

AbFlex® Nucleolin antibody (rAb)

Catalog Nos: 91231, 91232

RRID: AB_2793813

Isotype: IgG2a

Application(s): WB

Reactivity: Canine, Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 100 kDa

Background: AbFlex® antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports. AbFlex® Nucleolin antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

Nucleolin (NCL, C23) is an abundant and ubiquitously expressed protein that is a major component of the nucleolus. Nucleolin seems to be overexpressed in highly proliferative cells and is involved in many aspects of gene expression: chromatin remodeling, DNA recombination and replication, RNA transcription by RNA polymerase I and II (RNA pol II), rRNA processing, mRNA stabilization, cytokinesis and apoptosis. Its role as a histone chaperone has recently been identified. It is involved in the synthesis and maturation of ribosomes, and is located mainly in dense fibrillar regions of the nucleolus.

Immunogen: This antibody was raised against a salt-extracted nuclear matrix preparation from MDCK (canine kidney) cells.

Buffer: Purified IgG in 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, 30% glycerol. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

WB*: 0.1 - 1 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.

AbFlex® recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 39541.

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

The file cannot be found:
91231-00418001-1.jpg

AbFlex® Nucleolin antibody (rAb) tested by Western Blot. 25 µg of HeLa cell nuclear extract (Lane 1) or 30 µg K562 cell nuclear extract (Lane 2) was run on SDS-PAGE and probed with AbFlex Nucleolin antibody at 0.5 µg/ml.