

TIEG1 / KLF10 antibody (pAb)

Catalog Nos: 61117, 61118

RRID: AB_2615031

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 50 kDa

Background: TIEG1 / KLF10 (Transforming growth factor β -inducible early gene 1, Kruppel-like factor 10) is a transcriptional regulator involved in the control of cell growth. KLF10 binds to Sp1 / GC rich DNA sequences and can activate or repress the transcription of a number of genes. KLF10 is especially important for osteoblast function, including bone mineralization and support of osteoclast differentiation. KLF10 is rapidly induced by TGF β , bone morphogenetic protein, estrogen and epidermal growth factor. KLF10 plays a major role in TGF β -mediated inhibition of cell proliferation and inflammation and induction of apoptosis.

Immunogen: This TIEG1 / KLF10 antibody was raised against a peptide derived from human TIEG1 / KLF10.

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:

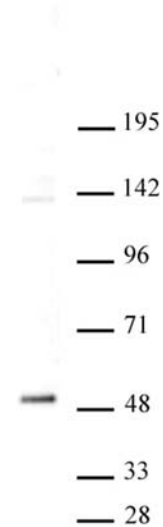
Applications Validated by Active Motif:

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

*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.



TIEG1 / KLF10 antibody (pAb) tested by Western blot

K-562 nuclear extract (20 μ g per lane) probed with the TIEG1 / KLF10 antibody (pAb) at a dilution of 1:500.