

PINK1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P33-35G

Lot # P1736-4

Product Description

Recombinant human PINK1 (114-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number for PINK1 is [BC028215](#).

Gene Aliases

BRPK, PARK6, FLJ27236

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.

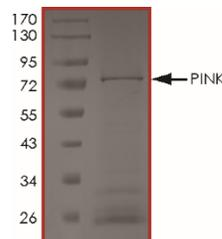
Scientific Background

PTEN-induced kinase 1 (PINK1) is a serine/threonine protein kinase that was identified initially in cancer cells as a gene up-regulated by overexpression of the major tumor suppressor, PTEN. It localizes to mitochondria, and regulates multiple aspects of mitochondrial biology. It is implicated in ubiquitin-dependent protein catabolic process to clear damaged mitochondria through selective autophagy (mitophagy). Mutations in the PINK1 gene are linked to early-onset recessive Parkinson disease.

References

1. Unoki M, Nakamura Y. Growth-suppressive effects of BPOZ and EGR2, two genes involved in the PTEN signaling pathway. *Oncogene*. 20:4457-652001, 2001.
2. Aerts L, et al. PINK1 kinase catalytic activity is regulated by phosphorylation on serines 228 and 402. *J Biol Chem*. 290:2798-811, 2015.

Purity



The purity of PINK1 was determined to be **>70%** by densitometry, PINK1 approx. MW **78kDa**.

PINK1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	P33-35G
Lot #	P1736-4
Purity	>70%
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.