

RBPJ antibody (mAb)

Catalog Nos: 61505, 61506

RRID: AB_2614972

Clone: 1F1

Isotype: IgG2b

Application(s): ChIP, EMSA, ICC, IF

Reactivity: Human, Mouse

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Rat

Concentration: 1 µg/µl

Molecular Weight: 68 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 a full-length recombinant protein corresponding to human RBPJ.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

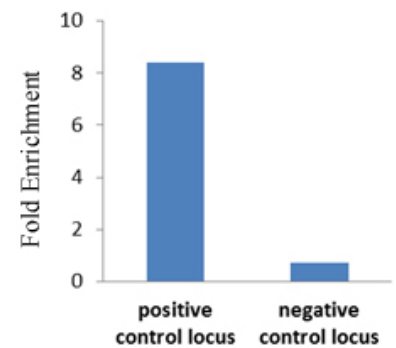
Application Notes:

Applications Validated by Active Motif:

ICC/IF: 0.5 µg/ml dilution

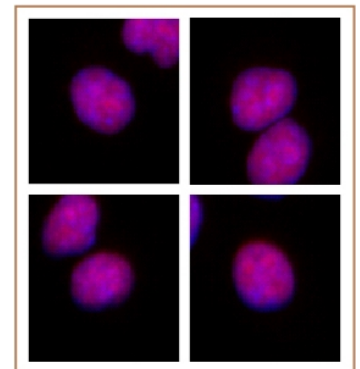
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



RBPJ antibody (mAb) tested by ChIP.

Chromatin IP was performed using chromatin of an Epstein-Barr virus infected lymphoblastoid cell line (2 x 10⁶ cell equivalents per ChIP) and RBPJ antibody or the equivalent amount of IgG negative control. Real time, quantitative PCR (RT-qPCR) was performed on DNA purified from each of the ChIP reactions using primer pairs for a positive and negative control region. Data are presented as Fold Enrichment of the ChIP



RBPJ antibody (mAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with RBPJ antibody at a 0.5 µg/ml dilution.