

AID antibody (mAb)

Catalog Nos: 39885, 39886**RRID:** AB_2793380**Clone:** 328.8b**Isotype:** IgG2a**Application(s):** ChIP, WB**Reactivity:** Human, Mouse**Quantities:** 100 µg, 10 µg**Purification:** Protein G Chromatography**Host:** Rat**Concentration:** 1 µg/µl**Molecular Weight:** 24 kDa

Background: AID (Activation-induced cytidine deaminase, AICDA) is a member of the cytidine and deoxycytidylate deaminase family. It acts as a single-strand DNA deaminase involved in somatic hypermutation, gene conversion and class-switch recombination in developing B-cells and is necessary for efficient antibody responses. It has also been implicated in active DNA demethylation, as deamination of methylcytosine yields thymine, which is read as a mismatch and is replaced by DNA damage repair enzymes.

Immunogen: This AID antibody was raised against a peptide corresponding to amino acids 178-198 of human AID.

Buffer: Purified IgG in 50 mM Tris-HCl (pH 7.2), 30% glycerol, and 0.035% sodium azide. Sodium azide is highly toxic.

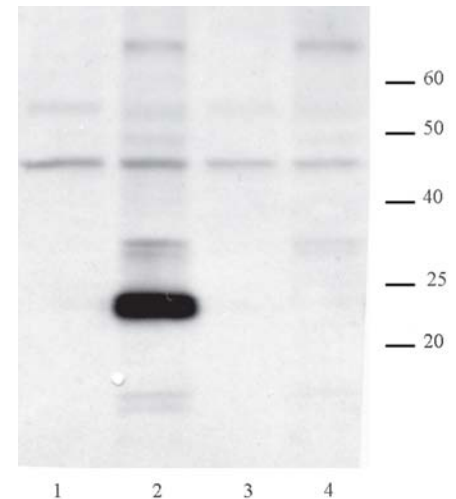
Application Notes:

Applications Validated by Active Motif:

WB: 1 - 3 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.

**AID antibody (mAb) tested by Western blot.**

Murine spleen cells were probed with AID antibody at a dilution of 3 µg/ml.

- Lane 1: Unstimulated spleen cells from wild type mice.
- Lane 2: Treated (LPS & IL4 for 96 hrs) spleen cells from wild type mice.
- Lane 3: Unstimulated spleen cells from AID knock-out mice.
- Lane 4: Treated (LPS & IL4 for 96 hrs) spleen cells from AID knock-out mice.