

EKLF antibody (mAb)

Catalog Nos: 61233, 61234

RRID: AB_2615069

Clone: 7B2

Isotype: IgG3

Application(s): ChIP, IP, WB

Reactivity: Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 42 kDa

Background: The **EKLF** (Erythroid krueppel like transcription factor, Erythroid transcription factor, KLF1) gene encodes a hematopoietic-specific transcription factor that induces high-level expression of adult beta-globin and other erythroid genes. The zinc-finger protein binds to the DNA sequence CCACACCCT found in the beta hemoglobin promoter. Furthermore, it binds to the BCL11A promoter and activates expression of BCL11A, which in turn represses the HBG1 and HBG2 genes.

Immunogen: This EKLF antibody was raised against a full-length recombinant protein to mouse EKLF (Kruppel-like factor 1 (erythroid)).

Buffer: Purified IgG in 12 mM Tris, pH 8.0, 60 mM Na Citrate, 120 mM NaCl with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.