

### Canine KIM-1 ELISA Kit **ab205084**

★★★★★ [1 Abreviews](#) [1 Image](#)

#### Overview

**Product name** Canine KIM-1 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** Urine, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 0.034 ng/ml

**Range** 0.031 ng/ml - 1 ng/ml

**Recovery** > 85 %

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	0.031ng/ml - 1ng/ml

**Assay time** 3h 30m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Dog

**Product overview** Kidney Injury Molecule-1 in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Kidney Injury Molecule-1 protein in dog biological fluids.

**Notes** In this assay the KIM-1 present in samples reacts with the anti-KIM-1 antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, the Detection Antibody, biotin conjugated anti-KIM-1, is added and complexes are formed. Following a wash step, the horseradish peroxidase (HRP) conjugated Streptavidin is

added and complexes are formed. After another washing step, the complexes are 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 KIM-1 in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of KIM-1 in the test sample. The quantity of KIM-1 in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

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

## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X Detection Antibody Conjugate	1 x 150µl
100X HRP Streptavidin	1 x 150µl
1X Diluent Concentrate	1 x 60ml
20X Wash Buffer Concentrate	1 x 50ml
Anti-Dog KIM-1 ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Dog KIM-1 Calibrator (Lyophilized)	1 vial
Stop Solution	1 x 12ml

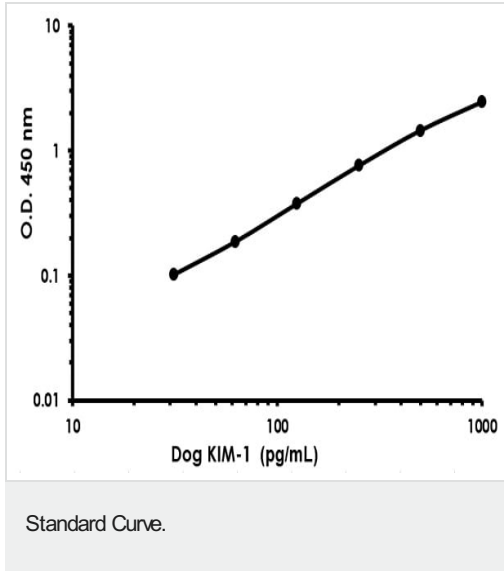
**Function** May play a role in T-helper cell development and the regulation of asthma and allergic diseases. Receptor for TIMD4 (By similarity). In case of human hepatitis A virus (HHAV) infection, functions as a cell-surface receptor for the virus. May play a role in kidney injury and repair.

**Tissue specificity** Widely expressed, with highest levels in kidney and testis. Expressed by activated CD4+ T-cells during the development of helper T-cells responses.

**Sequence similarities** Belongs to the immunoglobulin superfamily. TIM family. Contains 1 Ig-like V-type (immunoglobulin-like) domain.

**Cellular localization** Membrane.

## Images



Representative standard curve using ab205084 Dog KIM-1 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