

Parkin Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # P215-30G

Lot # J461-3

Product Description

Recombinant full-length human Parkin was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC022014](#).

Gene Aliases

PARK2, AR-JP, PDJ, PRKN, LPRS2

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

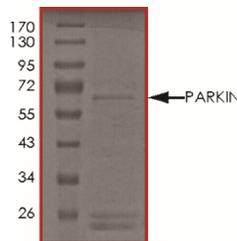
Scientific Background

Parkin or parkinson protein 2 is a RING domain-containing E3 ubiquitin protein ligase involved in proteasome-dependent degradation of proteins. Mutations in Parkin can lead to Parkinson's disease. Parkin is important for mitochondrial quality control by lysosome-dependent degradation of damaged mitochondria through autophagy, or mitophagy (1). Parkin plays an important role in directly supporting mitochondrial function and protecting mitochondrial genomic integrity from oxidative stress (2). Parkin also has essential role in regulating normal respiratory function in mitochondria and in the protection of cells from oxidative stress.

References

1. Yoshii, S. R. et.al: Parkin mediates proteasome-dependent protein degradation and rupture of the outer mitochondrial membrane. *J. Biol. Chem.* 286: 19630-19640, 2011.
2. Rothfuss, O. et.al: Parkin promotes mitochondrial genome integrity and supports mitochondrial DNA repair. *Hum. Molec. Genet.* 18: 3832-3850, 2009.

Purity



The purity of Parkin was determined to be **>70%** by densitometry. Approx. MW **68 kDa**.

Parkin Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	P215-30G
Lot #	J461-3
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.