

## Talin VHH Antibody

Z-N008

Support: echelon@echelon-inc.com



### Description:

Talin VHH Antibody was raised in llama and purified from *E. coli* as a VHH single domain antibody with a COOH-terminal HA epitope tag.

### Applications:

Pull-down (PD), Immunoprecipitation (IP), and Enzyme-Linked Immunoassay (ELISA).

Not recommended for Western Blot (WB).

Other in vitro, in vivo, cellular, & tissue applications are possible, but have not been verified by Echelon Biosciences.

### Properties:

**Form** – liquid

**Storage instructions** – Store at -20 °C upon arrival. Aliquot and store at -80 °C if longer storage is necessary. Avoid repeated freeze/thaw cycles.

**Storage buffer** – 20 mM Tris-HCl pH 8.0, 150 mM NaCl, 1mM DTT, 60 % glycerol

**Concentration** – 1.0 mg/mL

**Purity** – >80%

**Immunogen** – purified protein fragment encompassing mouse Talin amino acids 2364-2541

**Clonality** – monoclonal, Nb19

**Isotype** – VHH single domain antibody

**Tag** – C-Terminal HA

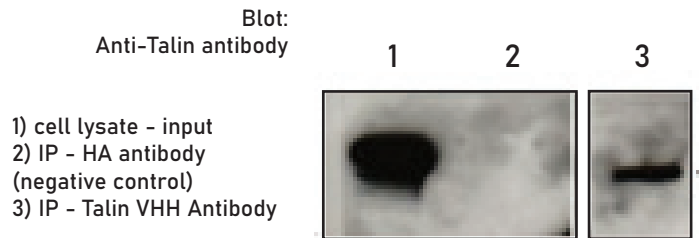
### Specificity:

Z-N008 reacts with mouse and human Talin. Reactivity with other species has not been tested.

### Background:

Talin is a ~270 kDa mechanosensitive cytoskeletal protein that links the actin cytoskeleton to integrins. It is concentrated in cellular regions that contact the extracellular matrix or other cells such as focal adhesions, immune synapses, invadopodia, and podosomes.

### Data: Immunoprecipitation



MDA-MB-231 cell lysate was incubated with HA antibody alone (2) or Talin VHH Antibody (3) followed by SDS-PAGE and WB with a rabbit polyclonal Talin antibody.

### References:

1. Van Audenhove I. et al. (2014) Stratifying fascin and cortactin function in invadopodia formation using inhibitory nanobodies and targeted subcellular fractionation. *FASEB J*; 28(4):1805-1818
2. Van Audenhove I. et al. (2015) Fascin actin bundling controls podosome turnover and disassembly while cortactin is involved in podosome assembly by its SH3 domain in THP-1 macrophages and dendritic cells. *Biochem Biophys Acta*; 1853(5):940-952

### Related Products:

Products	Catalog Number
Assays and Reagents	
PTEN Activity ELISA Kit	K-4700
GloPIP Biotin PI(4,5)P2	C-45B6a
Nanobodies	
Gelsolin VHH Antibody	Z-N005
C-Src VHH Antibody	Z-N007
Vinculin VHH Antibody	Z-N011

Technical Data Sheet Rev. 1a, 08-20-21 - For research use only. Not intended or approved for diagnostic or therapeutic use.