

MycK323ac polyclonal antibody

Cat. No. C15410346

Type: Polyclonal	Specificity: Human: positive. Other species: not tested.
Size: 50 µg/63 µl	Isotype: NA
Concentration: 0.8 µg/µl	Source: Rabbit
Lot No.: A506-0041	Purity: Affinity purified polyclonal antibody.
Storage buffer: PBS containing 0.05% azide and 0.05% ProClin 300.	Storage conditions: 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: May 3, 2018

Description

Other names: v-Myc, c-Myc, MRTL, BHLHE39

Polyclonal antibody raised in rabbit against Myc (v-myc myelocytomatosis viral oncogene homolog) acetylated at lysine 323 (MycK323ac) using a KLH-conjugated synthetic peptide.

Applications

Applications	Suggested dilution	References
ELISA	1:100 - 1:500	Fig 1
IF	1:500	Fig 2

Target Description

Myc (UniProtKB/Swiss-Prot entry P01106) participates in the regulation of gene transcription. Myc seems to activate the transcription of growth-related genes. It specifically recognizes the core sequence 5'-CAC[GA]TG-3' but also binds DNA in a non-specific manner. Overexpression of MYC is implicated in the etiology of a variety of hematopoietic tumors.

Validation data

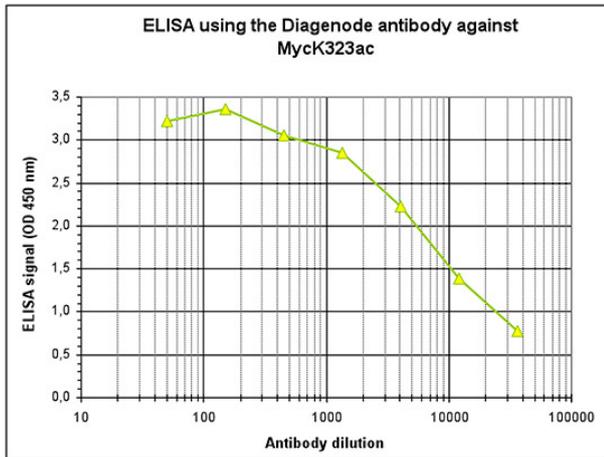


Figure 1. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human MycK323ac (Cat. No. C15410346). The plates were coated with the peptide 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 1:8,000.

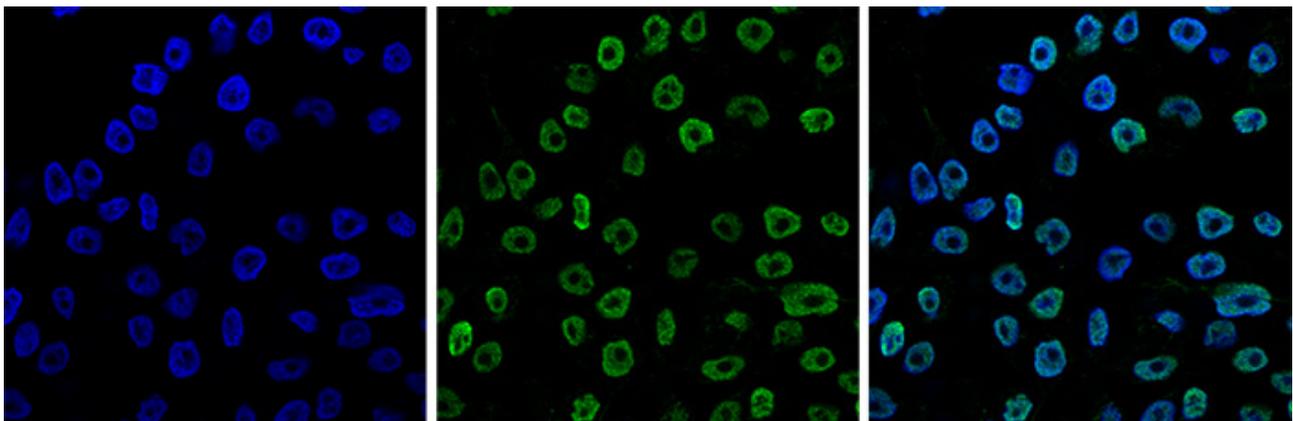


Figure 2. Immunofluorescence using the Diagenode antibody directed against MycK323ac

HeLa cells were stained with the Diagenode antibody against MycK323ac (cat. No. C15410346) and with DAPI. Cells were fixed with 4% formaldehyde for 10' and blocked with PBS/TX-100 containing 1% BSA. The cells were immunofluorescently labeled with the MycK323ac antibody (middle) diluted 1:500 in blocking solution followed by an anti-rabbit antibody conjugated to Alexa488. The left panel shows staining of the nuclei with DAPI. A merge of the two stainings is shown on the right.