

## Product Information

<b>Description:</b>	Recombinant human p53 (cellular tumor antigen p53), encompassing amino acids 2-394(end). This construct contains an N-terminal FLAG-tag followed by a Precision cleavage site. This protein was affinity purified.
<b>Background:</b>	p53 is a transcription factor and tumor suppressor very frequently mutated in human cancer and often termed “guardian of the genome”. Activated by DNA damage, oxidative stress, or deregulated oncogene expression, p53 binds to a specific site in the promoter region of target genes and leads to the transcriptional activation of downstream genes involved in DNA repair, cell cycle arrest, senescence, and apoptosis. Inactivation of p53 promotes genome instability and directly contributes to cell transformation.
<b>Species:</b>	Human
<b>Construct:</b>	p53 (FLAG-2-394(end))
<b>Concentration:</b>	0.36 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, and 80 µg/ml FLAG peptide
<b>MW:</b>	46 kDa
<b>Genbank Accession:</b>	NM_000546
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Applications:</b>	Useful for SDS-PAGE.

## Quality Control Data

## 4-20% SDS-PAGE Coomassie Staining

