

Data Sheet

c-CBL, GST-Tag

Human, Recombinant, N-terminal GST-Tag
Catalog #: 100370
Lot #: 190430 **Conc.:** 0.40 mg/ml

Description:

Human c-CBL, also known as CBL-2, Cbl Proto-Oncogene, Signal Transduction Protein CBL, and Cbl Proto-Oncogene, E3 Ubiquitin Protein Ligase, Genbank Accession No. NM_005188, a.a. 49-434 with N-terminal GST-tag, expressed in an E. coli expression system. MW = 71 kDa.

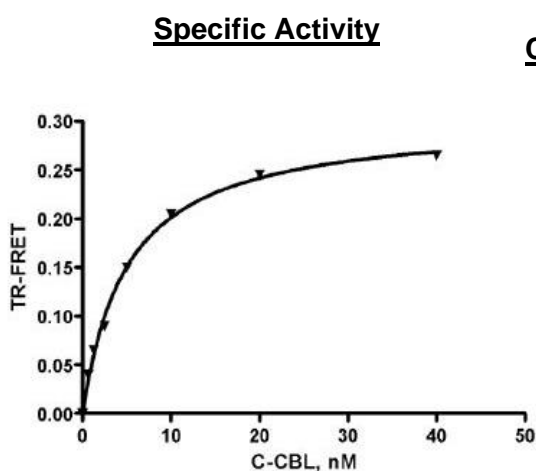
Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 0.04% Tween20, 20% glycerol, 16 mM glutathione, 3 mM DTT.

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

References:

1. Keane, M.M., *Oncogene*, 1999 June; **18**: 3365-3375.
2. Griffiths, E.K., *Molec. and Cellular Bio.* 2003; Oct. **23(21)**:7708-7718

Quality Assurance

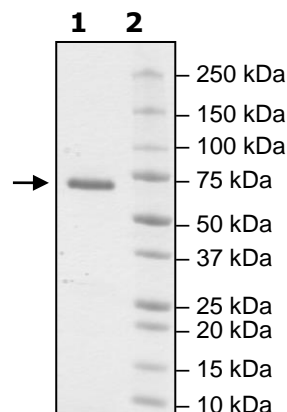


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
c-CBL
Lane 2:
Protein Marker

Purity: $\geq 90\%$

MW: 71 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