

CBL Protein

Recombinant human protein expressed in E. coli cells

Catalog # C272-381G

Lot # D2399-5

Product Description

Recombinant human CBL (1-375) was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM_005188](#).

Gene Aliases

C-CBL; CBL2; FRA11B; NSLL; RNF55

Formulation

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

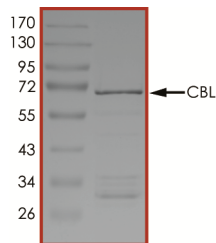
Scientific Background

CBL proto-oncogene acts as a RING finger E3 ubiquitin ligase, transfers ubiquitin from specific E2 ubiquitin-conjugating enzymes to targeting substrates for degradation by the proteasome(1). The N-terminal phosphotyrosine binding domain of CBL allows the interaction with the substrates containing phosphorylated tyrosine including the activated receptor tyrosine kinases, which plays the negative regulation of these pathways (1). Mutation or translocation of CBL is involved in acute myeloid leukemia, Jacobsen syndrome and Noonan syndrome-like disorder (2, 3).

References

1. Howlett C.J., et.al. The proto-oncogene p120 (Cbl) is a downstream substrate of the Hck protein-tyrosine kinase. *Biochem. Biophys. Res. Commun.* 257:129-138, 1999.
2. Muraoka M, et.al. Adults with germline CBL mutation complicated with juvenile myelomonocytic leukemia at infancy. *J Hum Genet. Jun; 61(6): 523-6, 2016.*
3. Martinelli S., et.al. Heterozygous germline mutations in the CBL tumor-suppressor gene cause a Noonan syndrome-like phenotype. *Am. J. Hum. Genet.* 87:250-257, 2010.

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **69kDa**.

CBL Protein

Recombinant human protein expressed in E. coli cells

Catalog #	C272-381G
Lot #	D2399-5
Purity	>85%
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.