

Mouse Myoglobin ELISA Kit ab157722

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

Overview

Product name Mouse Myoglobin 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 4.48 ng/ml

Range 9.375 ng/ml - 300 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 Myoglobin Mouse ELISA kit is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of Myoglobin in mouse samples.

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

Get better reproducibility in only 90 minutes with Mouse Myoglobin ELISA Kit ([ab210965](#)) from our SimpleStep ELISA® range.

Platform

Microplate

Properties

Storage instructions

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

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

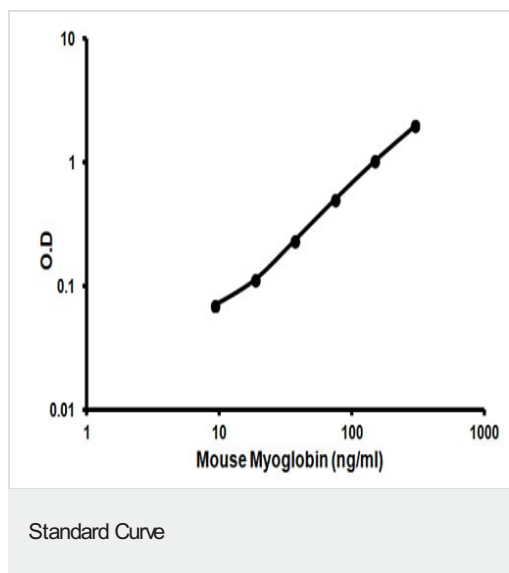
Function

Serves as a reserve supply of oxygen and facilitates the movement of oxygen within muscles.

Sequence similarities

Belongs to the globin family.

Images



Representative standard curve using ab157722 Myoglobin 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