

Akt2 (human), (recombinant) (GST- tag)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Application Notes Useful for kinetic, substrate, inhibitor studies and drug screening.

MW 68 kDa

Source Produced in a baculovirus expression system. The Δ PH domain of human Akt2 (PKB β , Protein Kinase B β) (aa 119-481) is fused at the N-terminus to a GST-tag.

UniProt ID P31751

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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