

# N-cadherin polyclonal antibody

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

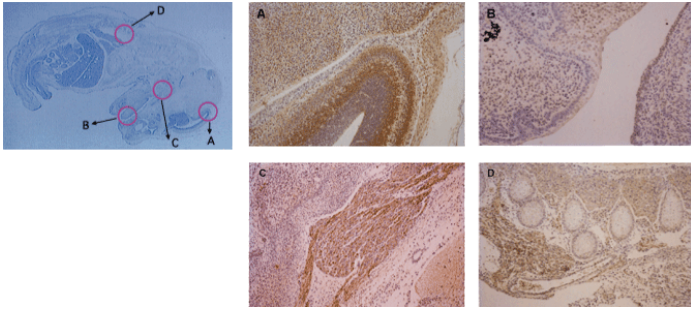
## Ordering Information

[Order Online »](#)

ALX-210-618-C400	400µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Immunohistochemistry of mouse embryo (15 days) paraffin embedded tissues using PAb to N-cadherin (Prod. No. ALX-210-618).

**Method:**Antigen : microwave treatment (3 min.×10 times) (10 mM citrate buffer, pH 6.0)

Reaction of the primary antibody : 10 µg/ml, 1 hr

Detection : Envision plus (Dako)

Colorimetric substrate : DAB

# Handling & Storage

Use/Stability	Stable for at least 2 years after receipt when stored at +4°C. Reconstituted antibody is stable for 6 months when stored at +4°C and for 1 year when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	CD325, Cadherin-2, Neural cadherin
Application	IHC (FS), IHC (PS), WB
Application Notes	Does not inhibit N-cadherin dependent cell-cell contact.
Crossreactivity	Does not cross-react with human E- or P-cadherin.
Formulation	Lyophilized from 10mM PBS, pH 7.4, containing 1% BSA. Does not contain preservatives.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 808-827 of human N-cadherin.
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (frozen and paraffin sections, 5-10µg/ml)Western Blot (5-10µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Reconstitution**

Reconstitute with 200µl distilled water. TBS should be used for dilution as PBS may cause precipitation. Dilute with 20mM TBS, pH 7.4, containing 1% BSA and 0.1% sodium azide and 10mM calcium chloride. In case of immunohistochemistry on paraffin or frozen sections, Ca<sup>2+</sup> is not necessary.

**Species Reactivity**

Bovine, Chicken, Human, Mouse

**UniProt ID**

P19022

**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)