

VEGFR-2 polyclonal antibody

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

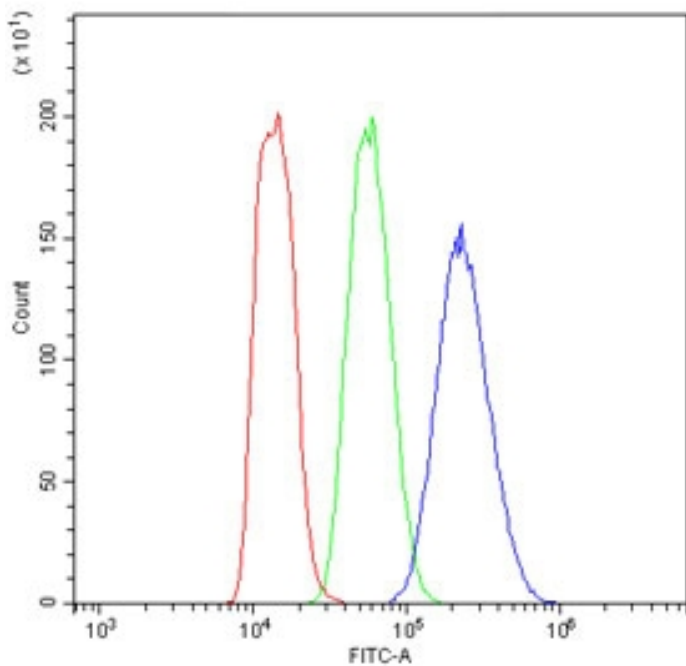
Ordering Information

[Order Online »](#)

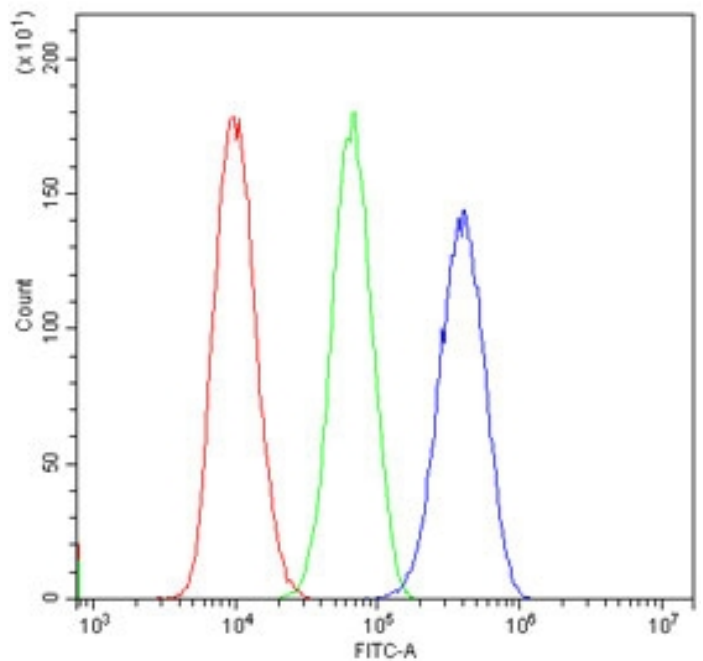
ENZ-ABS927-0100	100µg
-----------------	-------

Manuals, SDS & CofA

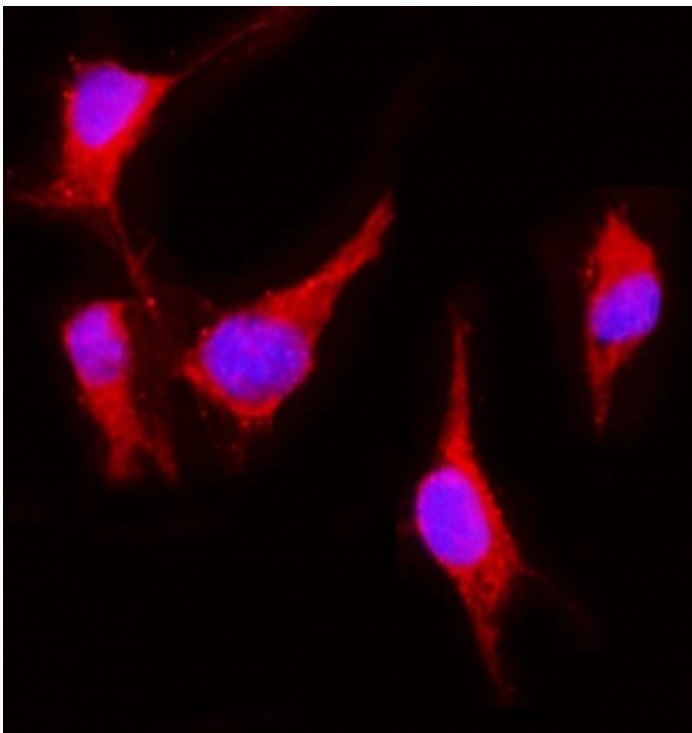
[View Online »](#)



