

Product Information

Description:	Recombinant human p53(tumor protein p53), full length, encompassing amino acids 2-393(end). In this construct, the N-terminal end of the protein is fused to GST via a short peptide linker placed between GST consisting of an optimal TEV protease target sequence that allows cleavage of GST (TEV: tobacco etch virus cysteine protease BPS Bioscience #50308). The construct also contains mutation of interest (Y220C). The recombinant protein was affinity purified
Species	Human
Construct:	p53 (Y220C) (GST-TEV-2-393(end))
Mutation:	Y220C
Concentration:	0.27 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	70 kDa
Genbank Accession:	NM_000546
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

