

PRODUCT NAME BCL7C polyclonal antibody		
Cat. No. C15310124 (CS-124-100)	Type: Polyclonal	Size: 100 µl
Lot #: A384-004	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against human BCL7C (B-cell CLL/lymphoma 7C), using two KLH-conjugated synthetic peptides containing a sequence from the central and from the C-terminal part of the protein, respectively.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:1,000	Fig 1
Western blotting	1:1,000	Fig 2

Purity: Whole antiserum from rabbit containing 0.05% azide.

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.

Last data sheet update: March 19, 2010

Target description

BCL7C (UniProt/Swiss-Prot entry Q8WUZ0) is homologous to the N-terminal region of BCL7A which was shown to be directly involved in a three-way gene translocation in a Burkitt lymphoma cell line. The function of BCL7A has not yet been elucidated, but it may play an anti-apoptotic role.

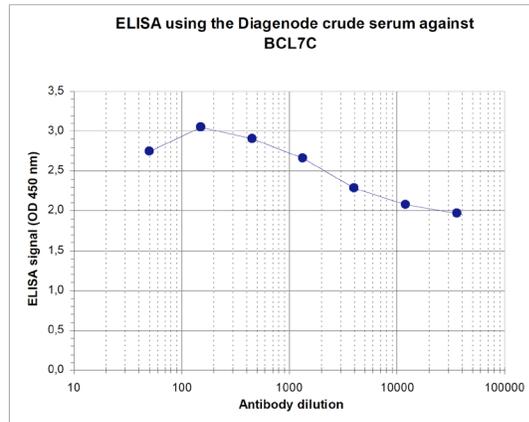


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human BCL7C (Cat. No. CS-124-100). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:209,000.

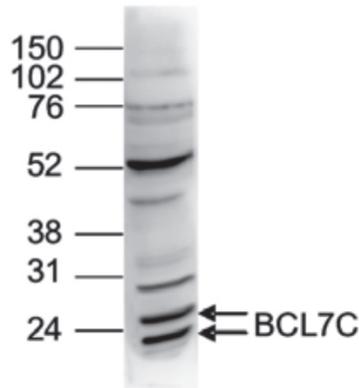


Figure 2
Western blot analysis using the Diagenode antibody directed against BCL7C

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against BCL7C (Cat. No. CS-124-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left. The antibody detects both known isoforms of the protein (expected sizes: 23.4 and 26.4 kDa).