

HER3 polyclonal antibody

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

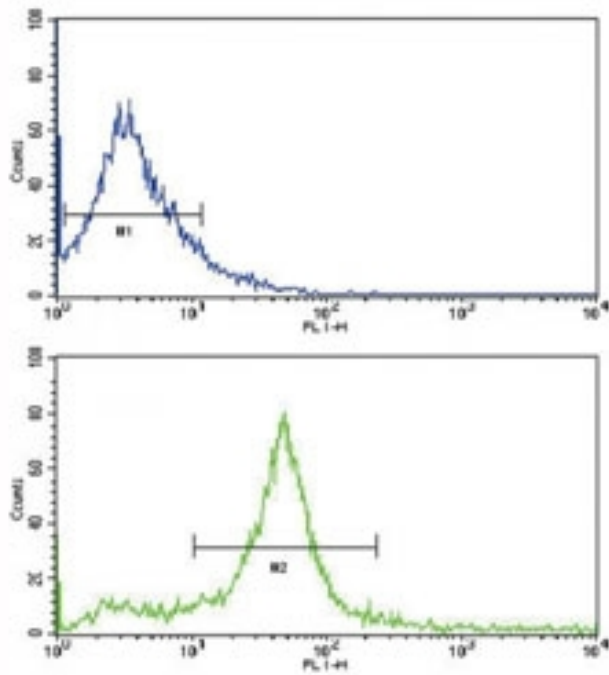
Ordering Information

[Order Online »](#)

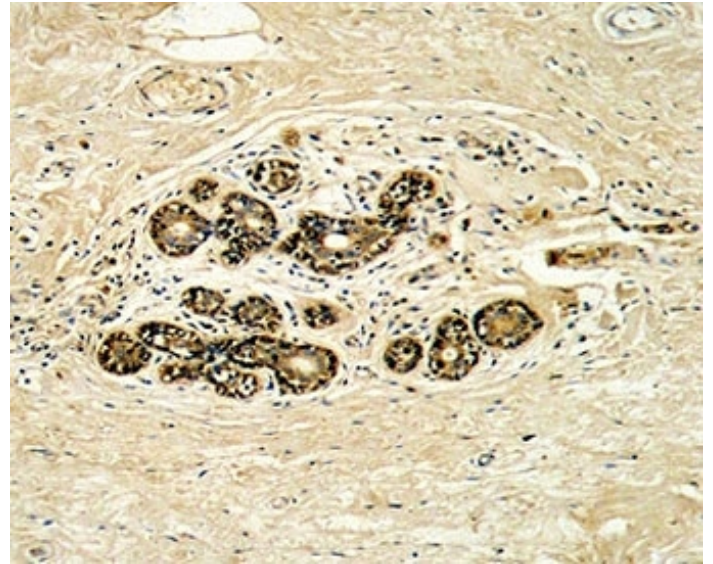
ENZ-ABS893-0400	0.4ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of 293 cells using HER3 antibody (bottom histogram) compared to a negative control (top histogram). FITC-conjugated goat-anti-rabbit secondary Ab was used for the analysis.



IHC analysis of FFPE human breast carcinoma with HER3 antibody



Western blot analysis of HER3 antibody and mouse brain lysate. Predicted molecular weight: 140-180 kDa.

Handling & Storage

Handling	Aliquot the HER3 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	Receptor tyrosine-protein kinase erbB-3
Application	ELISA, Flow Cytometry, IHC, WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human HER3 (aa 24-55).
Isotype	Ig
Purity Detail	Purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:10-1:50)Immunohistochemistry (paraffin sections) (1:10-1:50)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	P21860
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-
fr@enzolifesciences.com](mailto:info-fr@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)