

Abl1 (mouse), (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.

Concentration 0.1mg/ml

Formulation Liquid. In 50mM TRIS-HCl, pH 7.5, containing 150mM sodium chloride, 0.25mM DTT, 0.1mM EGTA, 0.1mM EDTA, 0.1mM PMSF and 25% glycerol.

MW ~135kDa

Purity ≥70% (SDS-PAGE)

Sequence Mouse recombinant Abl1 (aa 27-1123) with N-terminal histidine tag. Mouse Abl1 is 89% identical (91% similar) to human Abl1.

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

Specific Activity 820 nmol/min/mg using a peptide substrate, EAIYAAPFAKKK, in 25mM MOPS, pH 7.2, containing 12.5mM β-glycerol-phosphate, 25mM MgCl₂, 5mM EGTA, 2mM EDTA. Add 0.25mM DTT to assay buffer prior to use.

UniProt ID P00520



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989

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)

Last modified: May 29, 2024