

Echelon Biosciences Inc.

c-Src VHH Antibody

Z-N007

Support: echelon@echelon-inc.com



Description:

c-Src 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 human c-Src amino acids 1-252

Clonality – monoclonal, Nb76

Isotype – VHH single domain antibody

Tag – C-Terminal HA

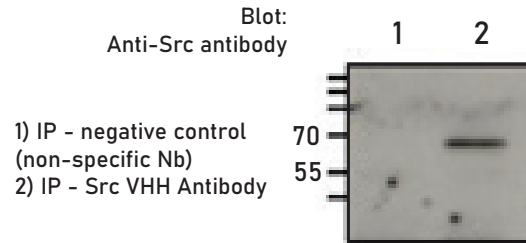
Specificity:

Z-N007 reacts with human c-Src. Reactivity with other species has not been tested.

Background:

Src is a 60 kDa proto-oncogene that codes for a non receptor tyrosine kinase. c-Src phosphorylates numerous substrates controlling their activity or localization. Src overactivation promotes tumor invasion, metastasis, cell cycle progression, and cytoskeletal reorganization.

Data: Immunoprecipitation



HEK293T cell lysate was incubated with either a non-specific nanobody clone or Src-specific VHH antibody followed by SDS-PAGE and WB with polyclonal Src antibody.

References:

1. Beghein E. et al. (2018) Cortactin and fascin-1 regulate extracellular vesicle trafficking or invadopodia formation and function. *Sci Rep*; 8(1):15606
2. Bertier L. et al. (2017) Inhibitory cortactin nanobodies delineate the role of NTA and SH3 domain-specific functions during invadopodium formation and cancer cell invasion. *FASEB J*; 31(6):2460-2476

Related Products:

Products	Catalog Number
Assays and Reagents	
PTEN Activity ELISA Kit	K-4700
LY294002 (PI3-K Inhibitor)	B-0294
Nanobodies	
Annexin VHH Antibody	Z-N002
FAK VHH Antibody	Z-N004
Tensin 4 VHH Antibody	Z-N010

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