

## PARK7 (DJ-1) Protein

Recombinant protein expressed in E.coli cells

**Catalog # P219-31H**

Lot # F495-3

### Product Description

Recombinant human PARK7 (DJ-1) (19-end) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM\\_007262](#).

### Gene Aliases

PARK7; DJ-1; DJ1

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . 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.

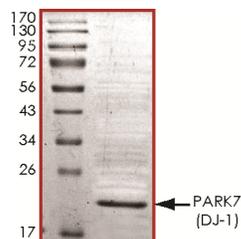
### Scientific Background

PARK7 or parkinson protein 7 belongs to the peptidase C56 family of proteins which acts as a positive regulator of androgen receptor-dependent transcription. PARK7 also functions as a redox-sensitive chaperone, as a sensor for oxidative stress, and it apparently protects neurons against oxidative stress and cell death. PARK7 mutations that impair transcriptional co-activator function can render dopaminergic neurons vulnerable to apoptosis and may contribute to the pathogenesis of Parkinson disease (1). PARK7 is an atypical peroxiredoxin-like peroxidase that scavenges hydrogen peroxide through oxidation of cys106 (2).

### References

- Xu, J.et.al: The Parkinson's disease-associated DJ-1 protein is a transcriptional co-activator that protects against neuronal apoptosis. *Hum. Molec. Genet.* 14: 1231-1241, 2005.
- Andres-Mateos.et.al: DJ-1 gene deletion reveals that DJ-1 is an atypical peroxiredoxin-like peroxidase. *Proc. Nat. Acad. Sci.* 104: 14807-14812, 2007.

### Purity



The purity of PARK7 (DJ-1) was determined to be **>85%** by densitometry. Approx. MW **22 kDa**.

## PARK7 (DJ-1) Protein

Recombinant protein expressed in E. coli cells

Catalog Number	P219-31H
Specific Lot Number	F495-3
Purity	>85%
Concentration	0.2µg/µl
Stability	1yr At $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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.**