

Rat CKM ELISA Kit ab187396

SimpleStep ELISA®

7 References 7 Images

Overview

|                    |   |   |           |    |            |
|--------------------|---|---|-----------|----|------------|
| Product name       | Rat CKM ELISA Kit   |   |           |    |            |
| Detection method   | Colorimetric  |   |           |    |            |
| Precision          | Intra-assay   |   |           |    |            |
|                    | Sample  | n | Mean      | SD | CV%        |
|                    | Serum   | 8 |           |    | 4%         |
|                    | Inter-assay   |   |           |    |            |
|                    | Sample  | n | Mean      | SD | CV%        |
|                    | Serum   | 3 |           |    | 5%         |
| Sample type        | Serum, Tissue Extracts, Cit plasma  |   |           |    |            |
| Assay type         | Sandwich (quantitative)   |   |           |    |            |
| Sensitivity        | 5.83 ng/ml  |   |           |    |            |
| Range              | 8.8 ng/ml - 100 ng/ml   |   |           |    |            |
| Recovery           | Sample specific recovery  |   |           |    |            |
|                    | Sample type   |   | Average % |    | Range      |
|                    | Serum   |   | 94        |    | 89% - 97%  |
|                    | Extraction Buffer   |   | 96        |    | 87% - 101% |
| Assay time         | 1h 30m  |   |           |    |            |
| Assay duration     | One step assay  |   |           |    |            |
| Species reactivity | Reacts with: Rat  |   |           |    |            |
| Product overview   | Rat CKM ELISA Kit (ab187396) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CKM protein in serum, tissue extracts, and cit plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Rat CKM with 5.83 ng/ml sensitivity. |   |           |    |            |

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

## Notes

Creatine kinase (CK), also known as creatine phosphokinase (CPK) or phospho-creatine kinase, is an enzyme that catalyzes the reversible transfer of a high-energy phosphate group between phosphocreatine and ATP. There are five known CK isozymes, three located in the cytosol (CKMM, CKMB, CKBB) and two in the mitochondria (Mi<sub>S</sub>CK and Mi<sub>L</sub>CK). Structurally, cytosolic CK is a 86 kDa dimeric enzyme of two identical or non-identical chains, CKM and/or CKB whereas mitochondrial CK is an octameric enzyme associated with mitochondrial membranes. The compartmentation of CK isoenzymes allows for direct association with ATP-providing or consuming processes both linked via metabolite channeling through the creatine phosphate shuttle. Thus CK isoenzymes allow for the generation of a large pool of phosphocreatine, considered to be a temporal and spatial ATP buffering system in tissues with large, fluctuating energy demands, such as skeletal muscle, heart, brain and spermatozoa.

Expression of cytosolic isoenzyme patterns differ between tissues, with skeletal muscle expressing high levels of CK-MM (98%) and low levels of CK-MB (1%), the myocardium (heart muscle) expressing CK-MM at 70% and CK-MB at 25–30% and the brain and neural tissue expressing mainly CK-BB. CKM is also found in serum from normal animals and levels vary based on muscle mass, age, physical activity and underlying pathologies.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

## Platform

Microplate

## Properties

### Storage instructions

Please refer to protocols.

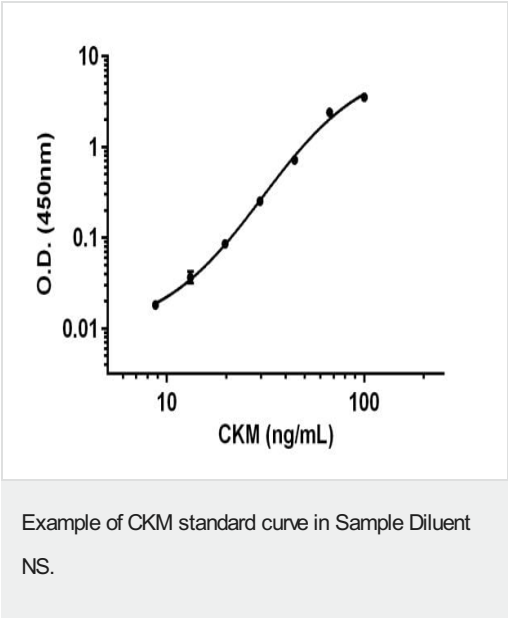
| Components                    | 1 x 96 tests |
|-------------------------------|--------------|
| 10X CKM Capture Antibody      | 1 x 600µl    |
| 10X CKM Detector Antibody     | 1 x 600µl    |
| 10X Wash Buffer PT (ab206977) | 1 x 20ml     |
|                               |              |

| Components  | 1 x 96 tests |
|---|--------------|
| 50X Cell Extraction Enhancer Solution (ab193971)    | 1 x 1ml      |
| 5X Cell Extraction Buffer PTR (ab193970)            | 1 x 10ml     |
| Antibody Diluent 5BI                                | 1 x 6ml      |
| Plate Seals   | 1 unit       |
| Rat CKM Lyophilized Purified Protein                | 2 vials      |
| Sample Diluent NS (ab193972)                        | 1 x 50ml     |
| SimpleStep Pre-Coated 96-Well Microplate (ab206978) | 1 unit       |
| Stop Solution                                       | 1 x 12ml     |
| TMB Development Solution                            | 1 x 12ml     |

