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 »

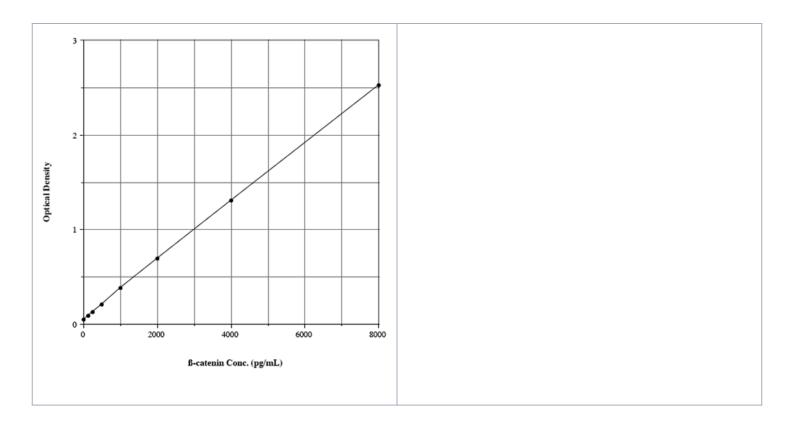
ADI-900-135

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 β -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

