

## p73 $\beta$ Protein

Full-length recombinant protein expressed in Sf9 cells

**Catalog # P08-30BG**

Lot # E273-1

### Product Description

Full-length recombinant human p73 $\beta$  was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The protein accession number is [O15350-2](#).

### Gene Aliases

TP73, TAp73beta

### 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.

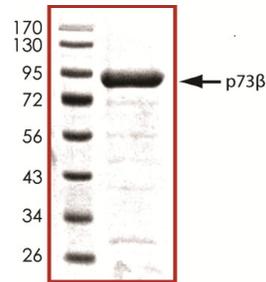
### Scientific Background

p73 $\beta$  is a member of the p53 family of transcription factors involved in cellular responses to stress and development. The p73 protein is expressed at very low levels in normal tissues and is differentially expressed in a number of tumors. P73 $\beta$  is strongly involved in malignancy acquisition and maintenance (1). 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). p73 is a component of a mismatch repair-dependent apoptosis pathway, which contributes to cisplatin-induced cytotoxicity. The regulation of p73 by c-Abl in response to DNA damage was also demonstrated by a failure of ionizing radiation-induced apoptosis after disruption of c-Abl-p73 interaction.

### References

1. Jost, C. A. et.al: p73 is a human p53-related protein that can induce apoptosis. Nature 389: 191-194, 1997.
2. Gong, J. et.al: The tyrosine kinase c-Abl regulates p73 in apoptotic response to cisplatin-induced DNA damage. Nature 399: 806-809, 1999.

### Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **88 kDa**.

## p73 $\beta$ Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	P08-30BG
Specific Lot Number	E273-1
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.**