

## ADCK3 Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # A37-11G**

Lot # Z1457-3

### Product Description

Recombinant human ADCK3 (156-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM\\_020247](#).

### Gene Aliases

ARCA2; CAB1; COQ10D4; COQ8; SCAR9

### 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.

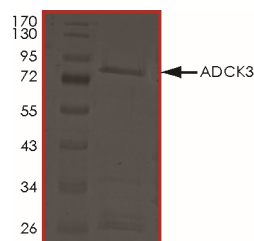
### Scientific Background

ADCK3 or aarF domain containing kinase 3 is a mitochondrial protein similar to yeast ABC1, which functions in an electron-transferring membrane protein complex in the respiratory chain. ADCK3 is involved in coenzyme Q10 synthesis which is essential for proper functioning of the mitochondrial respiratory chain (1). Expression of ADCK3 is induced by the tumor suppressor p53 and in response to DNA damage, and inhibiting its expression partially suppresses p53-induced apoptosis (2).

### References

1. Lagier-Tourenne, C. et.al: ADCK3, an ancestral kinase, is mutated in a form of recessive ataxia associated with coenzyme Q(10) deficiency. Am. J. Hum. Genet. 82: 661-672, 2008.
2. Iizumi, M. et.al: Isolation of a novel gene, CAB1, encoding a mitochondrial protein that is highly homologous to yeast activity of bc1 complex. Cancer Res. 62: 1246-1250, 2002.

### Purity



The purity of ADCK3 protein was determined to be **>75%** by densitometry. Approx. MW **83 kDa**.

## ADCK3 Protein

Recombinant human protein expressed in Sf9 cells

|                    |   |
|--------------------|---|
| Catalog #          | A37-11G   |
| Lot #              | Z1457-3   |
| Purity             | >75%  |
| 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.**