

KSR2, Unactive

Recombinant full-length protein expressed in Sf9 cells

Catalog # K08-15G

Lot # Y959-3

Product Description

Recombinant human KSR2 (554-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The KSR2 protein accession number is [Q6VAB6](#).

Gene Aliases

(None)

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.

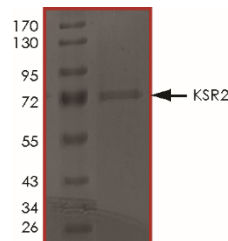
Scientific Background

KSR2 or kinase suppressor of Ras 2 specifically regulates the activity of MEKK3 and COT but not other members of the MAP3K family (1). KSR interacts with a regulatory Raf molecule in cis to induce a conformational switch of MEK, facilitating MEK's phosphorylation by a separate catalytic Raf molecule in trans (2).

References

- Brennan, D. F. et al: A Raf-induced allosteric transition of KSR stimulates phosphorylation of MEK. *Nature* 472: 366-369, 2011.
- Channavajhala, P. L. et al: hKSR-2 inhibits MEKK3-activated MAP kinase and NF-kappa-B pathways in inflammation. *Biochem. Biophys. Res. Commun.* 334: 1214-1218, 2005.

Purity



The purity of KSR2 was determined to be **>80%** by densitometry, approx. MW **72kDa**.

KSR2, Unactive

Recombinant human protein expressed in Sf9 cells

Catalog #	K08-15G
Lot #	Y959-3
Purity	>80%
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.