

BRPF1 antibody (pAb)

Catalog Nos: 61541, 61542

RRID: AB_2614967

Isotype: IgG

Application(s): WB

Reactivity: Mouse

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.727 µg/µl

Molecular Weight: 150 kDa

Background: **BRPF1** (bromodomain and PHD finger containing, 1) is a component of the MOZ/MORF complex which has a histone H3 acetyltransferase activity. The bromodomain recognizes acetylated histones and the PWWP domain has been reported to interact with trimethylated Lys36 of Histone H3 (H3K36me3). Positively regulates the transcription of RUNX1 and RUNX2.

Immunogen: This antibody was raised against a peptide within the C-terminal region of mouse BRPF1.

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

Application Notes:

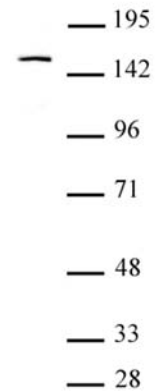
Applications Validated by Active Motif:

WB: 1:1,000 - 1:2,500 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.

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.



BRPF1 antibody (pAb) tested by Western blot. BRPF1 detection by Western blot. The analysis was performed using 20 µg mouse testis nuclear cell extract and BRPF1 antibody at a 1:1,000 dilution.