

## PEAK1, Unactive

Recombinant human protein expressed in Sf9 cells

**Catalog # P18-14G**

Lot # F576-3

### Product Description

Recombinant human PEAK1 (1127-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PTK9 gene accession number is [NM\\_024776](#).

### Gene Aliases

SGK269

### 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.

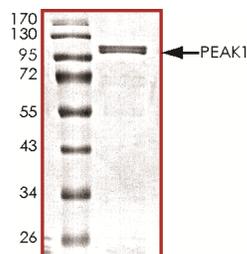
### Scientific Background

PEAK1 or Pseudopodium-enriched atypical kinase 1 is a protein tyrosine kinase which belongs to NKF3 kinase family and regulates the cytoskeleton and has been shown to be involved in cancer progression (1). PEAK1 promotes anchorage-independent growth in the MDA-MB-435 human cancer cell line and promotes tumor progression of XPA-1 human pancreatic cancer cells when transplanted into the pancreas of nude mice. PEAK1 is highly expressed in spleen, followed by heart, ovary, brain, lung, kidney, and fetal brain (2).

### References

1. Wang, Y. et.al: Pseudopodium-enriched atypical kinase 1 regulates the cytoskeleton and cancer progression. Proc. Nat. Acad. Sci. 107: 10920-10925, 2010.
2. Ohara, O. et.al: Characterization of size-fractionated cDNA libraries generated by the in vitro recombination-assisted method. DNA Res. 9: 47-57, 2002.

### Purity



The purity of PEAK1 was determined to be **>95%** by densitometry. Approx. MW **100 kDa**.

## PEAK1, Unactive

Recombinant human protein expressed in Sf9 cells

|                     |   |
|---------------------|---|
| Catalog Number      | P18-14G   |
| Specific Lot Number | F576-3  |
| 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**