

CoREST antibody (pAb)

Catalog Nos: 39955, 39956**RRID:** AB_2793407**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 66 kDa

Background: CoREST (REST corepressor 1, RCOR1) is a corepressor protein recruited by the REST, a neuronal gene-specific transcriptional repressor, to help repress neuron specific gene expression in non-neuronal cell lineages. CoREST is an essential component of the BHC repressor complex that also contains HDAC1, HDAC2, HMG20B/BRAF35, LSD1 (a lysine-4 specific histone demethylase) and PHF21A/BHC80. The Herpes Simplex virus ICP0 protein blocks the repression of viral genes by disrupting the interaction between CoREST and HDAC1/2.

Immunogen: This CoREST antibody was raised against a synthetic peptide derived from the C-terminal region of human CoREST.

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:

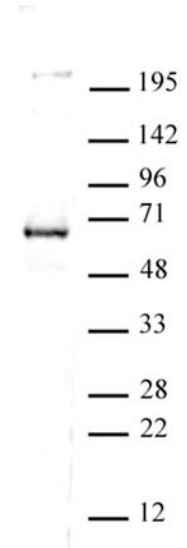
Applications Validated by Active Motif:

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

*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.

**CoREST antibody (pAb) tested by Western blot.**

Nuclear extract of HEK293 cells (25 µg) probed with CoREST antibody (1:500 dilution).