

Recombinant VRK1 protein

Catalog No: 31243

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.5 µg/µl in 20 mM Tris-HCl, pH 7.3, 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Background: VRK1, Vaccinia-related Kinase 1 belongs to a family of serine/threonine protein kinases. It phosphorylates histone, casein and transcription factors ATF2 and c-Jun. VRK-1 also phosphorylates Thr18 of p53/TP53 to reduce p53 ubiquitination. VRK1 is highly expressed in actively dividing cells, such as fetal liver, testis and thymus.

Protein Details: Recombinant VRK1 protein was expressed in *E. coli* as the full-length human protein (accession number NP_003375) with an N-terminal FLAG tag. The molecular weight of the protein is ~48 kDa.

Application Notes: Recombinant VRK1 protein is suitable for use in kinase assays. We recommended to start with a range of protein concentrations (*e.g.* 5 ng, 50 ng, and 500 ng) to determine the optimal amount needed for each assay.

Storage and Guarantee: This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.