

Swine Haptoglobin ELISA Kit ab223136

1 Image

Overview

Product name Swine Haptoglobin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Plasma				5.5%

Inter-assay

Sample	n	Mean	SD	CV%
Plasma				10.3%

Sample type Cell culture supernatant, Urine, Serum, Plasma, Tissue

Assay type Sandwich (quantitative)

Sensitivity 0.4 ng/ml

Range 2 ng/ml - 30 ng/ml

Recovery 97 %

Assay time 4h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Pig

Does not react with: Mouse, Rat, Rabbit, Horse, Cow, Human, Monkey

Product overview

The Swine Haptoglobin ELISA (Enzyme-Linked Immunosorbent Assay) Kit (ab223136) is designed for detection of swine haptoglobin in plasma, serum, urine, and cell culture samples.

This assay employs a quantitative sandwich enzyme immunoassay technique that measures swine haptoglobin in approximately 4 hours. A polyclonal antibody specific for swine haptoglobin has been pre-coated onto a 96-well microplate with removable strips. Swine haptoglobin in standards and samples is sandwiched by the immobilized antibody and a biotinylated polyclonal antibody specific for swine haptoglobin, which is recognized by a streptavidin-peroxidase conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Notes	Haptoglobin (HP, zonulin) is a plasma protein with hemoglobin-binding capacity and a plasma glycoprotein that forms a stable complex with hemoglobin to aid the recycling of heme iron.
Platform	Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Swine Haptoglobin Antibody	1 x 120µl
Anti-Swine Haptoglobin coated Microplate (12 x 8 wells)	1 unit
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml
Swine Haptoglobin Standard	1 vial

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.

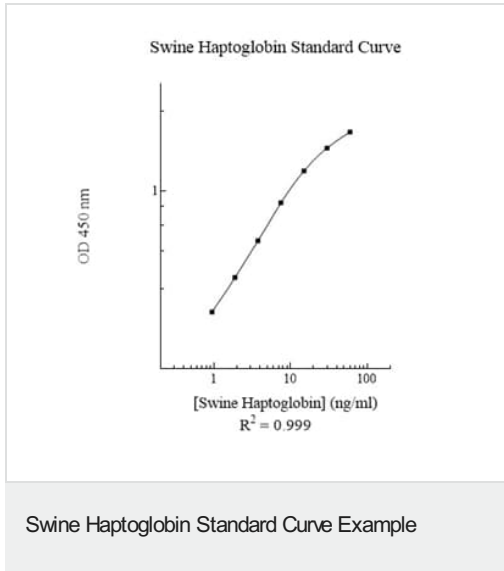
Tissue specificity Expressed by the liver and secreted in plasma.

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



Example of Swine Haptoglobin standard curve

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