

## Vinculin VHH Antibody

Z-N011

Support: echelon@echelon-inc.com



### Description:

Vinculin 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** – 0.5 mg/mL

**Purity** – >80%

**Immunogen** – purified C-terminal fragment of human Vinculin

**Clonality** – monoclonal, Nb85

**Isotype** – VHH single domain antibody

**Tag** – C-Terminal HA

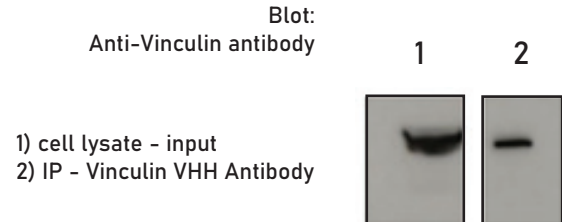
### Specificity:

Z-N011 reacts with human Vinculin. Reactivity with other species has not been tested.

### Background:

Vinculin is a large ~117 kDa cytoskeletal protein that links to integrins and other cytoskeletal proteins such as talin, actinin, and F-actin. Vinculin is a prominent regulator of focal adhesions, adherens junctions, and lamellipodium formation.

### Data: Immunoprecipitation



HEK293 cell lysate was immunoprecipitated with Vinculin VHH Antibody and blotted with a rabbit anti-Vinculin monoclonal antibody.

### References:

1. Wang F, et al. (2021) Enhanced non-toxic immunodetection of alternaria mycotoxin tenuazonic acid based on ferritin-displayed anti-idiotypic nanobody-nanoluciferase multimers. *J Agric Food Chem*; 69(16):4911-4917
2. Van Audenhove I, et al. (2016) Nanobodies as versatile tools to understand, diagnose, visualize, and treat cancer. *EBioMedicine*; 8:40-48

### Related Products:

Products	Catalog Number
Assays and Reagents	
PI(4,5)P2 Mass ELISA Kit	K-4500
Myosin Kinase Substrate	930-30
Nanobodies	
FAK VHH Antibody	Z-N004
C-Src VHH Antibody	Z-N007
Talin VHH Antibody	Z-N008

Technical Data Sheet Rev. 1a, 08-24-21 - For research use only. Not intended or approved for diagnostic or therapeutic use.