

# β-Catenin ELISA kit

The β-Catenin ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The β-Catenin EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours that measures pan β-Catenin. Absorbance is read at 450 nm. This kit has been validated in-house with human, mouse, and rat cell lysates and correlated Western blot. The β-Catenin EIA kit detects as little 33.8 pg/mL β-catenin surpassing semi-quantitative Western blot analysis and provides simple liquid color-coded reagents that reduce laboratory error.

β-catenin is a 781 amino acid adherens junction protein that plays an important role in two unrelated cellular functions. Adherens junctions are critical for establishment and maintenance of epithelial layers, such as those found lining the surface of organs. Involved in both cell-cell adhesion and in the wnt/wg signaling pathway, β-catenin can be found in the cytoplasm or the nucleus. In the cytoplasm, β-catenin is unstabilized by phosphorylation and association with E-cadherin. When β-catenin translocates to the nucleus, it has comparatively low levels of phosphorylation. Wnt/wg signaling results in β-catenin accumulation and transcriptional activation of specific target genes during development. Deregulation of β-catenin signaling is an important event in the evolution a number of malignancies, including colon cancer, melanoma, ovarian cancer, and prostate cancer. The homology between human, mouse and rat β-catenin is 99.9%.

Citations: 7

[View Online »](#)

## Ordering Information

[Order Online »](#)

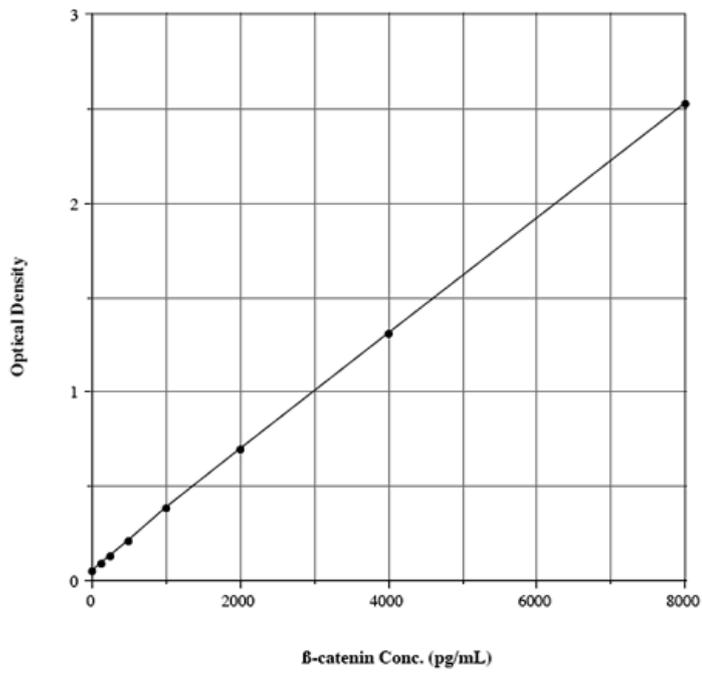
ADI-900-135

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Reliable - fully validated in-house with human, mouse, and rat cell lysates and Western blot correlation
- Sensitive - quantify less than 33.8 pg/ml β-catenin
- Simple - liquid color-coded reagents reduce error



## Handling & Storage

**Use/Stability** Store all components at 4°C, except standard at -20°C.

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cadherin-associated protein

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of  $\beta$ -Catenin in cell lysates of human, mouse, and rat origin.

**Assay Time** 3 hours

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Microtiter plate, Conjugate, Antibody, Assay buffer 13, RIPA Cell lysis buffer 2, Wash buffer concentrate, Standard, TMB Substrate, Stop solution 2

**Sensitivity** 33.8 pg/ml (range 125 – 8,000 pg/ml)

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** P35222 (human), Q02248 (mouse), Q9WU82 (rat)

**Wavelength** 450 nm

**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)