

p73 α Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # P08-30G

Lot # W188-3

Product Description

Full-length recombinant human p73 α was expressed by baculovirus in Sf9 cells using an N-terminal GST tag. The gene accession number is [NM_005427](#).

Gene Aliases

TP73, TAp73alpha

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°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.

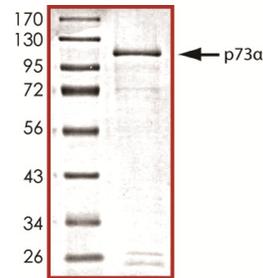
Scientific Background

p73 α is a member of the p53 family of transcription factors that are involved in cellular responses to stress and development. P73 has high sequence similarity to p53 and p63, which allows p63 and p73 to transactivate p53-responsive genes causing cell cycle arrest and apoptosis. P73 is a stress-response gene and a downstream effector in the p53 pathway (1). The p73 protein is expressed at very low levels in normal tissues and is differentially expressed in a number of tumors. p73 is the components of a mismatch repair-dependent apoptosis pathway which contributes to cisplatin-induced cytotoxicity (2).

References

1. Wang, J. et.al: p73 is a downstream target of p53 in controlling the cellular defense against stress. *J. Biol. Chem.* 282: 29152-29162, 2007.
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 **>90%** by densitometry. Approx. MW **110 kDa**.

p73 α Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	P08-30G
Specific Lot Number	W188-3
Purity	>90%
Concentration	0.2µg/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
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.