

## p63 Protein

Full-length recombinant protein expressed in Sf9 cells

**Catalog # P06-30G**

Lot # X561-3

### Product Description

Full-length recombinant human p63 was expressed by baculovirus in Sf9 cells using an N-terminal GST tag. The gene accession number is [NM\\_001114980](#).

### Gene Aliases

TP63; AIS; B(p51A); B(p51B); EEC3; KET; LMS; NBP; OFC8; p40; p51; p53CP; p73H; p73L; RHS; SHFM4; TP53CP; TP53L; TP73L

### 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  $-70^{\circ}\text{C}$ . 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.

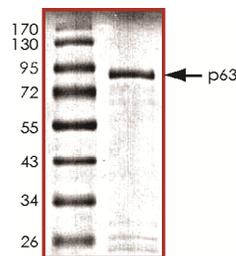
### Scientific Background

P63 is the primordial member of the p53 family that acts in a conserved process of monitoring the integrity of the female germline, whereas the functions of p53 are restricted to vertebrate somatic cells for tumor suppression. P63 protein plays an important role in the development and maintenance of stratified epithelial tissues (1). P63 is critical for maintaining the progenitor-cell populations that are necessary to sustain epithelial development and morphogenesis (2). Mutations in P63 are associated with ectodermal dysplasia, cleft lip/palate syndrome 3 (EEC3) and split-hand/foot malformation 4 (SHFM4).

### References

1. Yang, A., Schweitzer, R. et.al: p63 is essential for regenerative proliferation in limb, craniofacial and epithelial development. *Nature* 398: 714-718, 1999.
2. Suh, E.-K. et.al: p63 protects the female germ line during meiotic arrest. *Nature* 444: 624-628, 2006.

### Purity



The purity of p63 was determined to be **>95%** by densitometry. Approx. MW **92 kDa**.

## p63 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	P06-30G
Specific Lot Number	X561-3
Purity	>95%
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.**