# 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**

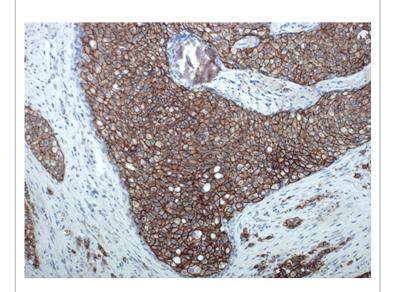
Order Online »

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

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  $\mu L$  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

UniProt ID P04626

Last modified: May 29, 2024



info-

eu@enzolifesciences.com