

PBRM1 antibody (pAb)

Catalog Nos: 61381, 61382

RRID: AB_2793612

Isotype: IgG

Application(s): ChIP-Seq, WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 180 kDa

Background: **PBRM1** (polybromo 1) is a component of the SWI/SNF-B (PBAF) ATP-dependent chromatin-remodeling complex. It is involved in transcriptional activation and repression of select genes by chromatin remodeling and is an integral component of complexes necessary for ligand-dependent transcriptional activation by nuclear hormone receptors. Defects in PBRM1 have been associated with primary clear cell renal cell carcinoma.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human PBRM1.

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:

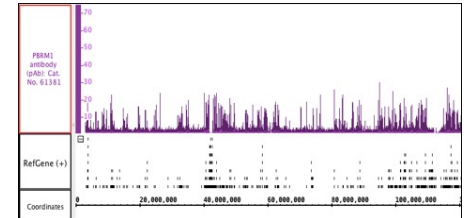
ChIP-Seq: 7 μ l (4 μ g) per ChIP

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

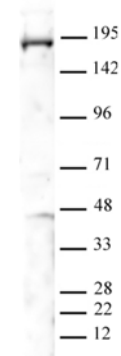
Addition of 0.05% Tween 20 in blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot.

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.



PBRM1 antibody (pAb) tested by ChIP-Seq. Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 μ g of chromatin from mouse liver and 4 μ g PBRM1 antibody. ChIP DNA was sequenced on the Illumina HiSeq and 3.9 million sequence tags were mapped to identify PBRM1 binding sites.



PBRM1 antibody (pAb) tested by Western blot.

Detection of PBRM1 by Western blot. The analysis was performed using 40 μ g of A-431 nuclear extract (Cat. No. 36004) and PBRM1 (pAb) at a 1:1000 dilution.