

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 »](#)

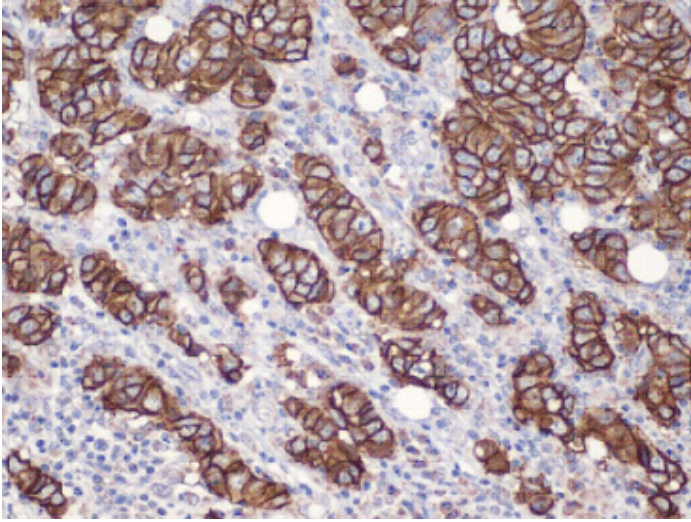


Figure 1: Formalin-fixed and paraffin-embedded human breast ductal carcinoma tissue (4 μ m) stained with C-erbB-2 (human), cAb (A24-V) (Prod. No. ALX-810-227) shows strong membranous immunostaining of neoplastic cells.(Kindly performed and provided by Katarína Poliaková, MD and Lubomír Straka, MD, Ph.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



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)