

KIM-1 (human) monoclonal antibody (3F4)

Kidney Injury Molecule-1 (KIM-1) is a type 1 transmembrane protein, with an immunoglobulin and mucin domain, whose expression is markedly up-regulated in the proximal tubule in the post-ischemic rat kidney. KIM-1 is a potential biomarker for renal injury or renal inflammation. It is a traditional blood and urine marker for the diagnosis of various renal diseases.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Ordering Information

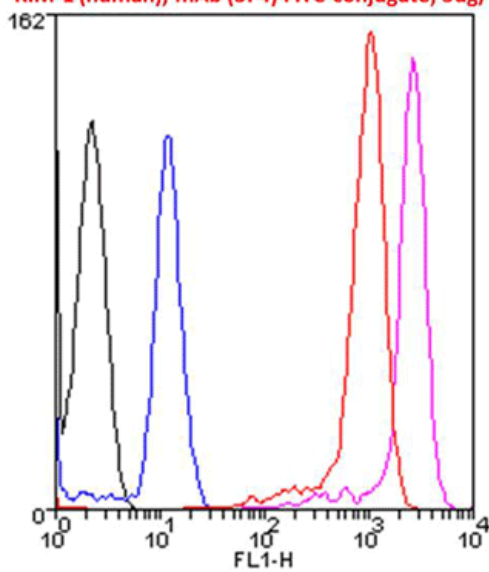
[Order Online »](#)

ADI-905-906-0100	100µg
------------------	-------

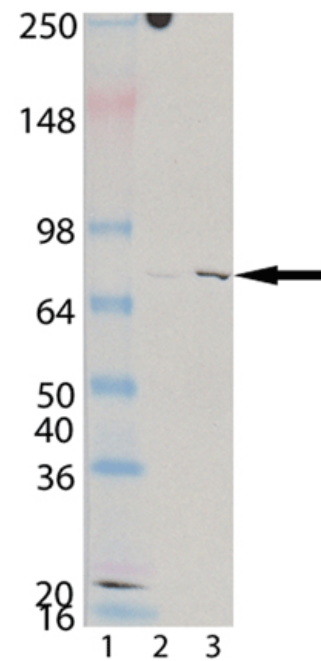
Manuals, SDS & CofA

[View Online »](#)

CaKi-2, Unstain
 ADI-SAB-600FI, IgG1 isotype control, 10ug/mL
 KIM-1 (human), mAb (3F4) FITC-conjugate, 25ug/mL
 KIM-1 (human), mAb (3F4) FITC-conjugate, 5ug/mL



Flow cytometry analysis of 10^6 CaKi-2 cells stained using KIM-1 (human), mAb (3F4), FITC Conjugate at concentrations of 25 μ g/ml and 5 μ g/ml.



Western blot analysis of KIM-1 (human), mAb (3F4) (Prod. No. ADI-905-906): Lane 1: MW marker, Lane 2: HeLa, (cell lysate) (Prod. No. ADI-LYC-HL100), Lane 3: CaKi-2 (cell lysate).

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Kidney Injury Molecule-1, HAVcr-1, TIMD-1, TIM-1, Hepatitis A virus cellular receptor 1 homolog, T cell immunoglobulin and mucin domain-containing protein 1, T-cell immunoglobulin mucin receptor, CD365

Application ELISA, Flow Cytometry, WB

Clone 3F4

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Host Mouse

Immunogen Recombinant human KIM-1.

Isotype IgG1

Purity Detail Protein G-affinity purified.

Source Purified from mouse ascites.

Species Reactivity Human

UniProt ID Q96D42.2

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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-be@enzolifesciences.com)

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

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

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

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