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 upregulated 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

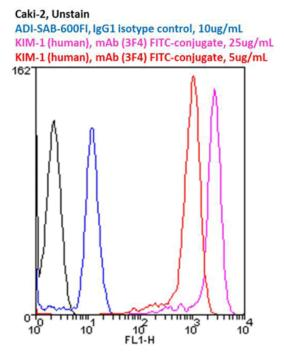
Order Online »

ADI-905-906-0100

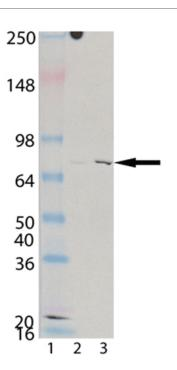
100µg

Manuals, SDS & CofA

View Online »



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



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.

lsotype lgG1

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com