

## Recombinant ABL1 protein

**Catalog No:** 31160

**Lot No:** 02606001

**Expressed In:** Baculovirus

**Quantity:** 10 µg

**Concentration:** 0.206 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg of Recombinant ABL1 protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol.

**Protein Details:** Recombinant human ABL1 protein was produced using baculovirus infected Sf9 cells. The protein was against amino acids P118-S535, accession number M14752 and N-terminally fused to GST-Factor Xa cleavage site. Purified by GST-agarose affinity purification.

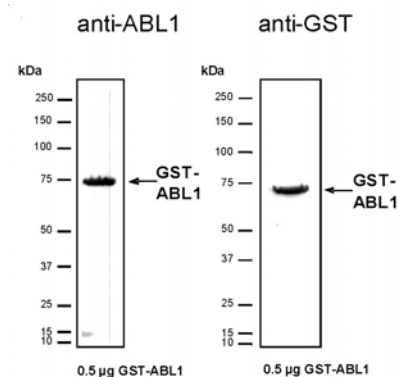
**Application Notes:** Recombinant ABL1 is suitable for kinase assays and Western blot. The molecular weight of the protein is ~74.14 kDa. The activity of the protein is ~ 33 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: PolyAEKY, 5 µg/50 µl, Recombinant ABL1: 50 ng/50 µl.

Kinase activity may vary depending on the substrate and reaction conditions used.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### Recombinant ABL1 protein tested by SDS-PAGE gel.

0.5 µg of ABL1 protein was analyzed using SDS/PAGE followed by Western blot with either an anti-ABL1 or anti-GST antibody.