

## GLYR1 antibody (pAb)

**Catalog Nos:** 65693, 65694**RRID:** AB\_3216387**Application(s):** WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 65 kDa

**Background:** GLYR1 / NDF (nucleosome-destabilizing factor) that is recruited to gene bodies during transcriptional activation and facilitates Pol II transcription through nucleosomes, where it is proposed to “unzip” chromatin during transcription, pushing the polymerase past pause points. It was discovered as a factor that allows p300 to acetylate Histone H3K56, which is normally an inaccessible residue for this modification. GLYR1 is shown to associate with a subset of highly transcribed genes tending to have longer mRNAs rather than shorter, and is also at a significant subset of loci containing the Histone H3K36me3. GLYR1 is expressed at high levels in most tissue types, including stem cells, and has been found to be overexpressed in breast cancer.

**Immunogen:** This antibody was raised against a peptide within the C-terminal region of human GLYR1.

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

**Application Notes:**

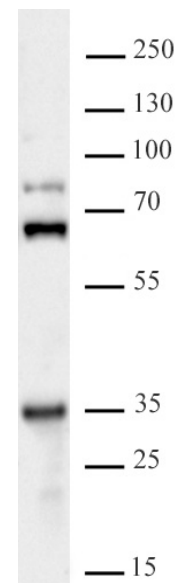
Applications Validated by Active Motif:

WB\*: 0.5 - 2 µg/ml

\*Note: Many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for 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.

**GLYR1 antibody (pAb) tested by Western blot**

20 µg of HEK293 nuclear extract was run on SDS-PAGE and probed with antibody at a dilution of 1:500.