

Recombinant AKT1 protein

Catalog No: 31511

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.22 µg/µl

Source: Human

Buffer Contents: Full length recombinant AKT1 protein expressed in Sf9 cells at a concentration of 0.22 µg/µl in 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol.

Background: **AKT1** is a serine-threonine protein kinase that is activated by Platelet-Derived Growth Factor (PDGF) and PI 3-kinase signaling pathways. AKT has been shown to be a critical mediator of survival and apoptosis signaling events. Activation of AKT1 involves (and requires) phosphorylation of serine 473. AKT is also referred to as Protein Kinase B (PKB) and RAC.

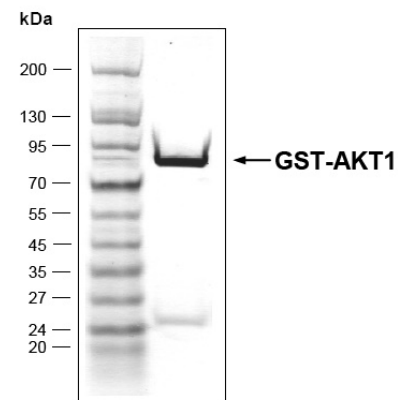
Protein Details: Recombinant AKT1 (accession number NP_005154.2) activated with PDK1, was expressed in Sf9 cells as an N-terminal GST-HIS6 fusion protein with a 3C cleavage site.

Application Notes: Recombinant AKT1 is suitable for kinase assays. The molecular weight of the fusion protein is ~ 81 kDa.

Specific activity: 55 pmol/µg×min ATP-KM: 5.0 µM

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant AKT1 protein gel. AKT1 protein was run on an SDS-PAGE gel and stained with Coomassie blue.