

PHLPP2 Protein

Recombinant human protein expressed in E. coli cells

Catalog # P72-21G

Lot # F469-4

Product Description

Recombinant human PHLPP2 (766-1043) was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_015020](#).

Gene Aliases

KIAA0931; PHLPL

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

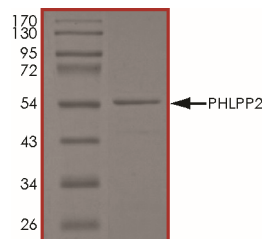
Scientific Background

PHLPP2 or PH domain and leucine rich repeat protein phosphatase 2 is the important regulators of Akt serine-threonine kinases and conventional/novel protein kinase C (PKC) isoforms. PHLPP1 and PHLPP2 differentially regulated AKT signaling by selectively dephosphorylating the hydrophobic motifs of AKT2 or AKT3. PHLPP may act as a tumor suppressor in several types of cancer due to its ability to block growth factor-induced signaling in cancer cells.

References

1. Brognard, J. et.al: PHLPP and a second isoform, PHLPP2, differentially attenuate the amplitude of Akt signaling by regulating distinct Akt isoforms. *Molec. Cell* 25: 917-931, 2007.

Purity



The purity of PHLPP2 was determined to be **>95%** by densitometry. Approx. MW **56 kDa**.

PHLPP2 Protein

Recombinant human protein expressed in E. coli cells

Catalog #	P72-21G
Lot #	F469-4
Purity	>95%
Concentration	0.1µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.