

Kaiso antibody (mAb)

Catalog Nos: 39365, 39366

RRID: AB_2793229

Clone: 6F1D

Isotype: IgG

Application(s): WB

Reactivity: Mouse

Quantities: 200 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 95 kDa

Background: Kaiso (Clone 6F1D) is a zinc-finger transcription factor protein that binds to methylated and non-methylated GTGCNA motifs, and is also known as ZNF-Kaiso or ZBTB33. Kaiso interacts with the cell adhesion cofactor p120 catenin and may also recruit N-CoR repressor complex to promote histone deacetylation and repress transcriptional activity.

Immunogen: This Kaiso antibody was raised against a His-Tagged fusion protein corresponding to amino acids 103-499 of mouse Kaiso.

Buffer: PBS containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

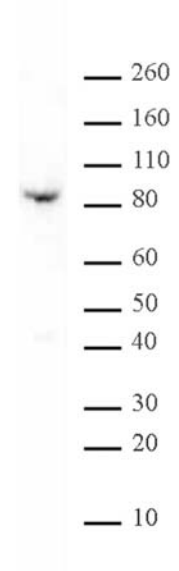
Application Notes:

Applications Validated by Active Motif:

WB: 0.5 - 4 µg/ml 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.



Kaiso mAb (Clone 6F1D) tested by Western blot.

Nuclear extract of CSML0 cells (20 µg) was probed with Kaiso mAb (Clone 6F1D) (1 µg/ml dilution).