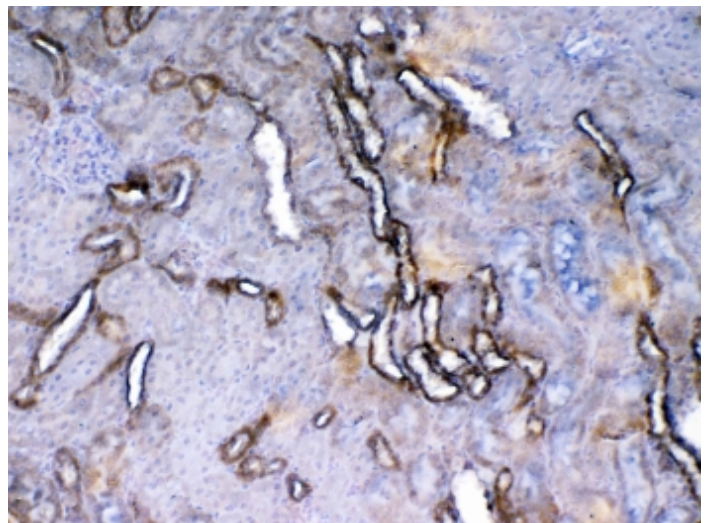
Flow cytometry testing of human MCF7 cells with Vegf receptor 2 antibody at 1 μ g/million cells (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= Vegf receptor 2 antibody.



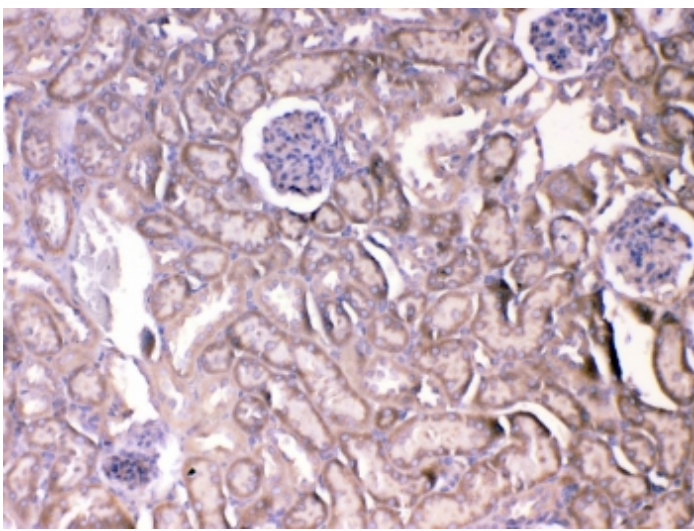
Flow cytometry testing of human LLC1 cells with Vegf receptor 2 antibody at 1 μ g/million cells (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= Vegf receptor 2 antibody.



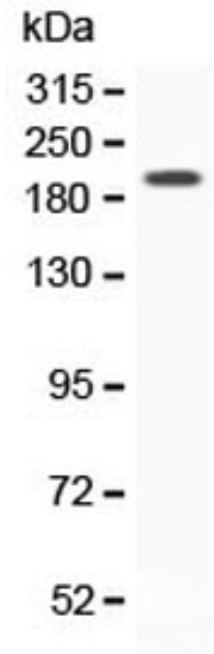
Immunofluorescent staining of FFPE human NIH 3T3 cells with Vegf receptor 2 antibody (red) and DAPI nuclear stain (blue). HIER: steam section in pH6 citrate buffer for 20 min.



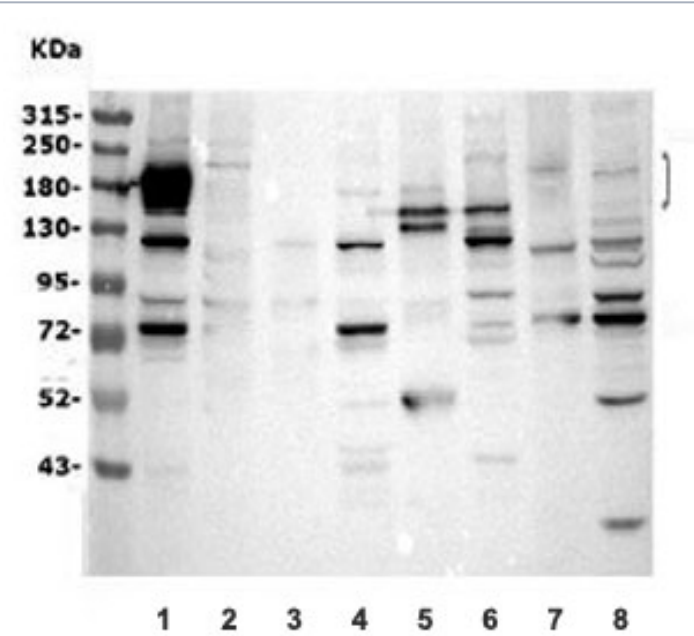
IHC testing of FFPE rat kidney tissue with Vegf receptor 2 antibody at 1 μ g/ml. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to testing.



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



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-
be@enzolifesciences.com](mailto:info-be@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)