

## RBPJ antibody (mAb)

**Catalog Nos:** 61073, 61074**RRID:** AB\_2793496**Clone:** 7A11**Isotype:** IgG2b**Application(s):** WB**Reactivity:** Human**Quantities:** 100 µg, 10 µg**Purification:** Protein G Chromatography**Host:** Rat**Concentration:** 1 µg/µl**Molecular Weight:** 55 kDa

**Background:** RBPJ (Recombination Signal Binding Protein For Immunoglobulin Kappa J Region) is a transcriptional regulator that plays a central role in Notch signaling, a signaling pathway involved in cell-cell communication that regulates a broad spectrum of cell-fate determinations. Acts as a transcriptional repressor when it is not associated with Notch proteins. When associated with some NICD product of Notch proteins (Notch intracellular domain), it acts as a transcriptional activator that activates transcription of Notch target genes. Probably represses or activates transcription via the recruitment of chromatin remodeling complexes containing histone deacetylase or histone acetylase proteins, respectively. Specifically binds to the immunoglobulin kappa-type J segment recombination signal sequence. Binds specifically to methylated DNA.

**Immunogen:** This antibody was raised against full-length recombinant human RBPJ protein.

**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.

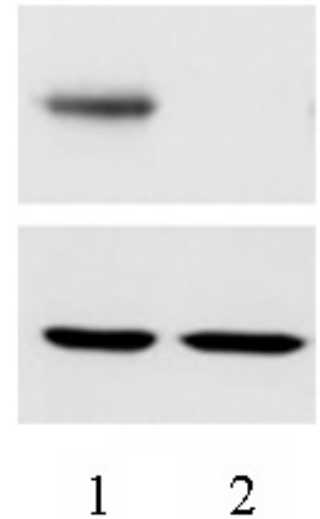
**Application Notes:**

Applications Validated by Active Motif:

WB: 2 - 4 µg/ml 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.

**Western blot of RBPJ antibody.**

Wild type (lane 1) or RBPJ KO (lane 2) DG75 cell lysate was probed with RBPJ antibody (top row) or a GAPDH antibody (bottom row) to show equal loading.