

p73ΔN antibody (mAb)

Catalog No: 40973

RRID: AB_2793466

Clone: 38C674

Application(s): WB

Reactivity: Human

Quantity: 100 µg**Purification:** Affinity Purified**Host:** Mouse**Isotype:** IgG1**Concentration:** 0.5 µg/µl**Molecular Weight:** 73 kDa

Background: p73ΔN is a member of the p53 protein family which includes p53 as well as p63. Whereas p53 transactivates the retinoblastoma gene, p63 and p73 induce the cyclin-dependent kinase inhibitor p57 to maintain RB1 in an active, hypophosphorylated state. p73ΔN inhibits the functions of the p53 family in differentiation control, prevents myogenic differentiation, and enables cooperating oncogenes to transform myoblasts to tumorigenicity. p73ΔN is frequently overexpressed in rhabdomyosarcoma and is essential for tumor progression *in vivo*. These findings establish differentiation control as a key tumor suppressor activity of the p53 family.

Immunogen: This p73DN antibody was raised against a peptide corresponding to amino acid residues 2-13 of human p73ΔN. This antibody does not cross react with full-length p73.

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

Application Notes:

Applications Validated by Active Motif:

WB: 1 - 2 µg/ml 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.