

Recombinant p53 protein

Catalog No: 31103

Expressed In: Baculovirus

Quantity: 5 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: 5 µg of Recombinant p53 protein in Dilution Buffer AM1 (20 mM Tris-Cl (pH 7.5), 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA).

Background: p53 is the most important tumor suppressor in the genome. It is responsive to numerous genotoxic stresses, which activates its transcription factor activity, in turn causing cell-cycle arrest by activating expression of p21 Cip/WAF. Mutant p53 that has lost its DNA-binding function interferes with the activity of native p53 and leads to oncogenic transformation. Alternatively, transformation may be caused by overexpression of Mdm2/Hdm2, an ubiquitin ligase specific for p53, which causes its destabilization. Inactivation of p53 is often coincident with hyperactivation of NFκB (NFκB p50 and NFκB p65), both of which serve to inhibit apoptosis.

Protein Details: Full-length recombinant p53 protein was expressed in a baculovirus system with a His-Tag and purified by an affinity column in combination with FPLC chromatography.

Active Motif also offers Recombinant p53 protein (Catalog No. 31465) that was expressed in *E. coli*.

Application Notes: Recombinant p53 is suitable for TransAM®, protein-protein and DNA protein assays. 1 ng is sufficient for a gelshift assay. The standard curve for TransAM p53 was generated using the range of 40 – 0.625 ng of protein.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

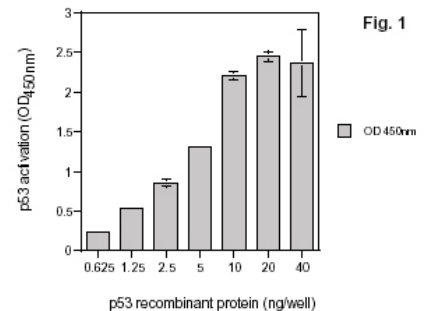
References:

This product was used in the following publications:

Miki, T., *et. al.* (2013). "p53 regulates Period2 expression and the circadian clock." *Nat. Commun.* 4:2444. PMID: 24051492.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



TransAM® standard curve generated using Recombinant p53 protein. The standard curve for TransAM® was generated using a range of 40 – 0.625 ng of protein and run on the TransAM® p53 ELISA Kit.