

# KIM-1 (human) ELISA kit

**Ultra sensitive (1.279 pg/ml) ELISA kit, enabling reduced input sample and matrix interference, for the quantification of KIM-1, an early biomarker for kidney injury or disease.**

The KIM-1 (human) ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of human KIM-1 in urine with results in <2 hours.

Chronic kidney disease (CKD) is a major public health problem, as one in nine American adults have CKD. There have been numerous studies focused on this disease over the years but there remains a need for improved therapeutics and identifying and/or predicting patient outcomes. One protein that has been identified as playing an important role in kidney disease is Kidney Injury Molecule-1 (KIM-1). KIM-1 is a 30KDa, type 1 membrane protein with an ectodomain that contains immunoglobulin and highly O-glycosylated mucin subdomains as well as multiple N-glycosylation sites. It is the most highly upregulated protein in the proximal tubule of the injured kidney. It exists in very low levels in normal kidneys but when upregulated during injury, it is detectable in urine in a wide variety of human diseases. KIM-1 is a potential biomarker for renal injury, which would suggest it has great importance in various kidney diseases and disorders, such as chronic kidney disease (as mentioned above), as well as acute tubular necrosis and acute kidney failure.

Citations: 26

[View Online »](#)

## Ordering Information

[Order Online »](#)

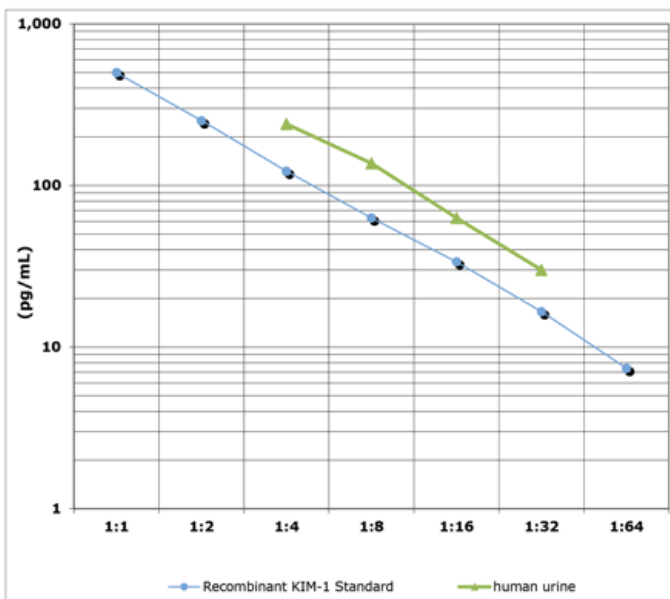
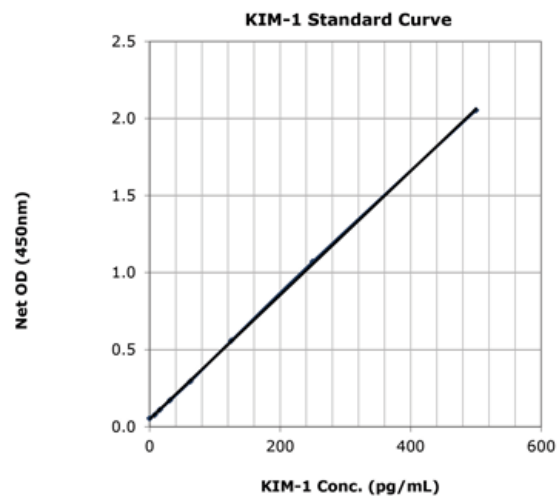
ADI-900-226-0001

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive – reliably measure as little as 1.279 pg/ml of KIM-1
- High Throughput – results in <2 hours from up to 38 samples in duplicate
- Specific – low cross-reactivity to similar proteins
- Quantitative – fully quantitative results for both normal and disease state samples that surpass semi-quantitative Western blot analysis



## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

### Alternative Name

Kidney injury molecule-1, T-cell immunoglobulin and mucin domain-containing protein 1, TIMD-1, TIM-1, HAVCR1, Hepatitis A virus cellular receptor 1, T-cell immunoglobulin mucin receptor 1, CD365

### Application

Colorimetric detection, ELISA

### Application Notes

For the quantitative determination of human KIM-1 in urine.

### Compatibility

This product is compatible with the [Absorbance 96 Plate Reader](#).

### Contents

Microtiter Plate, Assay Buffer, Standard, Antibody-conjugate, Streptavidin-HRP, TMB Substrate, Stop Solution, Wash Buffer Concentrate

### Crossreactivity

No cross reactivity to similar molecules (< 0.02% TIM-3, TIM-4).

### Sensitivity

1.279 pg/ml (range 7.813 – 500 pg/ml)

### Species Reactivity

Human

### Technical Info / Product Notes

This product is licensed from the General Hospital Corporation and is protected by US patents 6,664,385 B1, 7,041,290 B2 and 7,696,321.

### UniProt ID

Q96D42

### Wavelength

450 nm



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.