

# $\beta$ -Catenin (NT/exon 2) monoclonal antibody (7D11)

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

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-0600	100 $\mu$ g
--------------	-------------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable at -80°C up to 1 year, at 4°C up to 3 months.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Application</b>	ELISA, IHC, IP, WB
<b>Clone</b>	7D11
<b>Formulation</b>	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human $\beta$ -catenin.
<b>Isotype</b>	IgG2a
<b>Positive Control</b>	Included. (Prod. No. ALX-840-027)
<b>Purity Detail</b>	Thiophilic adsorption and size exclusion chromatography purified.
<b>Recommendation Dilutions/Conditions</b>	Western Blot (0.5 $\mu$ g/ml for HRPO/ECL detection)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Reconstitution</b>	Reconstitute with 1ml water (15 minutes at room temperature).
<b>Species Reactivity</b>	Dog, Human, Mouse

**Specificity**

Recognizes the N-terminus (exon 2) of human, mouse, rat (weak) and dog  $\beta$ -catenin.

**UniProt ID**

P35222

**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)