

C17orf96 phospho Ser15 antibody (pAb)

Catalog Nos: 61745, 61746

RRID: AB_2793754

Isotype: IgG

Application(s): DB

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 50 kDa

Background: C17orf96 has been shown to be highly expressed in mammalian ES cells and has been implicated as an interacting protein of the PRC2 complex (polycomb repressive complex 2). It may function as a CpG island binding protein and thus a potential regulator of gene transcription.

Phosphorylation of C17orf96 may be involved in regulation of the interaction of C17orf96 with PRC2.

Immunogen: This antibody was raised against a synthetic peptide containing phospho-serine 15 of human C17orf96.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

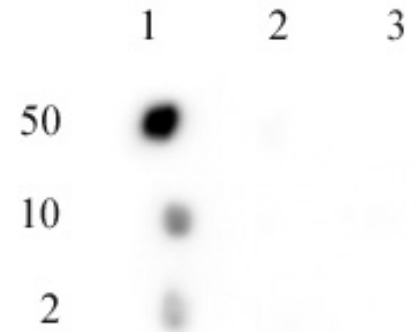
Application Notes:

Applications Validated by Active Motif:

DB: 1:1,000 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.



C17orf96 phospho Ser15 antibody (pAb) tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of C17orf96 phospho Ser15 antibody for phosphorylation of serine 15 of C17orf96. Peptides corresponding to the immunogen and related sequences derived were spotted onto PVDF and probed with the antibody at 1:1,000. The amount of peptide (picomoles) spotted is indicated next to each row.

Lane 1: Phospho serine 15 peptide.

Lane 2: Unmodified serine 15 peptide.