

FANCL Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog # F289-30G

Lot # J691-2

Product Description

Recombinant full-length human FANCL was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_018062](#).

Gene Aliases

FAAP43; PHF9; POG

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

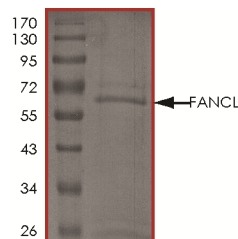
Scientific Background

FANCL or Fanconi anemia complementation group L is a member of the Fanconi anemia complementation group. Fanconi anemia is a genetically heterogeneous recessive disorder characterized by cytogenetic instability, hypersensitivity to DNA crosslinking agents, increased chromosomal breakage, and defective DNA repair. FANCL is crucial in the FA pathway as the catalytic subunit required for monoubiquitination of FANCD2 (1). UBE2W interacts with FANCL and regulates the monoubiquitination of Fanconi anemia protein FANCD2 (2).

References

- Meetei.et.al: A novel ubiquitin ligase is deficient in Fanconi anemia. *Nature Genet.* 35: 165-170, 2003.
- Zhang, Y.et.al: UBE2W interacts with FANCL and regulates the monoubiquitination of Fanconi anemia protein FANCD2. *Molec. Cells* 31: 113-122, 2011.

Purity



The purity of FANCL was determined to be **>75%** by densitometry. Approx. MW **67 kDa**.

FANCL Protein

Full length human recombinant protein expressed in Sf9 cells

Catalog Number	F289-30G
Specific Lot Number	J691-2
Purity	>75%
Concentration	0.1 µg/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.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.