

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.

Ordering Information

[Order Online »](#)

BML-SE562-0005	5µg
----------------	-----

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)