

## Catenin α Protein

Full-length recombinant protein expressed in Sf9 cells

**Catalog # C05-30G**

Lot # U003-2

### Product Description

Recombinant full-length human Catenin α was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_001903](#).

### Gene Aliases

CTNNA1, CAP102, FLJ36832

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

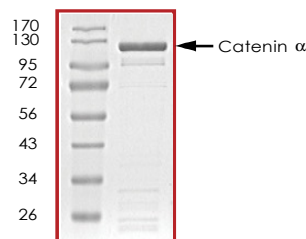
### Scientific Background

Catenin α is a novel actin-binding and bundling protein. Catenin α is responsible for organizing and tethering actin filaments at the zones of E-cadherin-mediated cell-cell contact (1). Monomeric Catenin α can bind strongly to E-Cadherin-β-Catenin, whereas the dimer preferentially binds actin filaments. Different molecular conformations are associated with these different binding states, indicating that Catenin α is an allosteric protein. α-Catenin directly regulates actin-filament organization by suppressing Arp2/3-mediated actin polymerization, likely by competing with the Arp2/3 complex for binding to actin filaments (2)

### References

1. Rimm, D L. et al: Alpha 1(E)-catenin is an actin-binding and -bundling protein mediating the attachment of F-actin to the membrane adhesion complex. Proc Natl Acad Sci U S A. 1995 Sep 12;92(19):8813-7.
2. Drees, F. et al: Alpha-catenin is a molecular switch that binds E-cadherin-beta-catenin and regulates actin-filament assembly. Cell. 2005 Dec 2;123(5):903-15.

### Purity



The purity of Catenin α was determined to be **>95%** by densitometry. Approx. MW **123kDa**.

## Catenin α Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	C05-30G
Specific Lot Number	U003-2
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**