

ACTR2 Protein

Full-length recombinant human protein expressed in sf9 cells

Catalog # A39-30G

Lot # U1763-7

Product Description

Full-length recombinant human ACTR2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC014546](#).

Gene Aliases

ARP2

Formulation

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

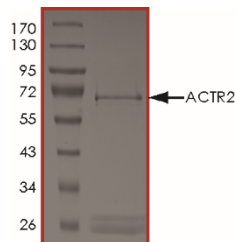
Scientific Background

Belongs to the actin-related protein (Arp) superfamily whose primary sequences exhibited similarity, but not identity, to conventional actin. Along with Arp3 and 5 other proteins, ACTR2 makes up the evolutionary conserved Arp2/3 complex that was first identified in *Acanthamoeba* using profiling affinity chromatography and is an essential and ubiquitous component of the actin filament network. Functions as ATP-binding component of the Arp2/3 complex which is involved in regulation of actin polymerization and together with an activating nucleation-promoting factor (NPF) mediates the formation of branched actin networks.

References

1. Machesky LM, et al: 1994. Purification of a cortical actin complex containing two unconventional actins from *Acanthamoeba* by affinity chromatography on profilin-agarose. *J. Cell Biol.* 127:107-15. 1994.
2. Schafer DA, et al: Actin-related proteins. *Annu Rev Cell Dev Biol.* 15:341-63. 1999.

Purity



The purity of ACTR2 protein was determined to be **>70%** by densitometry, ACTR2 approx. MW **70kDa**.

ACTR2 Protein

Full-length recombinant human protein expressed in sf9 cells

Catalog #	A39-30G
Lot #	U1763-7
Purity	>70%
Concentration	0.05 µg/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.