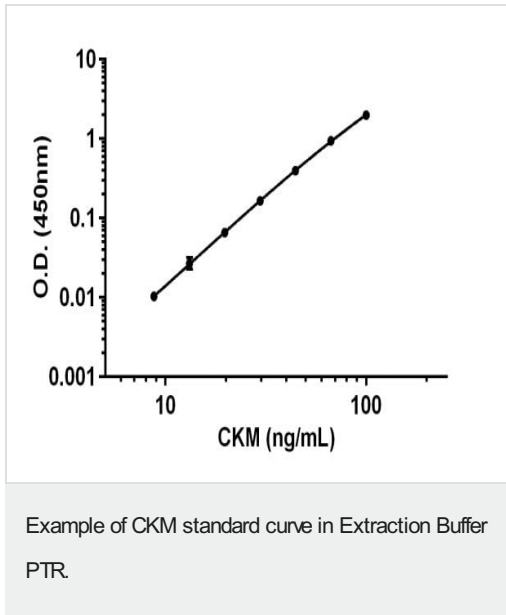
Cellular localization

Cytoplasmic

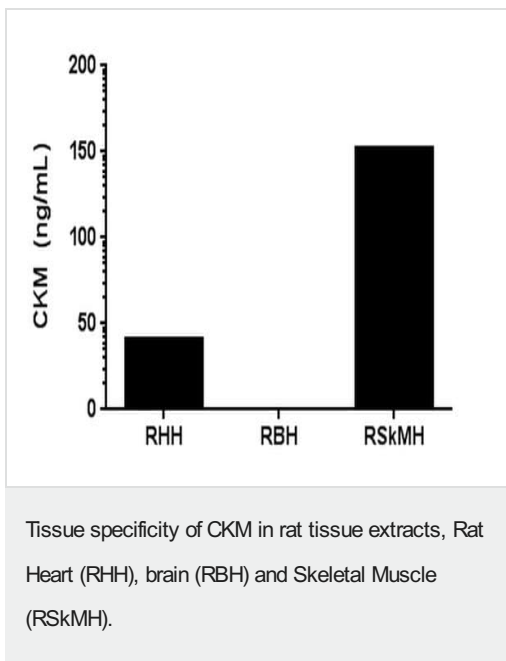
Images



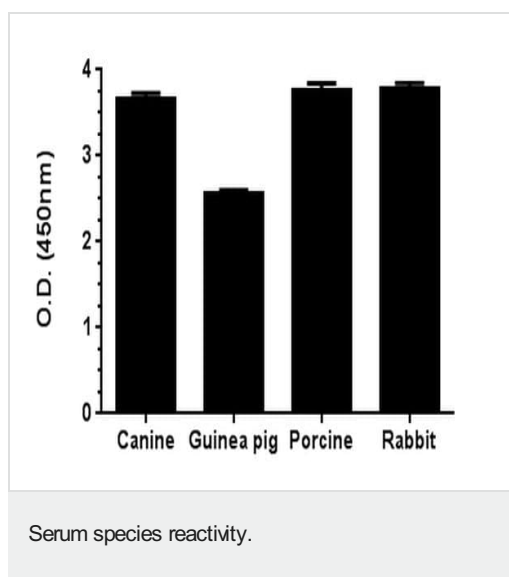
Background-subtracted data values (mean +/- SD) are graphed.



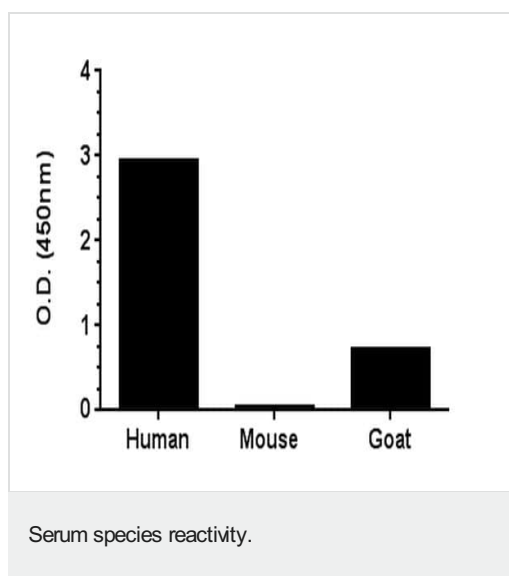
Background-subtracted data values (mean +/- SD) are graphed.



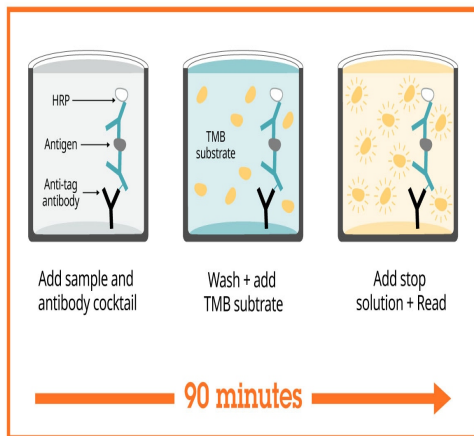
Levels of CKM in 10  $\mu$ g/mL extract were interpolated from a standard curve in 1X Cell Extraction Buffer PTR.



Canine, Guinea pig, Porcine and Rabbit serum was loaded in the assay at 2.5% in Sample Diluent NS. Absorbance levels after background subtraction are displayed.



Human, Mouse and Goat serum was loaded in the assay at 1% in Sample Diluent NS. Absorbance levels after background subtraction are displayed.



Sandwich ELISA - Rat CKM ELISA Kit (ab187396)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

#### Get more done with SimpleStep ELISA



**Easy to use**  
Single-wash 90-minute protocol



**Flexible**  
Matched antibody pairs available



**Precision antibodies**  
High sensitivity, specificity and reproducibility



**Scalable**  
Now in 10-pack and 384-well formats

Sandwich ELISA - Rat CKM ELISA Kit (ab187396)

To learn more about the advantages of SimpleStep ELISA® kits see [here](#).

**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