

KLF6 antibody (pAb)

Catalog Nos: 61243, 61244

RRID: AB_2793565

Isotype: IgG

Application(s): WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 40 kDa

Background: KLF6 (Kruppel-like factor 6) is member of the Kruppel-like family of transcription factors and is a zinc finger protein that acts as a transcriptional activator and functions as a tumor suppressor. Could play a role in B-cell growth and development. Multiple transcript variants encoding different isoforms have been found for this gene, some of which are implicated in carcinogenesis.

Immunogen: This antibody was raised against a peptide within the internal region of human KLF6.

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:

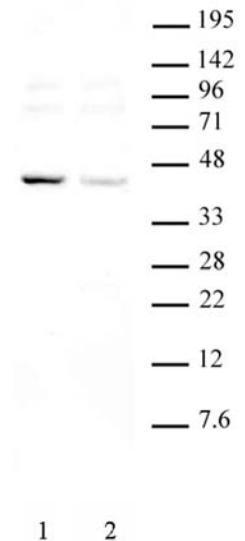
WB*: 0.5 - 2 µg/ml dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.



KLF6 pAb tested by Western blot.

3T3-L1 nuclear extract extract following stimulation with a defined hormonal mixture (DMSO, dexamethasone, insulin, and IBMX) at 30 µg per lane was probed with KLF6 pAb at a dilution of 2 µg/ml .

Lane 1: Day 0.

Lane 2: Day 10 - fully differentiated adipocytes.