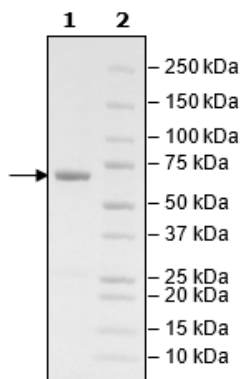


## Product Information

<b>Construct:</b>	CBL-B (Y363E) (GST-39-426)
<b>Mutation:</b>	Y393E
<b>Concentration:</b>	1.14 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 0.2 mM TCEP, and 8 mM glutathione
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_170662
<b>MW:</b>	72 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Assay was performed according to the protocol in CBL-B TR-FRET Assay Kit (BPS Bioscience #79575).
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CBL-B Activity

