

p73 antibody (mAb)

Catalog No: 40972

RRID: AB_2793465

Clone: 5B429

Application(s): WB

Reactivity: Human

Quantity: 100 µg**Purification:** Protein G Chromatography**Host:** Mouse**Isotype:** IgG1, κ**Concentration:** 0.5 µg/µl**Molecular Weight:** 73 kDa

Background: p73 is a p53-related tumor protein, with some structural resemblance to p53, and is therefore considered a tumor-suppressor protein. p73 has numerous isoforms arising from alternative splicing, resulting in 6 different C-termini. Overexpression of p73 promotes growth arrest and/or apoptosis in a manner similar to that seen with p53.

Immunogen: This p73 antibody was raised against a full-length protein corresponding to human p73. This p73 antibody reacts with an epitope that is located on the N-terminal region of human p73 and does not cross react with p53. The p73 protein may run at 63 kDa in some cell lines due to post-transcriptional modifications.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

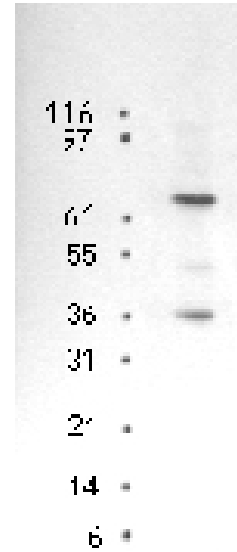
Applications Validated by Active Motif:

WB: 1:500 dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



p73 mAb tested by Western blot.
p73 monoclonal antibody used at a 1:500 dilution against 10 µg of transfected cell lysate.