

CRKL Protein

Full length recombinant protein expressed in insect cells

Catalog # C53-30H

Lot # N212-2

Product Description

Recombinant full length human CRKL was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_005207](#).

Gene Aliases

crk-like protein

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

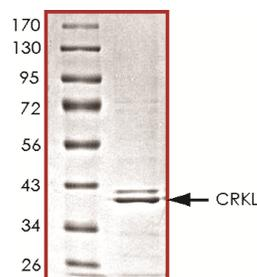
Scientific Background

CRKL or v-crkl sarcoma virus CT10 oncogene homolog (avian)-like is a protein kinase containing SH2 and SH3 (src homology) domains that activate the RAS and JUN kinase signaling pathways and transform fibroblasts in a RAS-dependent fashion (1). It is a substrate of the BCR-ABL tyrosine kinase which plays a role in fibroblast transformation by BCR-ABL, and may be oncogenic (2). CRKL is also required for normal cellular responses to Fgf8, including survival and migration, Erk activation, and target gene expression.

References

1. Ten Hoeve.J et.al: Isolation and chromosomal localization of CRKL, a human crk-like gene. *Oncogene* 8: 2469-2474, 1993.
2. Senechal, K. et.al: The CRKL adaptor protein transforms fibroblasts and functions in transformation by the BCR-ABL oncogene. *J. Biol. Chem.* 38: 23255-23261, 1996.

Purity



The purity was determined to be **>80%** by densitometry. Approx. MW **39 kDa**.

CRKL Protein

Full length recombinant protein expressed in insect cells

Catalog Number	C53-30H
Specific Lot Number	N212-2
Purity	>80%
Concentration	0.2µg/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.