

Rat Hemopexin ELISA Kit (HPX) ab157734

[1 References](#) [1 Image](#)

Overview

Product name Rat 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 1.1 ng/ml

Range 6.25 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:** Rat

Product overview

Abcam's Hemopexin (HPX) Rat ELISA Kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring Hemopexin (HX) in biological samples of rats.

In this assay the Hemopexin present in samples reacts with the anti-Hemopexin antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-HX antibodies conjugated with horseradish peroxidase (HRP) are added. These enzyme-labeled antibodies form complexes with the previously bound HX. 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 HX in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of HX in the test sample. The quantity of HX in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

ab157734 has been redeveloped with new goat capture and detector antibodies, Kits purchased prior to 2020 used chicken antibodies.

Platform Microplate

Properties

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

Components	1 x 96 tests
100X HRP-conjugated anti-rat Hemopexin antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Rat Hemopexin Calibrator (lyophilized)	1 vial
Rat IgA ELISA Microplate	1 x 12 units
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.

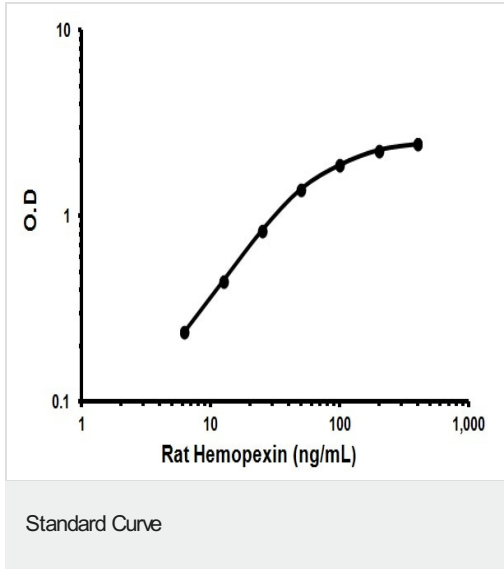
Tissue specificity Expressed by the liver and secreted in plasma.

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 ab157734 Hemopexin Rat 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