

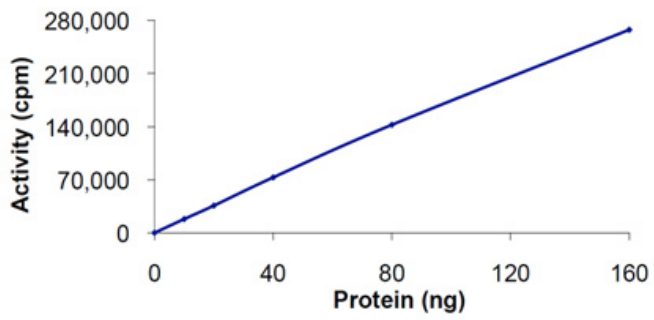
Akt1 (human), (recombinant) (GST- tag)

Citations: 1

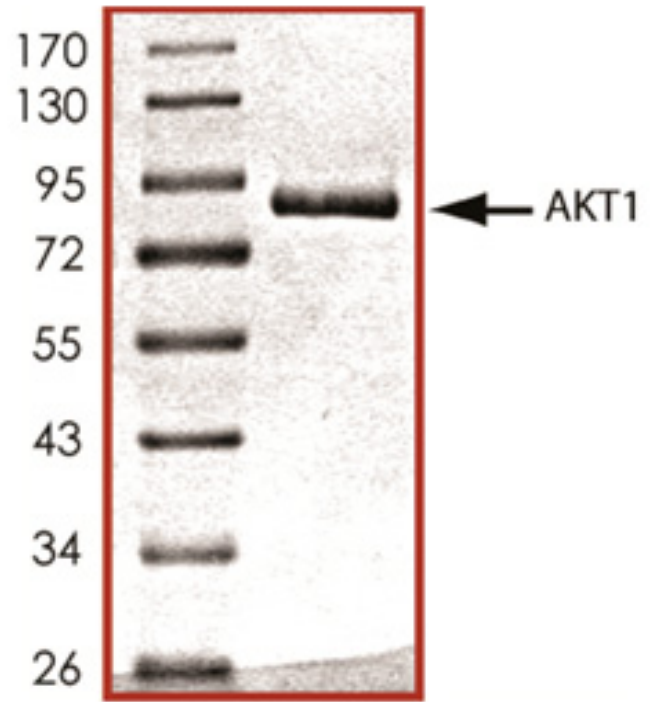
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Specific Activity: Sample Kinase Activity Plot. For specific information on a given lot, see related technical data sheet.



Purity: Sample Purity Data. For specific information on a given lot, see related technical data sheet.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Protein kinase B α , PKB α , RAC- α serine/threonine protein kinase

Application Notes Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.

MW 85 kDa

Source Produced in insect cells. Full length active enzyme.
Produced in a baculovirus expression system.

Technical Info / Product Notes Replacement for **ADI-PPK-410**

UniProt ID P31749

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)