

Mouse Haptoglobin ELISA Kit ab157714

5 References 1 Image

Overview

Product name	Mouse Haptoglobin ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
Sample type	Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.484 ng/ml				
Range	1.95 ng/ml - 125 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		> 85		% - %
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Mouse Haptoglobin ELISA kit is designed for the quantitative measurement of Haptoglobin in mouse serum, plasma and other biological samples.				

In this assay the Haptoglobin present in samples reacts with the anti-Haptoglobin antibodies which have been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti-Haptoglobin antibodies conjugated with horseradish

peroxidase (HRP) are added. These enzyme-labeled antibodies form complexes with the previously bound Haptoglobin. Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of Haptoglobin in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Haptoglobin in the test sample. The quantity of Haptoglobin in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Platform Microplate

Properties

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

Components	1 x 96 tests
100X HRP-conjugated anti-mouse Haptoglobin antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Mouse Haptoglobin Calibrator (lyophilized)	1 vial
Mouse Haptoglobin ELISA Microplate	1 unit
Stop Solution	1 x 12ml

Function As a result of hemolysis, hemoglobin is found to accumulate in the kidney and is secreted in the urine. Haptoglobin captures, and combines with free plasma hemoglobin to allow hepatic recycling of heme iron and to prevent kidney damage. Haptoglobin also acts as an Antimicrobial; Antioxidant, has antibacterial activity and plays a role in modulating many aspects of the acute phase response. Hemoglobin/haptoglobin complexes are rapidly cleared by the macrophage CD163 scavenger receptor expressed on the surface of liver Kupfer cells through an endocytic lysosomal degradation pathway. Uncleaved haptoglobin, also known as zonulin, plays a role in intestinal permeability, allowing intercellular tight junction disassembly, and controlling the equilibrium between tolerance and immunity to non-self antigens.

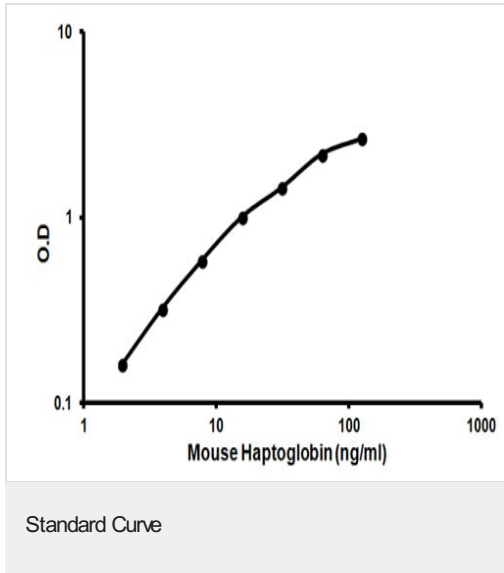
Tissue specificity Expressed by the liver and secreted in plasma.

Involvement in disease Anhaptoglobinemia

Sequence similarities Belongs to the peptidase S1 family.
Contains 1 peptidase S1 domain.
Contains 2 Sushi (CCP/SCR) domains.

Cellular localization Secreted.

Images



Representative standard curve using ab157714 Haptoglobin
Mouse ELISA kit

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