

HER2/Neu monoclonal antibody

Monoclonal antibody to HER2 for use with IHC applications

The HER2/Neu (c-erbB-2) proto-oncogene is a trans-membrane receptor tyrosine kinase that is clinically indicated in a number of carcinomas. Overexpression of the HER2/Neu protein has been associated with ductal breast cancer, as well as pulmonary and gastric adenocarcinomas. A correlation between HER2/Neu and p53 has also been documented, as overexpression of both proteins has been associated with early invasion and metastasis in bladder cancer.

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

Citations: 1

[View Online »](#)

Ordering Information

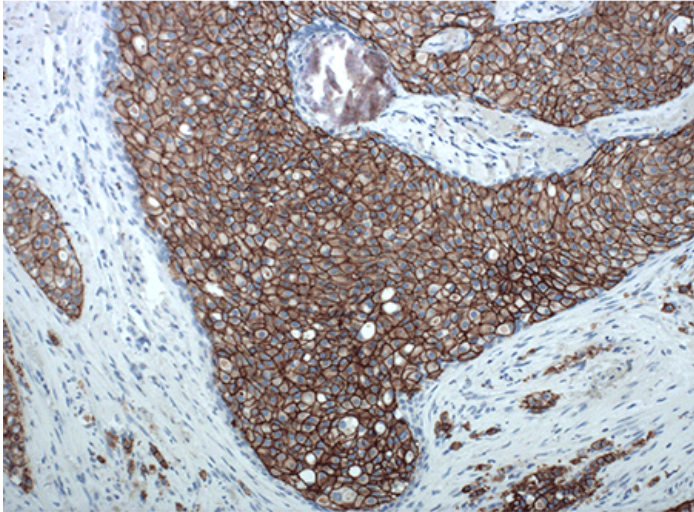
[Order Online »](#)

ENZ-ABS381-1000	1ml
ENZ-ABS381-0100	100µl

Manuals, SDS & CofA

[View Online »](#)

- Strong, clear staining with minimal background
- Cost-effective option compared to competitor's antibody



HER2+ breast cancer tissue incubated with HER2/Neu monoclonal antibody (ENZ-ABS381) and then stained with POLYVIEW® PLUS (anti-Mouse) HRP-DAB Kit (ENZ-KIT160).

Handling & Storage

Use/Stability	When stored correctly, the antibody is stable until the date indicated on the label. To ensure proper stability of the antibody after each run, replace the cap and immediately place the bottle in a refrigerator in an upright position.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Proto-oncogene c-ErbB-2, CD340
Application	IHC
Application Notes	Validated for use with formalin-fixed paraffin embedded (FFPE) samples. For IHC applications: Dilute the concentrate as appropriate in Antibody Blocking Buffer (ENZ-ACC108), and then apply 150 µL to the slide, as indicated in the user's IHC protocol.
Formulation	Liquid. In Tris buffer, pH 7.5, containing 1% BSA and <0.1% sodium azide.
Host	Mouse
Immunogen	Recombinant HER2.
Isotype	IgG
Purity Detail	Affinity-purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:100-1:200)Optimal conditions must be determined individually for each application.
Species Reactivity	Human



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

P04626

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Worry-free Guarantee

This antibody is covered by our **Worry-Free Guarantee.**