

Bcl6 antibody (mAb)

Catalog Nos: 61193, 61194

RRID: AB_2615054

Clone: IG191E

Isotype: IgG1

Application(s): IHC, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 80 kDa

Background: Bcl6 (B-cell lymphoma 6 protein, Bcl5, LAZ3, ZBTB27, ZNF51) is sequence-specific transcriptional repressor that is required for the formation of lymph node germinal centers (GC). Bcl6 is a member of the BTB-POZ family of transcriptional regulators and contains six Kruppel-type zinc fingers. The POZ domain in the amino terminus of Bcl6 is involved in the recruitment of corepressor proteins including NCoR, BCoR, SMRT, Mi-2/NURD and several HDACs. Bcl6 promotes proliferative expansion of B-cells in the germinal center through repression of genes regulating cell cycle and differentiation. Bcl6 also suppresses the DNA damage response, to prevent class-switch recombination being repaired. Unregulated expression of Bcl6 is involved in the development of many B-cell lymphomas.

Immunogen: This Bcl6 antibody was raised against a full-length protein corresponding to human Bcl6.

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

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91287.

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.