VEGFR-2 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

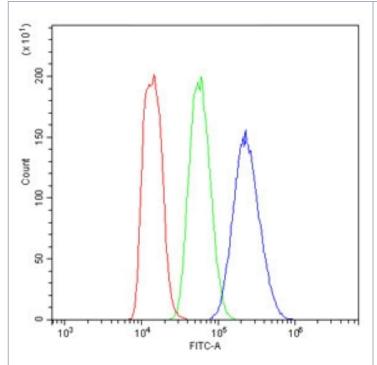
Order Online »

ENZ-ABS927-0100

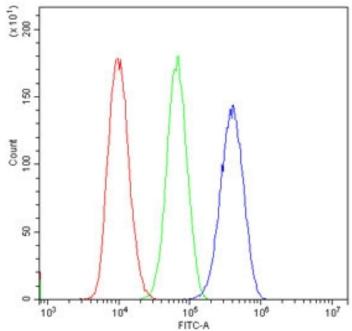
100µg

Manuals, SDS & CofA

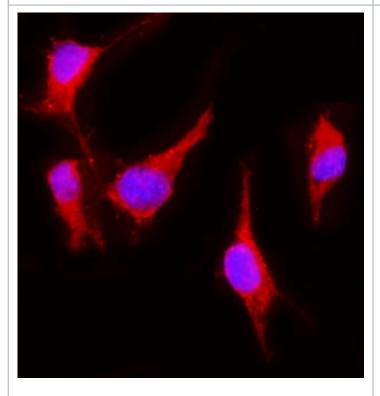
View Online »



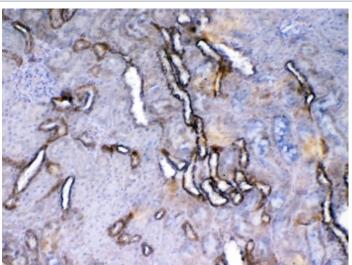
Flow cytometry testing of human MCF7 cells with Vegf receptor 2 antibody at 1µg/million cells (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= Vegf receptor 2 antibody.



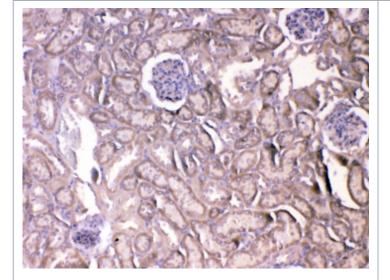
Flow cytometry testing of human LLC1 cells with Vegf receptor 2 antibody at 1µg/million cells (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= Vegf receptor 2 antibody.



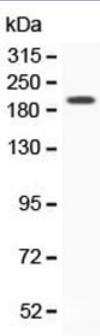
Immunofluorescent staining of FFPE human NIH 3T3 cells with Vegf receptor 2 antibody (red) and DAPI nuclear stain (blue). HIER: steam section in pH6 citrate buffer for 20 min.



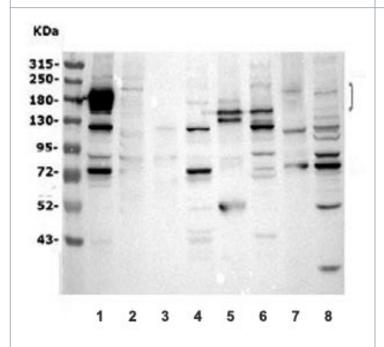
IHC testing of FFPE rat kidney tissue with Vegf receptor 2 antibody at $1\mu g/ml$. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to testing.



IHC testing of FFPE mouse kidney tissue with Vegf receptor 2 antibody at 1µg/ml. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to testing.



Western blot testing of recombinant mouse Vegfr2 protein with Vegf receptor 2 antibody at 0.5µg/ml.



Western blot testing of mouse 1) heart, 2) thymus, 3) spleen, 4) kidney, 5) brain, 6) lung, 7) HEPA1-6 and 8) NIH 3T3 lysate with Vegf receptor 2 antibody. Predicted molecular weight: ~152 (immature), 180-200 kDa (intermediate) and 220-230 kDa (mature).

Handling & Storage

Use/Stability After reconstitution, the Tissue Factor antibody can be stored for up to one month at

+4°C.

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Vascular endothelial growth factor receptor 2, Kdr, FLK-1,

Fetal liver kinase 1, CD309

Application Flow Cytometry, IF, IHC (PS), WB

Formulation Lyophilized from 1X PBS containing 2% Trehalose and

0.025% sodium azide.

Host Rabbit

Immunogen Recombinant mouse VEGFR-2 (aa 20-244).

Isotype IgG

Purity Detail Antigen affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1-3µg/million cells)Immunofluorescence

(FFPE) (2-4μg/ml)Immunohistochemistry (FFPE) (1-2μg/ml)Western blot (0.5-1μg/ml)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Mouse, Rat

UniProt ID P35918

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: July 12, 2024



eu@enzolifesciences.com