

ACKA Protein

Recombinant full-length bacterial protein expressed in E.coli cells

Catalog # A120-30G

Lot # V2491-5

Product Description

Full-length recombinant bacterial ACKA was expressed in E. coli cells using an N-terminal GST tag. The protein accession number is [NP_416799](#).

Gene Aliases

ECK2290; JW2293

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.25mM DTT, 0.1mM EGTA, 0.1mM EDTA, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

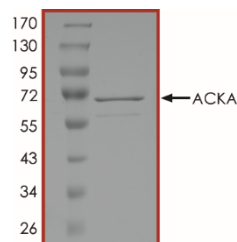
Scientific Background

ACKA belongs to the ASKHA superfamily of phosphotransferases, and is the prototypic carboxylate kinase. Widely distributed in the bacteria and archaea domains. Catalyzes the phosphorylation of acetate.

References

- Buss KA, et al: (2001) Urkinase: structure of acetate kinase, a member of the ASKHA superfamily of phosphotransferases. J Bacteriol. 183(2):680-686.
- <http://www.uniprot.org/uniprot/F3MTX6>

Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **71 kDa**.

ACKA Protein

Recombinant full-length bacterial protein expressed in E. coli

Catalog #	A120-30G
Lot #	V2491-5
Purity	>90%
Concentration	0.1µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.