

PDGF receptor α polyclonal antibody

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

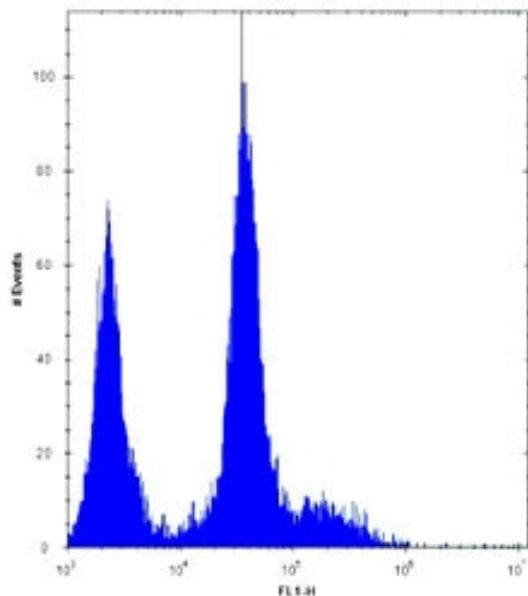
Ordering Information

[Order Online »](#)

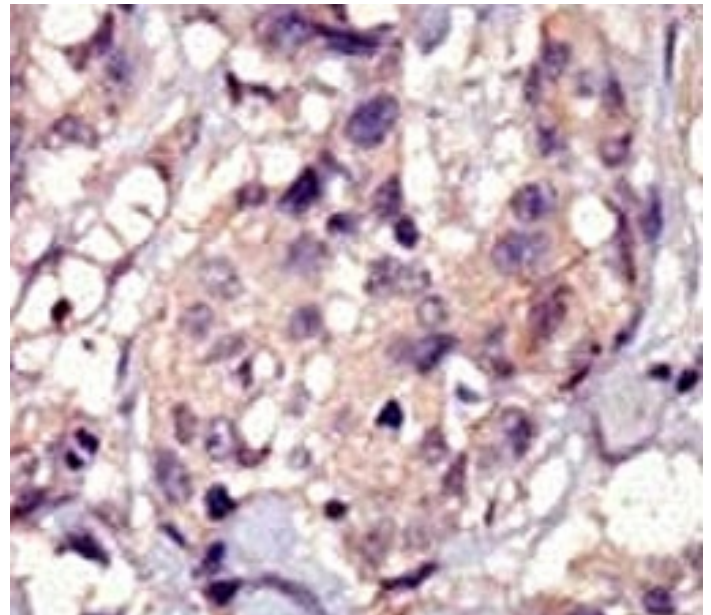
ENZ-ABS910-0400	0.4ml
-----------------	-------

Manuals, SDS & CofA

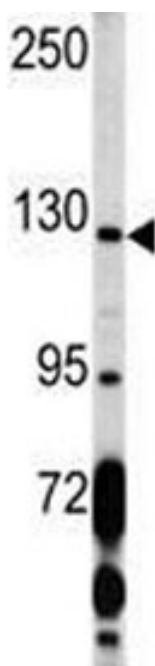
[View Online »](#)



PDGFRA antibody flow cytometric analysis of U251 cells (right histogram) compared to a negative control (left histogram). FITC-conjugated goat-anti-rabbit secondary Ab was used for the analysis.



IHC analysis of FFPE human breast carcinoma tissue stained with PDGFRA antibody



Western blot analysis of PDGFRA antibody and K562 lysate. Predicted molecular weight: 120-195 kDa.



Western blot analysis of PDGFRA antibody and mouse lung tissue lysate. Predicted molecular weight: 120-195 kDa.

Handling & Storage

Handling	Aliquot the PDGFRA antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PDGFRA, Platelet-derived growth factor receptor α , PDGFR-2, CD140a
Application	ELISA, Flow Cytometry, IHC, WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human PDGF receptor α (aa 1048-1077).
Isotype	Ig
Purity Detail	Purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:10-1:50)Immunohistochemistry (paraffin sections) (1:50-1:100)Western blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	P16234

Worry-free Guarantee

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

.



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

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

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

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

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

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

Last modified: July 12, 2024