

## p63 antibody (mAb)

---

**Catalog No:** 40974

**Clone:** 87C553

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human

**Purification:** Affinity Purified

**Host:** Mouse

**Concentration:** 0.5 µg/µl

**Quantity:** 100 µg

**Molecular Weight:** 63 kDa

---

**Background:** p63, also known as TP63 (for tumor protein p63), is a p53-related tumor protein with some structural resemblance to p53, and is therefore considered a tumor-suppressor protein. p63 has numerous isoforms arising from alternative splicing, resulting in 6 different C-termini. Overexpression of p63 promotes growth arrest and/or apoptosis in a manner similar to that seen with p53. It is associated with ectrodactyly-ectodermal dysplasia-cleft syndrome and Hay-Wells syndrome.

**Immunogen:** This p63 antibody was raised against a synthetic peptide corresponding to human p63.

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

**Application Notes:**

Validated Applications:

WB: 3 - 5 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all Blotting 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. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. 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.