

Product datasheet

Human CKMB ELISA Kit ab193696

[1 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.

This assay employs an antibody specific for Human CKMB coated on a 96-well plate. Standards

and samples are pipetted into the wells and the immobilized antibody captures CKMB present in 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.

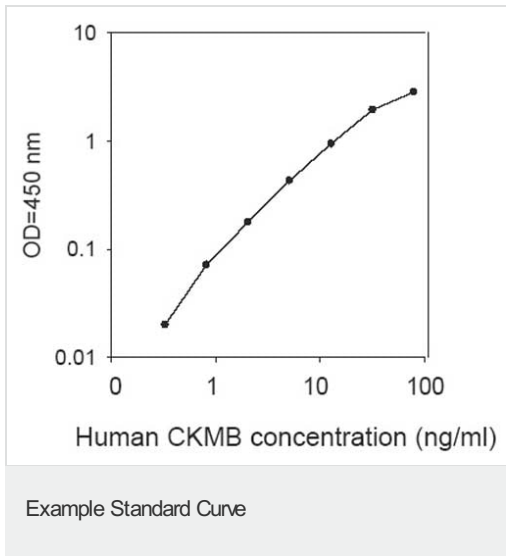
Components	1 x 96 tests
100X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent D	1 x 15ml
5X Assay Diluent B	1 x 15ml
Biotinylated Human CKMB detection antibody (lyophilized)	2 vials
Human CKMB Standards (lyophilized)	2 vials
Pre-coated Human CKMB Microplate (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 phosphagens (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