

Data Sheet

CBL-B, His-Avi-Tag

Human, Recombinant, N-terminal His-tag,
N-terminal Avi-tag

Catalog #: 80414

Lot #: 161003 **Conc:** 0.36 mg/ml

Formulated in: 40 mM HEPES pH 7.4, 120 mM NaCl, 20% glycerol.

Stability: At least 6 months at -80°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, 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. 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:

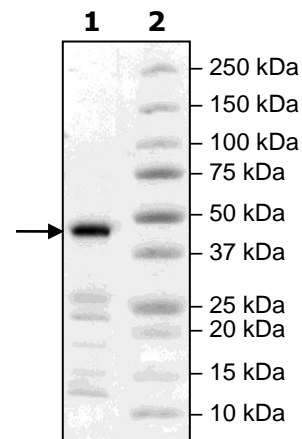
2 μg CBLB

Lane 2:

Protein Marker

MW: 48 kDa

Purity: $\geq 73\%$



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

Please visit our website at: www.bpsbioscience.com