

## VRK3 Protein

Recombinant full-length protein expressed in Sf9 cells

**Catalog # V03-30G**

Lot # F590-1

### Product Description

Recombinant full-length human VRK3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The VRK3 gene accession number is [NM\\_016440](#).

### 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

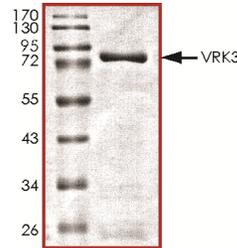
VRK3 is a member of the vaccinia-related kinase (VRK) family of serine/threonine protein kinases which has substitutions at several residues within the ATP binding motifs that in other kinases have been shown to be required for catalysis. Profound changes to the active site region underlie the loss of catalytic activity of VRK3 since it cannot bind ATP because of residue substitutions in the binding pocket (1). VRK3 is widely expressed in human tissues and the protein localizes to the nucleus. VRK3 regulates several transcription factors, nuclear envelope assembly, and chromatin condensation and is also required for cell cycle progression (2).

### References

1. Scheeff, E D. et al: Structure of the pseudokinase VRK3 reveals a degraded catalytic site, a highly conserved kinase fold, and a putative regulatory binding site. Structure. 2009 Jan 14;17(1):128-38.
2. Sanz-García M, et.al: Proteomics identification of nuclear Ran GTPase as an inhibitor of human VRK1 and VRK2 (vaccinia-related kinase) activities. Mol Cell Proteomics. 2008 Nov;7(11):2199-214.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

### Purity



The purity of VRK3 was determined to be **>95%** by densitometry, approx. MW **80kDa**.

## VRK3 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog Number	V03-30G
Specific Lot Number	F590-1
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.