

Mouse Hemopexin ELISA Kit (HPX) ab157716

★★★★★ 1 Abreviews 4 References 1 Image

Overview

Product name Mouse Hemopexin ELISA Kit (HPX)

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, Other biological fluids

Assay type Sandwich (quantitative)

Sensitivity 4.5092 ng/ml

Range 12.5 ng/ml - 400 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 Abcam's Hemopexin (HPX) Mouse ELISA kit is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of Hemopexin in biological samples of mice.

In this assay the Hemopexin present in samples reacts with the anti-Hemopexin antibodies which have been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti- Hemopexin antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound Hemopexin. 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 Hemopexin in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Hemopexin in the test sample. The quantity of Hemopexin 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 Hemopexin antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Mouse Hemopexin Calibrator (lyophilized)	1 vial
Mouse Hemopexin ELISA Microplate	1 unit
Stop Solution	1 x 12ml

**Function** Binds heme and transports it to the liver for breakdown and iron recovery, after which the free hemopexin returns to the circulation.

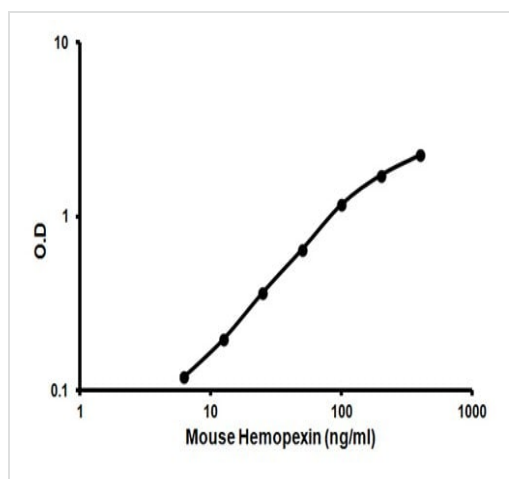
**Tissue specificity** Expressed by the liver and secreted in plasma.

**Sequence similarities** Belongs to the hemopexin family.  
Contains 5 hemopexin-like domains.

**Post-translational modifications** N- and O-glycosylated. O-glycosylated with core 1 or possibly core 8 glycans.

**Cellular localization** Secreted.

## Images



Representative standard curve using ab157716 Hemopexin Mouse ELISA Kit.

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