

Tensin 4 VHH Antibody

Z-N010

Support: echelon@echelon-inc.com



Description:

Tensin 4 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 protein fragment, residues 428-714 of human Tensin 4

Clonality – monoclonal, Nb12

Isotype – VHH single domain antibody

Tag – C-Terminal HA

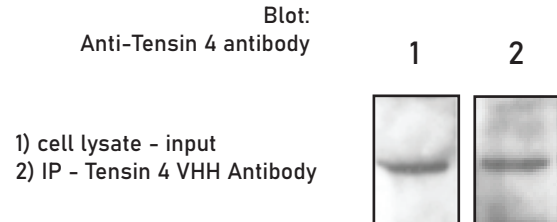
Specificity:

Z-N010 reacts with human Tensin 4. Reactivity with other species has not been tested.

Background:

Tensin 4 (TNS4, CTEN) is a component of the cytoskeleton that connects integrins with actin filaments. TNS4 regulates cell motility, including metastasis in cancer cells, and is aberrantly expressed in prostate, lung and breast cancer.

Data: Immunoprecipitation



HeLa cell lysate was immunoprecipitated with Tensin 4 VHH Antibody and blotted with a mouse anti-Tensin monoclonal antibody.

References:

- Bertier L, et al. (2018) Nanobodies targeting cortactin proline rich, helical, and actin binding regions downregulate invadopodium formation and matrix degradation in SCC-61 cancer cells. *Biomed Pharmacother*; 102:230-241
- 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
SF1670 (PTEN Inhibitor)	B-0350
Nanobodies	
Gelsolin VHH Antibody	Z-N005
C-Src VHH Antibody	Z-N007
Vinculin VHH Antibody	Z-N011

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