

Human CKMB ELISA Kit ab193696

3 References 1 Image

Overview

Product name	Human CKMB ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Serum, Plasma, Cell culture media				
Assay type	Sandwich (quantitative)				
Sensitivity	0.3 ng/ml				
Range	0.3 ng/ml - 80 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		97.17		71% - 123%
	Plasma		87.57		78% - 108%
	Cell culture media		86.45		67% - 106%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's CKMB Human ELISA Kit (ab193696) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of human CKMB in serum, plasma, and cell culture supernatants.				

the samples. The wells are washed and biotinylated anti-Human CKMB antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of CKMB bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

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

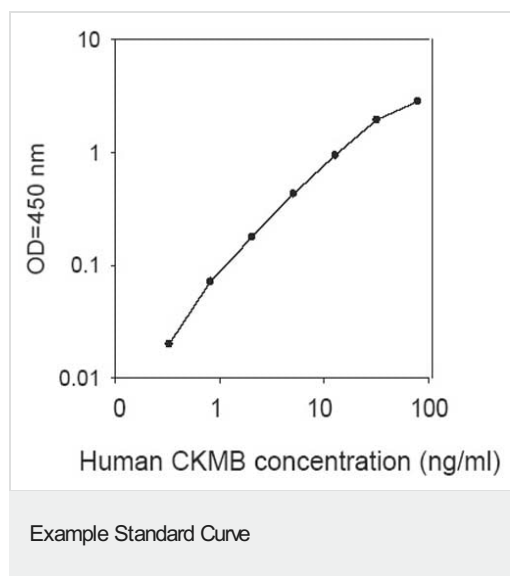
Components	1 x 96 tests
100X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent D	1 x 15ml
5X Assay Diluent B	1 x 15ml
Biotinylated Human CKMB detection antibody	2 vials
Human CKMB Standards (lyophilized)	2 vials
Pre-coated Human CKMB Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Reversibly catalyzes the transfer of phosphate between ATP and various phosphogens (e.g. creatine phosphate). Creatine kinase isoenzymes play a central role in energy transduction in tissues with large, fluctuating energy demands, such as skeletal muscle, heart, brain and spermatozoa.

**Sequence similarities** Belongs to the ATP:guanido phosphotransferase family.  
Contains 1 phosphagen kinase C-terminal domain.  
Contains 1 phosphagen kinase N-terminal domain.

**Cellular localization** Cytoplasm.

## Images



**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