

Echelon Biosciences Inc.

Focal Adhesion Kinase VHH Antibody

Z-N004

Support: echelon@echelon-inc.com



Description:

Recombinant Focal Adhesion Kinase 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 suitable 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 – >90%

Immunogen – purified protein fragment encompassing mouse FAK amino acids 750-1053

Clonality – monoclonal, Nb54

Isotype – VHH single domain antibody

Tag – C-Terminal HA

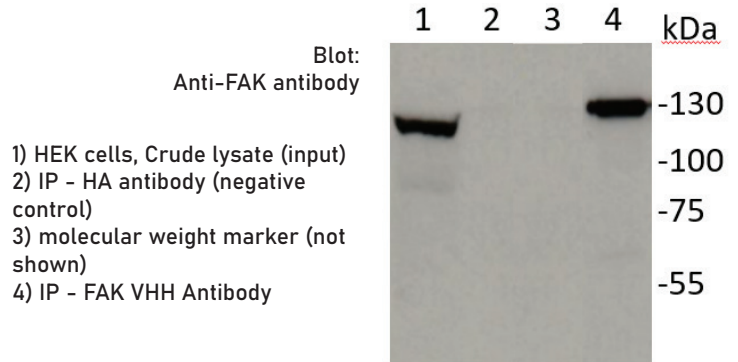
Specificity:

Z-N004 reacts with human and mouse Focal Adhesion Kinase. Other species have not been tested.

Background:

Focal adhesion kinase (FAK) is a 125 kDa cytosolic (non-receptor) tyrosine kinase. FAK is found in most cells promoting cell migration by inducing turn-over of cell contacts with the underlying matrix. During apoptosis, FAK is cleaved by caspase 3 contributing to loss of focal contacts, cell rounding and formation of blebs.

Data: Immunoprecipitation



HEK293T cell lysate was incubated with control antibody or Focal Adhesion Kinase VHH Antibody followed by SDS-PAGE and WB with a rabbit monoclonal FAK antibody.

References:

1. Van Audenhove I, Debeuf N, Boucherie C, Gettemans J. (2015) Fascin actin bundling controls podosome turnover and disassembly while cort-actin is involved in podosome assembly by its SH3 domain in THP-1 macrophages and dendritic cells. *Biochim Biophys Acta*; 1853(5):940-952.

Related Products:

Products	Catalog Number
Assays and Reagents	
PTEN Activity ELISA Kit	K-4700
PI3-Kinase Activity ELISA	K-1000S
Nanobodies	
Annexin VHH Antibody	Z-N002
C-Src VHH Antibody	Z-N007
Tensin 4 VHH Antibody	Z-N010

Technical Data Sheet Rev. 1a, 08-13-21 - For research use only. Not intended or approved for diagnostic or therapeutic use.