

Data Sheet

CBL-B, His-Avi-Tag, Biotin-Labeled

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

Catalog #: 80412

Lot #: 170821-B

Conc: 1.08 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 3 mM DTT, and 150 mM Imidazole.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

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. This protein was enzymatically biotinylated using Avi-Tag™ technology.

Quality Assurance

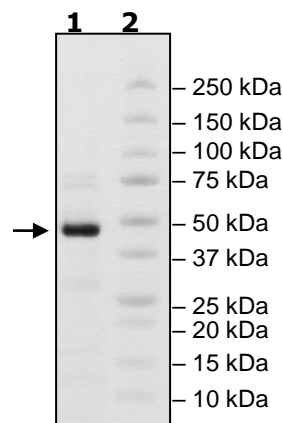
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
CBL-B(39-426)

Lane 2:
Protein Marker

Purity: $\geq 84\%$

MW: 48 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com