

## DNAPK Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # D08-35G**

Lot # Q2357-9

### Product Description

Recombinant human DNAPK (3746-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM\\_006904](#).

### Gene Aliases

PRKDC, HYRC, p350, DNA-PK, DNP1, HYRC1, XRCC7

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, and 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.

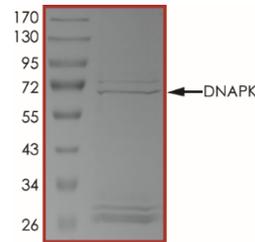
### Scientific Background

DNAPK is a member of the PI3/PI4-kinase family which encodes the catalytic subunit of a nuclear DNA-dependent serine/threonine protein kinase (DNA-PK), which is involved in DNA nonhomologous end-joining (NHEJ) during DNA double-strand break (DSB) repair and for V(D)J recombination during immune development. The second component of DNA-PK is Ku, which is required for proper activation of PRKDC (1). The initial events in retroviral integration are detected as DNA damage by the host cell, and that completion of the integration process requires the DNA-PK-mediated repair pathway (2).

### References

1. Van der Burg, M.et.al: A DNA-PKcs mutation in a radiosensitive T-B-SCID patient inhibits Artemis activation and nonhomologous end-joining. J. Clin. Invest. 119: 91-98, 2009.
2. Daniel, R.et.al: A role for DNA-PK in retroviral DNA integration. Science 284: 644-647, 1999.

### Purity



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

## DNAPK Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	D08-35G
Lot #	Q2357-9
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](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.**