C-erbB-2 (human) clonal antibody (A24-V)

This antibody is covered by our Worry-Free Guarantee.

Citations: 8

View Online »

Ordering Information

Order Online »

ALX-810-227-L001

1ml

Manuals, SDS & CofA

View Online »

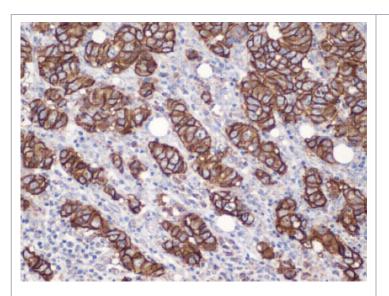


Figure 1: Formalin-fixed and paraffin-embedded human breast ductal carcinoma tissue (4μm) stained with C-erbB-2 (human), cAb (A24-V) (Prod. No. ALX-810-227) shows strong membranous immunostaining of neoplastic cells.(Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph. D. from Clinical Pathology Presov, Ltd., Presov, Slovakia)

Handling & Storage

Use/Stability Stable for at least 12 months after receipt when stored as recommended.

Handling After opening, prepare aliquots and store at +4°C.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ErbB2, HER-2, NEU proto-oncogene, CD340

Application IHC (FS), IHC (PS)

Clone A24-V

Formulation Liquid. In 20mM TRIS/HCl, pH 8.0, containing 20mg/ml

BSA and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the C-terminal region

of human C-erbB-2.

Isotype IgG

Purity Detail Major clone of IgGs obtained from immunoaffinity purified

immunoglobulins corresponding to immunogenic peptide.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 1:100-

1:300)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

Technical Info / Product Notes

Clonal Rabbit Antibody is a pure homogenous fraction of rabbit immunoglobulin (IgG) which corresponds to a unique linear epitope on the target protein. This epitope is designed after a detailed analysis of the antigen structure, regulatory post-translational modifications, protein/protein interaction domain and protein folding. The chosen sequence (peptide) is synthesized and the rabbit is immunized. The clone is then separated from the crude antiserum by "in vitro cloning technology". The resulting product is monospecific and characterized by superior specificity, affinity and avidity.

UniProt ID

P04626

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



eu@enzolifesciences.com