

## Data Sheet

### **CBL-B Y363F, His, Biotin-labeled**

Human, recombinant, CBL-B Y363F with N-terminal His-tag and biotin-tag

**Catalog #:** 80413

**Lot #:** 150430

**Conc:** 0.47 mg/ml

**Formulated in:** 40 mM HepesI, pH 7.4, 120 mM NaCl, 20% glycerol, 200 mM imidazole

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

**References:**

1. Jeon, M.S., *et al.*, *Immunity*. 2004; **21**(2): 167-177.
2. Chiang, Y.J. *et al.*, *Nature*. 2000; **403**: 216-220.

**Description:** Human Casitas B-lineage lymphoma proto-oncogene b (CBL-B), also known as E3 ubiquitin-protein ligase CBL-B, RING finger protein 56, and SH3-binding protein CBL-B with Y363F mutation GenBank Accession No. NM\_170662, a.a. 39 – 426 with N-terminal His-tag and N-terminal Avi-tag, expressed in an *E. coli* expression system. This protein was enzymatically biotinylated *in vivo* using AviTag™ technology. MW=48 kDa.

**Application:** Useful for the study of enzyme kinetics and regulation, to dephosphorylate target substrates and for screening inhibitors.

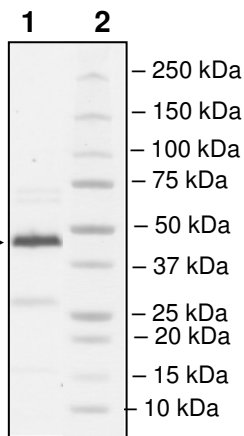
### Quality Assurance

**4-20% SDS-PAGE  
Coomassie staining**

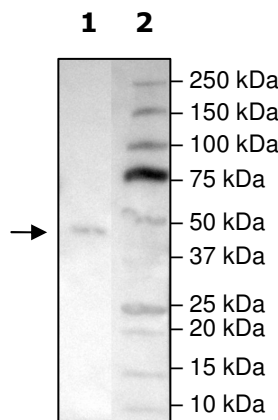
**Lane 1:**  
2 µg CBLB Mutant

**Lane 2:**  
Standard

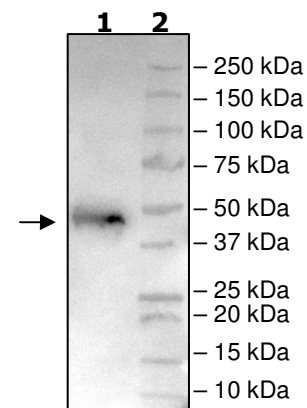
**Purity:** ≥85%  
**MW:** 48 kD



**4-20% SDS-PAGE  
Anti-His Western**



**4-20% SDS-PAGE  
Streptavidin western**



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Sequence

MGSSHHHHHSQDPNSSARLQVDGLNDIFEAQKIEWHEGGAAADRRTVEKTKWLMDKVVRLCQ  
NPKLQLKNSPPYILDILPDYQHLRLILSKYDDNQLAQLSENEYFKIYIDSLMKKSKRAIRLFK  
EGKERMYYEQSODRRNLTKLSLIFSHMLAEIKAIFFPNGQFQGDNFRIKADAAEFWRKFFGDKTI  
VPWKVFRQCLHEVHQISSGLEAMALKSTIDLTCNDYISVFEFDIFTRLFQPWGSILRNWNFLAVT  
HPGYMAFLTYDEVKARLQKYSTKPGSYIFRLSCTRLGQWAIQYVTGDGNILQTI PHNKPLFQALI  
DGSREGFYLYPDGRSYNPDLTGLCEPTPHDHIKVTQEYELFCEMGSTFQLCKICAENDKDVKIE  
PCGHLMCTSLTAWQESDGQGCFFCRCEIKGTEP IIVDPF

**Red:** His-Avi-tag

**Black:** CBLB (a.a. 39-426)

**Yellow:** Y363F

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)