

## p73 $\gamma$ Protein

Full-length recombinant protein expressed in Sf9 cells

**Catalog # P08-30CG**

Lot # E249-1

### Product Description

Full-length recombinant human p73 $\gamma$  was expressed by baculovirus in Sf9 cells using an N-terminal GST tag. The protein accession number is [O15350-3](#).

### Gene Aliases

TP73, TAp73gamma

### 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.

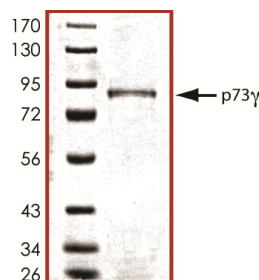
### Scientific Background

p73 $\gamma$  is a member of the p53 family of transcription factors which are involved in cellular responses to stress and development. There are four isoforms of p73 (p73 $\alpha$ , p73 $\beta$ , p73 $\delta$  and p73 $\gamma$ ) which differed in their abilities to form homodimeric and heterodimeric interactions with each other and with p53 (1). The p73 isoforms also differed in their abilities to activate transcription of the p21(Waf1) promoter and to inhibit colony formation. The p73 protein is expressed at very low levels in normal tissues and is differentially expressed in a number of tumors. p73 is a stress-response gene that activates transcription of p53-responsive genes and inhibits cell growth in a p53-like manner by inducing apoptosis (2).

### References

1. De Laurenzi, V. et al: Two new p73 splice variants, gamma and delta, with different transcriptional activity. J. Exp. Med. 188: 1763-1768, 1998.
2. Jost, C. A. et.al: p73 is a human p53-related protein that can induce apoptosis. Nature 389: 191-194, 1997.

### Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **85 kDa**.

## p73 $\gamma$ Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	P08-30CG
Specific Lot Number	E249-1
Purity	>90%
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.**