

α -Catenin monoclonal antibody (15D9)

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

Citations: 22

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-101-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Reconstituted antibody is stable for 1 year when stored at -80°C. Thaw aliquots at +37°C. Thawed aliquots may be stored at +4°C for up to 3 months.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ICC, IP, WB
Clone	15D9
Crossreactivity	Does not cross-react with β -catenins.
Formulation	Lyophilized from 1ml 2x PBS, 0.09% sodium azide, PEG and sucrose.
Host	Mouse
Immunogen	Recombinant human α -catenin.
Isotype	IgG1
Positive Control	Included. (Prod. No. ALX-840-026)
Purity Detail	Purified from serum-free cell culture supernatant by subsequent thiophilic adsorption and size exclusion chromatography.
Recommendation Dilutions/Conditions	Immunocytochemistry (0.1-1 μ g/ml; formaldehyde fixation is tolerated)Western Blot (0.5 μ g/ml for HRPO/ECL detection; recommended blocking buffer CPPT: 10mM TRIS-HCl, pH 7.4, 0.5% (w/v) casein, 1% (w/v) PEG 4,000, 1% (w/v) polyvinylpyrrolidone, 0.1% (v/v) Tween 20, 150mM sodium chloride)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Reconstitution	Reconstitute with 1ml distilled water (15 min., RT).

Species Reactivity Dog, Human, Mouse, Rat

UniProt ID P35221 (α 1), P26232 (α 2), Q9UI47 (α 3)

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



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)