

RIOK2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # R05-34G

Lot # Z1385-2

Product Description

Full-length recombinant human RIOK2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC000953](#).

Gene Aliases

FLJ11159

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.

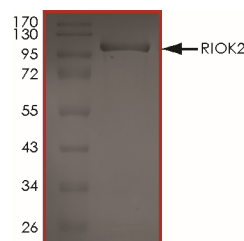
Scientific Background

RIOK2 or RIO kinase 2 is a novel cytoplasmic component of pre-40S pre-ribosomal particle(s) in human cells which is required late 40S maturation at several distinct steps (1). RIOK2 regulates metaphase-anaphase transition during mitotic progression (2).

References

1. Zemp I. et.al: Distinct cytoplasmic maturation steps of 40S ribosomal subunit precursors require hRio2. J Cell Biol. 2009 Jun 29;185(7):1167-80. doi: 10.1083/jcb.200904048.
2. Liu T. et.al: Phosphorylation of right open reading frame 2 (Rio2) protein kinase by polo-like kinase 1 regulates mitotic progression. J Biol Chem. 2011 Oct 21;286(42):36352-60. doi: 10.1074/jbc.M111.250175. Epub 2011 Aug 31.

Purity



The purity of RIOK2 protein was determined to be **>95%** by densitometry. Approx. MW **106 kDa**.

RIOK2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	R05-34G
Lot #	Z1385-2
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.