

## Ezrin VHH Antibody

Z-N003

Support: echelon@echelon-inc.com



### Description:

Recombinant anti-human Ezrin 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), Western Blot (WB), and Enzyme-Linked Immunoassay (ELISA).

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 – >90%

Immunogen – purified full-length human recombinant protein

Clonality – monoclonal, Nb22

Isotype – VHH single domain antibody

Tag – C-Terminal HA

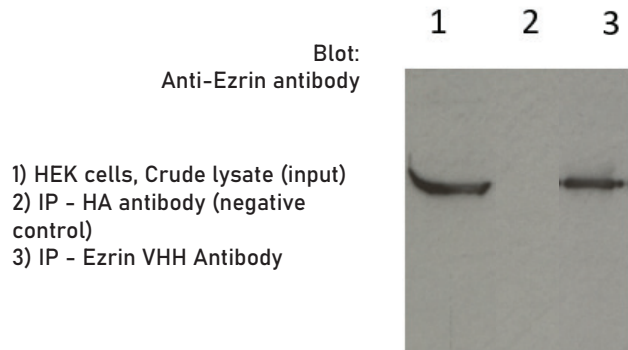
### Specificity:

Z-N003 reacts with human Ezrin.

### Background:

Ezrin (cytovillin, villin-2) is a member of the closely related family of ERM proteins (Ezrin, Radixin, Moesin). Ezrin interacts with NHERF, phospholipids, actin and CD44, with its N-terminal FERM domain anchoring the protein to the plasma membrane. Ezrin plays a role in wound healing, cell migration/invasion and metastasis.

### Data: Immunoprecipitation



HEK293T cell lysate was incubated with control antibody or Ezrin VHH Antibody followed by SDS-PAGE and WB with a rabbit polyclonal Ezrin antibody.

### References:

1. Beghein E, Gettemans J. (2017) Nanobody Technology: A Versatile Toolkit for Microscopic Imaging, Protein-Protein Interaction Analysis, and Protein Function Exploration. *Front Immunol*; 8:771

### Related Products:

Products	Catalog Number
Assays and Reagents	
PI(4,5)P2 Mass Assay	K-4500
BODIPY FL PI(4,5)P2	C-45F6
Nanobodies	
Pan-Akt VHH Antibody	Z-N001
Talin VHH Antibody	Z-N008
Tensin 4 VHH Antibody	Z-N010