
RNA Polymerase II Polyclonal Antibody

(Catalog A-2033)

Background

This gene encodes the largest subunit of RNA polymerase II, the polymerase responsible for synthesizing messenger RNA in eukaryotes. The product of this gene contains a carboxy-terminal domain composed of heptapeptide repeats that are essential for polymerase activity. These repeats contain serine and threonine residues that are phosphorylated in actively transcribing RNA polymerase. In addition, this subunit, in combination with several other polymerase subunits, forms the DNA binding domain of the polymerase, a groove in which the DNA template is transcribed into RNA.

Description

RNA Polymerase II Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thiomersal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P24928

Purification

Affinity purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1135-1311 of human POLR2A (NP_000928.1).

Storage

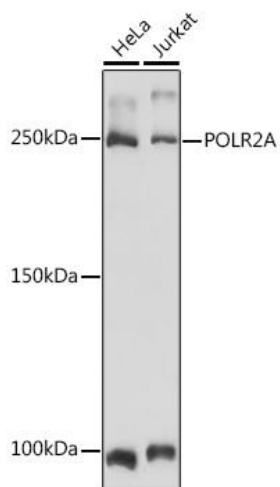
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

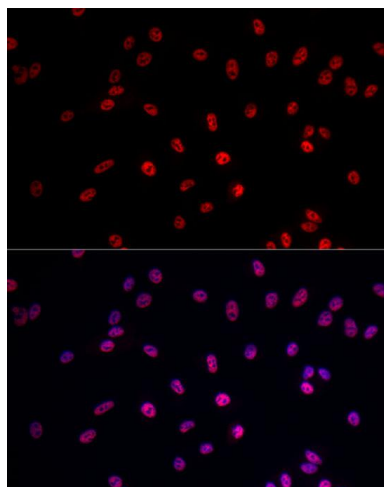
POLR2A; POLR2; POLRA; RPB1; RPBh1; RPO2; RPOL2; RpIILS; hRPB220; hsRPB1; Pol II, RNA Polymerase II

Application

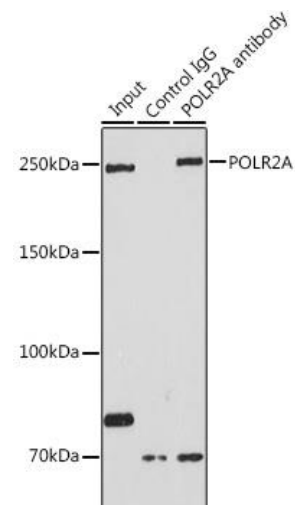
WB, IF, ICC, IP, ChIP; Recommended dilution: WB 1:500 - 1:2000, IF/ICC 1:50 - 1:200, IP 1:500 - 1:1000



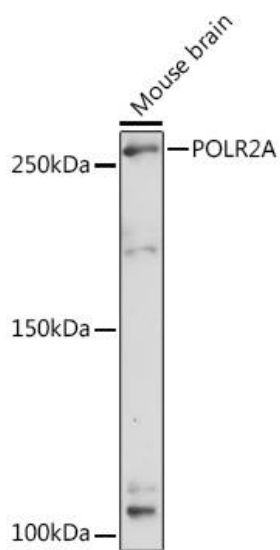
Western blot analysis of extracts of various cell lines, using RNA Polymerase II pAb at 1:3000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



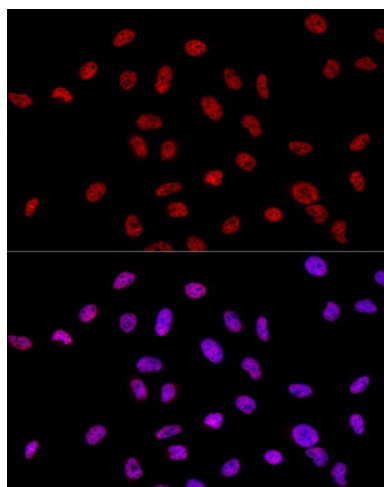
Immunofluorescence analysis of A-549 cells using RNA Polymerase II pAb at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Immunoprecipitation analysis of 200ug extracts of HeLa cells using 3ug RNA Polymerase II pAb. Western blot was performed from the immunoprecipitate using RNA Polymerase II pAb at a dilution of 1:1000.



Western blot analysis of extracts of Mouse brain, using RNA Polymerase II pAb at 1:3000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST.



Immunofluorescence analysis of HeLa cells using RNA Polymerase II pAb at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.