

# Abl1 (human), (recombinant) (His- tag)

The ABL1 protooncogene encodes a cytoplasmic and nuclear protein tyrosine kinase that has been implicated in cell differentiation, cell division, cell adhesion, and stress responses. The c-Abl protein is inhibited by its SH3 domain, and deletion of this domain turns ABL1 into an oncogene. A chromosomal translocation results in the head-to-tail fusion of the BCR and ABL1 genes. This translocation is present in many cases of chronic myelogenous leukemia. The DNA-binding activity of Abl1 is regulated by CDC2-mediated phosphorylation.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Abelson murine leukemia viral oncogene homolog 1

**Application Notes** Useful tool to study enzyme regulation and kinetics, phosphorylate target substrates, screen for inhibitors.

**Formulation** Liquid. In 50mM sodium phosphate, pH 7.0, containing 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

**MW** ~135kDa

**Purity** ≥70% (SDS-PAGE)

**Sequence** Human recombinant Abl1 (aa 27-1130) with N-terminal histidine tag.

**Source** Produced in insect cells. Produced in a baculovirus expression system.

**UniProt ID** P00519

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)