

## BCL6 polyclonal antibody - Classic

**Cat. No.** C15410221

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 40163

**Size:** 25 µl/100 µl

**Concentration:** 1 µg/µl

**Specificity:** Human, mouse.

**Purity:** Purified by antigen-affinity chromatography.

**Storage:** Store at -20°C; for long storage, store at -80°C.  
Avoid multiple freeze-thaw cycles.

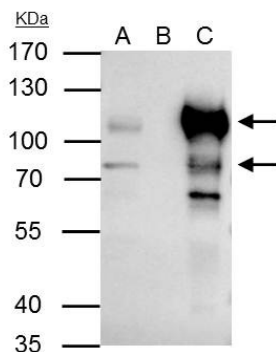
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

### Applications

|                     | Suggested dilution* | Results     |
|---------------------|---------------------|-------------|
| ChIP                | 5 µg/ChIP           | Figure 5    |
| IHC                 | 1:100-1:1,000       | Figure 4    |
| Immunoprecipitation | 1:100-1:500         | Figure 1    |
| Western blot        | 1:500-1:3,000       | Figure 2, 3 |

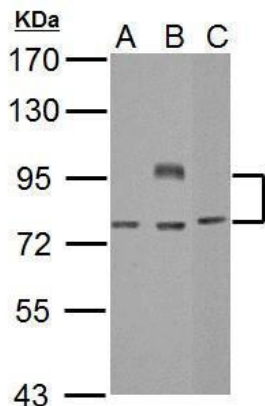
\*Optimal dilutions/concentrations should be determined by the researcher.

### Results



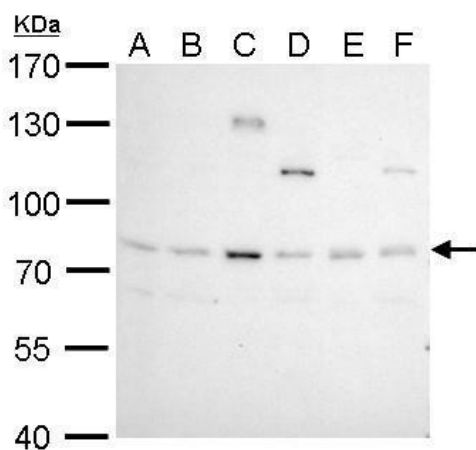
**Figure 1. Immunoprecipitation**

BCL6 antibody immunoprecipitates BCL6 protein in IP experiments. IP Sample: Raji whole cell extract A: 30 µg whole cell extract of BCL6 protein expressing Raji cells B: Control with 3 µg of pre-immune rabbit IgG C : Immunoprecipitation of BCL6 by 3 µg of BCL6 antibody. The immunoprecipitated BCL6 protein was detected by Western blot with the BCL6 antibody diluted 1:1,000.



**Figure 2. Western blot results**

Sample: 30 µg of whole cell lysate  
A: Jurkat  
B: Raji  
C: NCI-H929  
7.5% SDS PAGE  
BCL6 antibody diluted 1:1,000



**Figure 3. Western blot results**

BCL6 antibody detects BCL6 protein by Western blot analysis.

A. 30 µg Neuro2A whole cell extract

B. 30 µg GL261 whole cell extract

C. 30 µg C8D30 whole cell extract

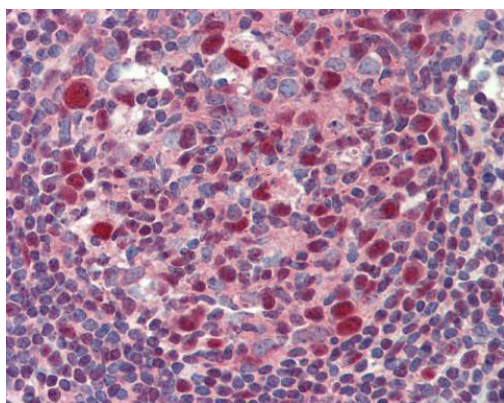
D. 30 µg NIH-3T3 whole cell extract

E. 30 µg Raw264.7 whole cell extract

F. 30 µg C2C12 whole cell extract

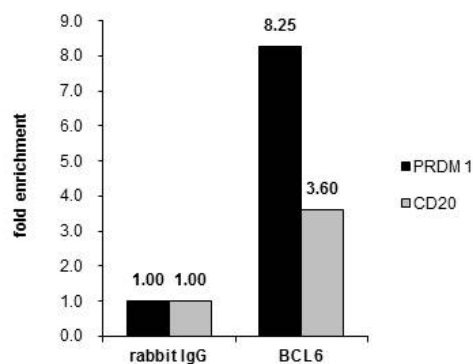
7.5 % SDS-PAGE

The BCL6 antibody (Cat. No. C15410221) was used at a dilution of 1:1,000



**Figure 4. IHC**

Immunohistochemical analysis of paraffin-embedded Tonsil, using BCL6 (C15410221) antibody (10 µg/ml).



**Figure 5. ChIP**

ChIP was performed with Raji chromatin extract and 5 µg of either control rabbit IgG or BCL6 antibody. The precipitated DNA was detected by PCR with primer set targeting to PRDM1 or CD20. Results are expressed as fold enrichment over the IgG negative control.