

Rat Myoglobin ELISA Kit ab157739

1 References 1 图像

概述

产品名称	大鼠Myoglobin ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Overall				< 10%
	批次间				
	样品	n	Mean	SD	CV%
	Overall				< 10%
样品类型	Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	3.365 ng/ml				
范围	6.25 ng/ml - 200 ng/ml				
回收率	特定样本回收率				
	样品类型	平均%		范围	
	Serum	> 85		% - %	
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Rat				
产品概述	Abcam's Myoglobin Rat ELISA Kit is a two-site enzyme-linked immunosorbent assay (ELISA) for				

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.

**平台** Microplate

**性能**

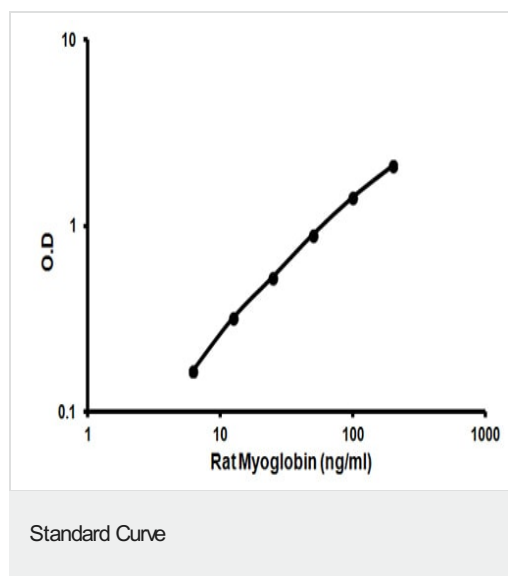
**存放说明** Please refer to protocols.

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

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

**序列相似性** Belongs to the globin family.

**图片**



Representative standard curve using ab157739 Myoglobin 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.cn/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors