

## Pan-Akt VHH Antibody

Z-N001

Support: echelon@echelon-inc.com



### **Description:**

Pan-Akt VHH Antibody was raised in llama and purified from E. coli cells 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.

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** – human, full length recombinant p53 purified from insect cells

DNA-binding domain (p53 DBD)

**Clonality** – monoclonal, Nb139

**Isotype** – VHH single domain antibody

**Tag** – C-Terminal HA

### **References:**

1. Merckaert, T., O. Zwaenepoel, K. Gevaert, and J. Gettemans (2020) Development and characterization of protein kinase B/AKT isoform-specific nanobodies. PLoS One. 15:e0240554.
2. Merckaert, T., O. Zwaenepoel, K. Gevaert, and J. Gettemans (2021) An AKT2-specific nanobody that targets the hydrophobic motif induces cell cycle arrest, autophagy and loss of focal adhesions in MDA-MB-231 cells.

### **Related Products:**

Products	Catalog Number
Assays and Reagents	
PTEN Activity ELISA	K-4700
PI(3,4)P2 diC8	P-3408
Nanobodies	
Annexin V VHH Antibody	Z-N002
FAK VHH Antibody	Z-N004
Tks5 VHH Antibody	Z-N009

### **Specificity:**

Z-N001 interacts with human Akt. Reactivity with other species has not been tested.

### **Background:**

Akt kinases are three highly homologous (Akt 1, 2, 3) isoforms of serine/threonine kinases. Akt, also known as protein kinase b (PKB) is frequently over-activated in cancer and considered a therapeutic target.

