effector pathway carried out by polymorphonuclear granulocytes (PMN). During PMN apoptosis, arginine is depleted in the microenvironment leading to suppressed T cell and natural killer cell proliferation and cytokine secretion. Arginase also plays a key role in human sexual response in smooth muscle tissue.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties**Storage instructions**

Store at +4°C. Please refer to protocols.

Components	1 x 96 tests	1 x 96 tests
10X Mouse Arginase1 Capture Antibody	1 x 600µl	1 x 600µl
10X Mouse Arginase1 Detector Antibody	1 x 600µl	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 20ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml	1 x 10ml
Antibody Diluent CPR2	1 x 6ml	1 x 6ml

Components	1 x 96 tests	1 x 96 tests
Mouse Arginase1 Lyophilized Recombinant Protein	2 vials	2 vials
Plate Seals	1 unit	1 unit
Sample Diluent NS (ab193972)	1 x 50ml	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	1 unit
Stop Solution	1 x 12ml	1 x 12ml
TMB Development Solution	1 x 12ml	1 x 12ml

Pathway

Nitrogen metabolism; urea cycle; L-ornithine and urea from L-arginine: step 1/1.

Involvement in disease

Defects in ARG1 are the cause of argininemia (ARGIN) [MIM:207800]; also known as hyperargininemia. Argininemia is a rare autosomal recessive disorder of the urea cycle. Arginine is elevated in the blood and cerebrospinal fluid, and periodic hyperammonemia occurs. Clinical manifestations include developmental delay, seizures, mental retardation, hypotonia, ataxia, progressive spastic quadriplegia.

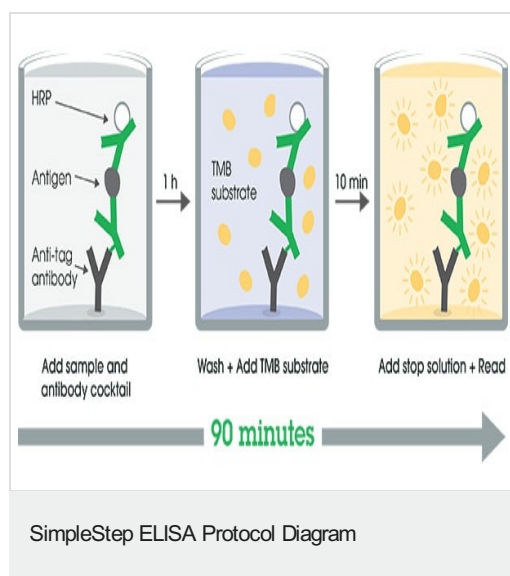
Sequence similarities

Belongs to the arginase family.

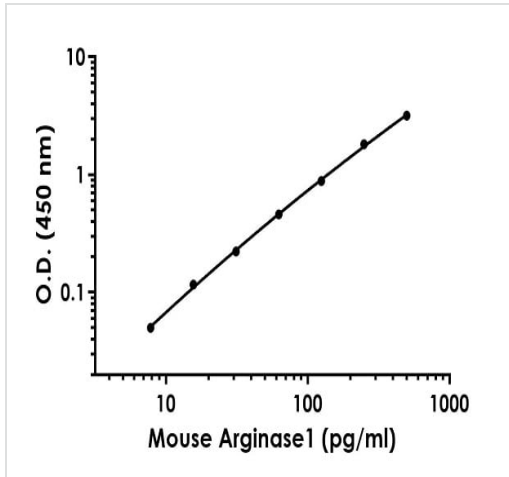
Cellular localization

Cytoplasm.

Images

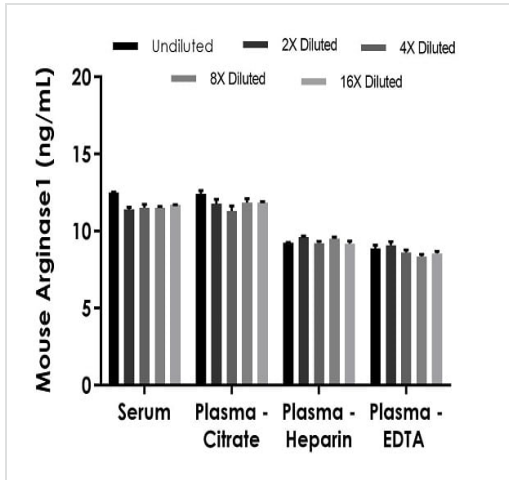


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



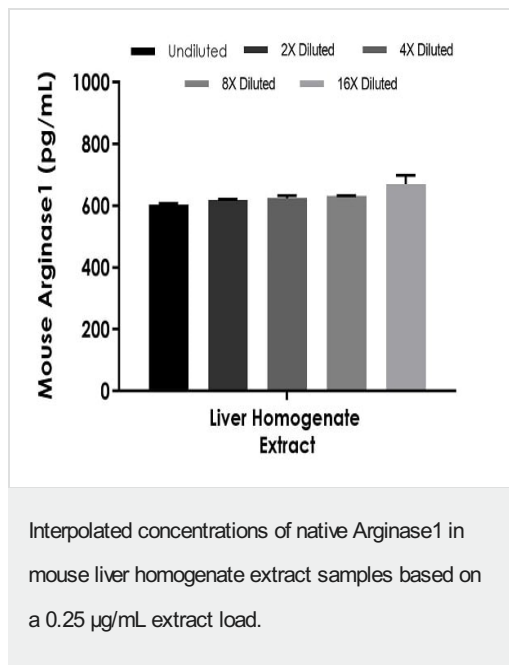
The Arginase1 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean \pm SD) are graphed.

Example of mouse Arginase1 standard curve in Sample Diluent NS.



The concentrations of Arginase1 were measured in duplicates, interpolated from the Arginase1 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1: 40, plasma (citrate) 1: 40, plasma (heparin) 1: 20, and plasma (EDTA) 1: 20. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean Arginase1 concentration was determined to be 11.61 ng/mL in serum, 11.86 ng/mL in plasma (citrate), 9.37 ng/mL in plasma (heparin), and 8.70 ng/mL in plasma (EDTA).

Interpolated concentrations of native Arginase1 in mouse serum and plasma samples.



The concentrations of Arginase1 were measured in duplicate and interpolated from the Arginase1 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Arginase1 concentration was determined to be 636.63 pg/mL in 0.25 µg/mL liver homogenate extract.

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 Arginase 1 ELISA Kit (ab269541)

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.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