

HGF/SF monoclonal antibody (7-2)

HGF/SF (Hepatocyte growth factor/Scatter factor) is a multifunctional heterodimeric polypeptide. It mediates the growth and scattering of various cell types, epithelial mesenchymal transition, the formation of tubules and lumens, and promotes angiogenesis. The ligand-receptor pair has also shown to be uniquely involved in most human solid tumors and to participate in tumor development, invasion, and metastasis.

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

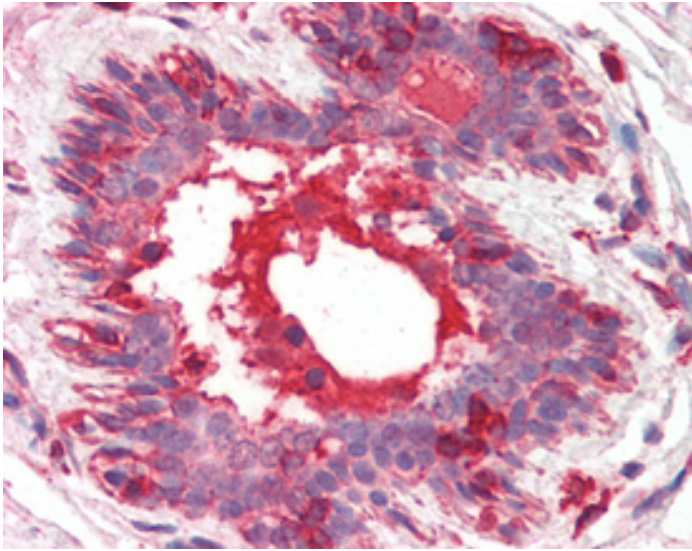
Ordering Information

[Order Online »](#)

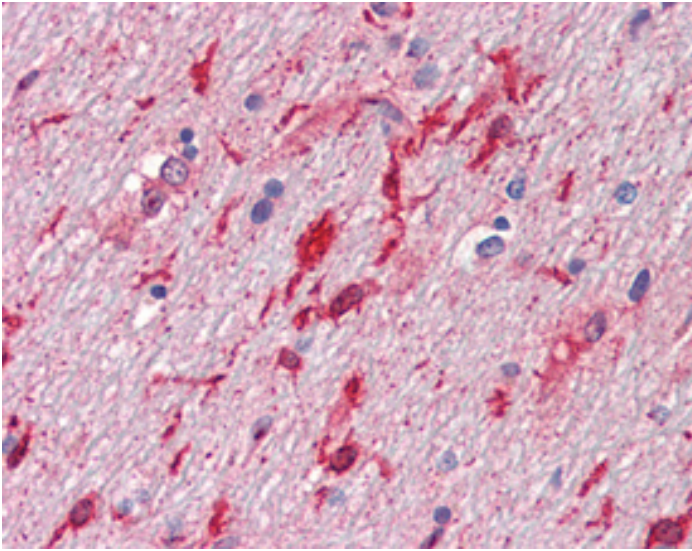
ADI-905-165-100	100µg
-----------------	-------

Manuals, SDS & CofA

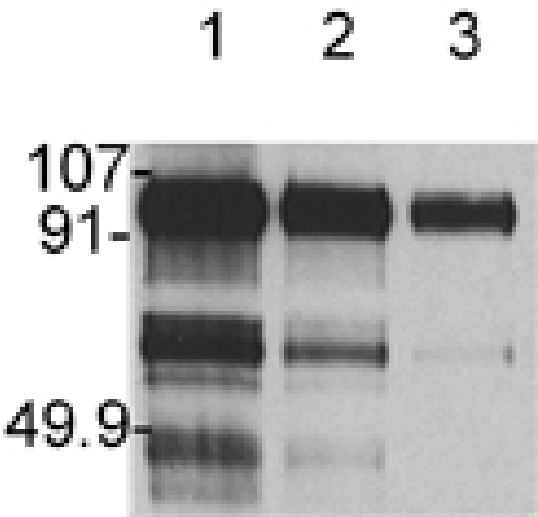
[View Online »](#)



IHC Analysis: Human breast tissue stained with HGF/SF, mAb (7-2) at 10µg/ml.



IHC Analysis: Human brain, cerebellum, white matter tissue stained with HGF/SF, mAb (7-2) at 10µg/ml.



Western blot analysis: Lane 1 hHGF 1µg, Lane 2 hHGF 0.5µg and Lane 3 hHGF 0.25µg.

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	HGFB, HPTA, hepatocyte growth factor
Application	IHC (PS), WB
Clone	7-2
Formulation	Liquid. In PBS containing 0.1% sodium azide.
Host	Mouse
Immunogen	Human Hepatocyte Growth Factor/Scatter Factor.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (2µg/ml colorimetric, 2µg/ml ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Mouse
UniProt ID	P14210
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-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)