

KLF4 antibody (pAb)

Catalog Nos: 39745, 39746**RRID:** AB_2614947**Isotype:** IgG**Application(s):** WB**Reactivity:** Human, Wide Range Predicted**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 55 kDa

Background: KLF4 (Kruppel-like factor 4) is a transcription factor that binds GC-rich DNA sequences having a consensus core DNA binding sequence of CACCC. It is a member of the Kruppel-like factor (KLF) family, highly expressed in the gastrointestinal tract, skin and other epithelial tissues. Kruppel-like factors are a family of zinc-finger transcription factors that play important roles in many fundamental biologic processes including development, proliferation, differentiation, and apoptosis. KLF4 can function both as an activator and repressor of transcription, depending on the gene targeted. Generally, KLF4 serves as a cell cycle regulator, and it may be a tumor suppressor. KLF4 also appears to function as an oncogene in some situations. Additionally, overexpression of KLF4 (along with Oct3/4, Sox2 and c-Myc) can transform mouse fibroblasts into a state resembling embryonic stem cells (ES cells), referred to as Induced Pluripotency.

Immunogen: This KLF4 antibody was raised against recombinant full-length human KLF4.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

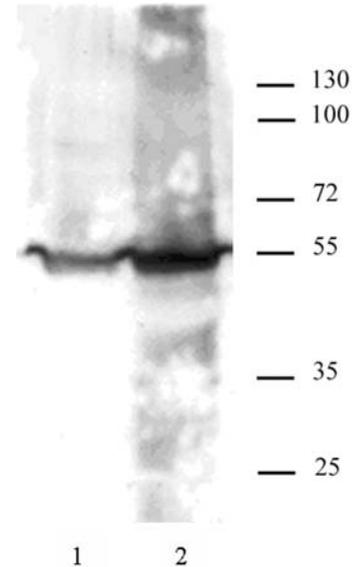
Application Notes:

Applications Validated by Active Motif:

WB: 1:2,000 - 1:5,000 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. 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.

**KLF4 antibody tested by Western blot.**

Lane 1: Nuclear extract (20 µg) of Hct116 cells.

Lane 2: Nuclear extract (20 µg) of primary human keratinocyte cells. Both probed with KLF4 antibody at a 1:5,000 dilution.