

## BCL11A antibody (mAb)

**Catalog Nos:** 61231, 61232

**RRID:** AB\_2793560

**Clone:** BCL11A/123

**Isotype:** IgG1

**Application(s):** ICC, IF, IHC, WB

**Reactivity:** Human

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 100 kDa

**Background:** **BCL11A** (B-cell CLL/lymphoma 11A) is a transcriptional regulator, a myeloid and B-cell proto-oncogene and plays important roles in leukemogenesis and hematopoiesis. BCL11A is involved in regulating hemoglobin gene expression. BCL11A regulates hemoglobin gene switching during fetal development and silencing of  $\gamma$ -globin expression in adults. BCL11A interacts with the Mi-2/NuRD chromatin remodeling complexes, as well as the erythroid transcription factors GATA1 and FOG1. Abundant expression of full-length forms of BCL11A is developmentally restricted to adult erythroid cells.

**Immunogen:** This BCL11A antibody was raised against a recombinant protein corresponding to amino acids 637-835 of BCL11A-XL (B-cell lymphoma/leukemia 11A extra long form).

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

**Application Notes:**

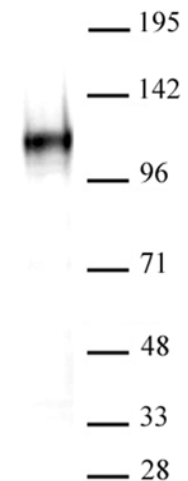
Applications Validated by Active Motif:

WB\*: 1 – 2 µg/ml dilution

\*Note: we recommend the addition of 0.05% Tween 20 to all blocking solutions to reduce background.

**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.



**BCL11A antibody (mAb) tested by Western blot.**

HeLa nuclear extract (35 µg per lane) probed with BCL11A antibody at 2 µg/ml dilution.