

# Cbl-b (casitas B-lineage lymphoma proto-oncogene-b), Full Length

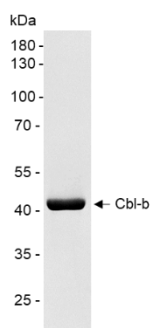
Cat. # UB308FL

<b>Background</b>	Cbl-b is an E3 ubiquitin ligase that regulates a wide range of signaling molecules in the cell. Members of the Cbl protein family share a highly homologous tyrosine kinase-binding (TKB) domain, followed by a conserved helical linker (L) domain and a RING (Really Interesting New Gene) finger (RF) domain, which interact with ubiquitin-conjugating enzymes (E2s). In contrast, the C-terminal regions of these proteins are less conserved. This entry refers to the full-length form of Cbl-b.
<b>Alternate Names</b>	Cbl Proto-Oncogene B, RING Finger Protein 56 (RNF56), Casitas B-Lineage Lymphoma Proto-Oncogene B, RING-Type E3 Ubiquitin Transferase CBL-B, E3 Ubiquitin-Protein Ligase CBL-B, Signal Transduction Protein CBL-B, Cas-Br-M (Murine) Ectropic Retroviral Transforming Sequence B 2
<b>Application(s)</b>	Ubiquitin ligation reactions

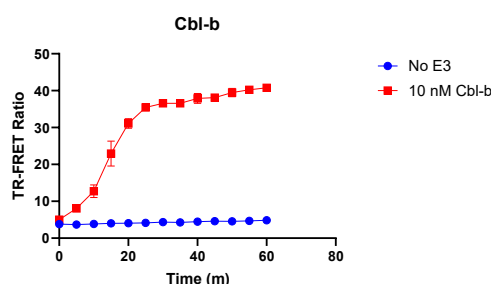
## Product Specifications

<b>Tag</b>	FLAG
<b>Purity</b>	> 95% by RP-HPLC
<b>Molecular Weight</b>	110 kDa
<b>Quantity</b>	25 µg
<b>Species</b>	Human
<b>Expression System</b>	E. coli
<b>Physical State</b>	Liquid
<b>Buffer</b>	50 mM Tris, 150 mM NaCl, 10 mM DTT, 10% glycerol
<b>Activity</b>	A typical enzyme concentration of 1-100 nM is used for <i>in vitro</i> conjugation, depending on experimental conditions
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

## Product QC



**SDS-Page Analysis of purified Cbl-b.** Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



**Activity Assay of Cbl-b.** 10 nM Cbl-b was tested in a TR-FRET assay showing robust E3 ligase activity.

## References

1. Augustin, RC., et al., J Immunother Cancer, 2023. 11(2):e006007.
2. Kumar, J., et al., J Immunother Cancer, 2021. 9(1):e001688.