

SMAD1 phospho Ser214 (rAb), 100 µl

Catalog No: 91695

Clone: 3K8

Application(s): WB

Reactivity: Human, Mouse

Volume: 100 µl

Purification: Protein A Chromatography

Host: Rabbit

Isotype: IgG

Molecular Weight: 52 kDa

Background: SMAD family member 1 (**SMAD1**) have been involved in metastatic progression of many cancer types. **SMAD1** can be induced by many tumor-stimulating cytokines such as the bone morphogenetic protein 2 (BMP2) and TNF α and plays important roles in cell invasion and metastasis. BMP2 signaling is initiated by binding to its specific receptors, which leads to in the phosphorylation and nuclear translocation of Smad1. Translocated **SMAD1** then modulates the expression of downstream osteogenic genes.

Immunogen: This information is proprietary to Active Motif and/or its suppliers.

Buffer: Purified IgG in PBS, pH 7.3, with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

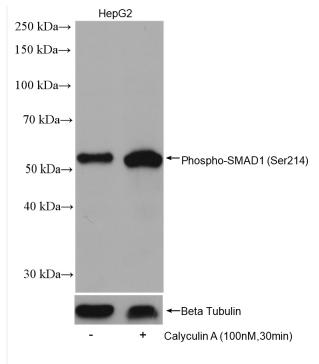
Application Notes:

Applications Validated by Active Motif:

WB: 1:2000-1:168000 dilution

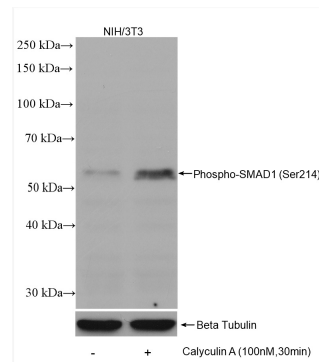
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



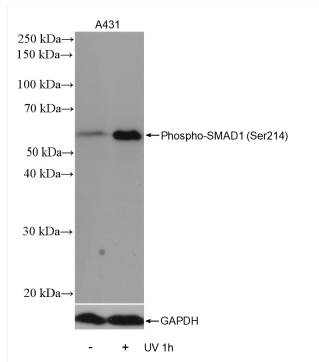
SMAD1 antibody (rAb) tested by Western blot.

Detection of SMAD1 by Western blot analysis using 30 µg of treated and non-treated A431 lysate and probed with SMAD1 antibody at 1:8000 dilution. The membrane was stripped and re-blotted with GAPDH antibody loading control.



SMAD1 antibody (rAb) tested by Western blot.

Detection of SMAD1 by Western blot analysis using 30 µg of treated and non-treated HepG2 lysate and probed with SMAD1 antibody at 1:8000 dilution. The membrane was stripped and re-blotted with GAPDH antibody loading control.



SMAD1 antibody (rAb) tested by Western blot.

Detection of SMAD1 by Western blot analysis using 30 µg of treated and non-treated NIH/3T3 lysate and probed with SMAD1 antibody at 1:8000 dilution. The membrane was stripped and re-blotted with GAPDH antibody loading control.