

## p63 antibody (pAb)

**Catalog No:** 39739, 39740

**Isotype:** IgG

**Application(s):** ChIP, ICC, IF, WB

**Reactivity:** Human

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 0.3 µg/µl

**Volume:** 100 µl, 10 µl

**Molecular Weight:** 70 kDa

**Background:** p63 is a sequence-specific DNA binding transcription factor that is related to the p53 protein. p63 binds to and regulates a subset of genes also regulated by p53. p63 has been postulated to have a critical role in maintaining the identity of epithelial cell layers by regulating the differentiation of epithelial stem cells. Defects in p63 are associated with a number of developmental disorders involving the limbs and palate, as well as a variety of head and neck cancers.

**Immunogen:** This p63 antibody was raised against a recombinant protein corresponding to full-length human p63.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Validated Applications:

ChIP: 2 - 10 µl per ChIP

ICC/IF: 1:500 - 1:1,000 dilution

WB: 1:500 - 1:2,000 dilution

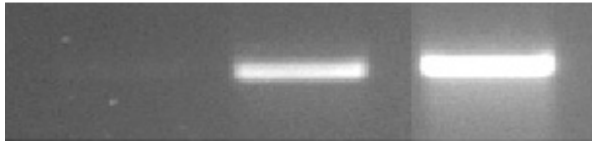
Published Applications:

ChIP:

**Storage and Guarantee:** Antibodies in solution can be stored at -20°C for 2 years. 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.

1 2 3



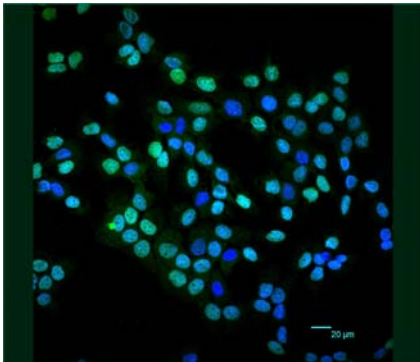
**p63 antibody (pAb) tested by ChIP.**

Chromatin IP was performed on chromatin from HaCaT cells. PCR primers were used to amplify a region of the FXDR gene promoter.

Lane 1: negative control IgG.

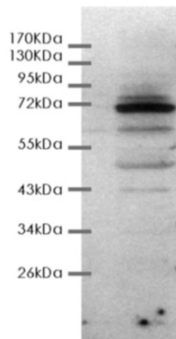
Lane 2: p63 antibody.

Lane 3: Input DNA control.



**p63 antibody (pAb) tested by immunofluorescence.**

HaCaT cells stained with p63 antibody (pAb) (1:500 dilution).



**p63 antibody (pAb) tested by Western blot.**

HaCaT cell nuclear extract (20 μg per lane) was probed with p63 antibody at a dilution of 1:2,000.