

RIOK1 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # R04-34G

Lot # A1159-3

Product Description

Full-length recombinant human RIOK1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC006104](#).

Gene Aliases

AD034; RRP10; bA288G3.1; FLJ30006; MGC12903

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

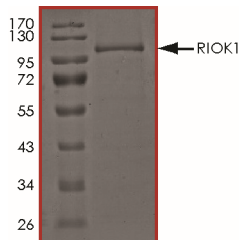
Scientific Background

RIOK1 or RIO kinase 1 is a novel cytoplasmic component of pre-40S pre-ribosomal particle(s) in human cells which is required in the final stages of cytoplasmic pre-40S maturation (1). RioK1 is a new interactor of protein arginine methyltransferase 5 (PRMT5) which competes with pICln for binding and modulates PRMT5 complex composition and substrate specificity (2). RIOK1 may reveal new insights into defects underlying glioblastoma and related cancers and may reveal new therapeutic opportunities for these cancers.

References

1. Widmann B. et.al: The kinase activity of human Rio1 is required for final steps of cytoplasmic maturation of 40S subunits. Mol Biol Cell. 2012 Jan;23(1):22-35. doi: 10.1091/mbc.E11-07-0639. Epub 2011 Nov 9.
2. Guderian G1. et.al: RioK1, a new interactor of protein arginine methyltransferase 5 (PRMT5), competes with pICln for binding and modulates PRMT5 complex composition and substrate specificity. J Biol Chem. 2011 Jan 21;286(3):1976-86. doi: 10.1074/jbc.M110.148486. Epub 2010 Nov 16.

Purity



The purity of RIOK1 protein was determined to be **>95%** by densitometry. Approx. MW **110 kDa**.

RIOK1 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # R04-34G

Lot # A1159-3

Purity >95%

Concentration 0.1 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.