

acrIIA4 Antibody

Cat. No. C15200248

Type: Monoclonal. IgG1 kappa.	Specificity: Listeria monocytogenes
Size: 50 µg/50 µl	Isotype: NA
Concentration: 1 µg/µl	Host: Mouse
Lot No.: 002	Purity: Protein A purified monoclonal antibody
Storage buffer: PBS containing 0.05 % Na-azide.	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: October 27, 2020

Description

Monoclonal antibody raised in mouse against acrIIA4 (anti CRISPR protein IIA4) using a recombinant protein.

Applications

Applications	Suggested dilution/amount	References
Western blotting	1:1,000	Fig 1

Target Description

CRISPR systems are adaptable immune mechanisms which are present in many bacteria to protect themselves from foreign nucleic acids, such as viruses, transposable elements or plasmids. Recently, the CRISPR/Cas9 (CRISPR-associated protein 9 nuclease) system from *S. pyogenes* has been adapted for inducing sequence-specific double stranded breaks and targeted genome editing. This system is unique and flexible due to its dependence on RNA as the moiety that targets the nuclease to a desired DNA sequence and can be used to induce indel mutations, specific sequence replacements or insertions and large deletions or genomic rearrangements at any desired location in the genome. In addition, Cas9 can also be used to mediate upregulation of specific endogenous genes or to alter histone modifications or DNA methylation. acrIIA4 is a protein that inhibits CRISPR/Cas9 activity.

Validation data

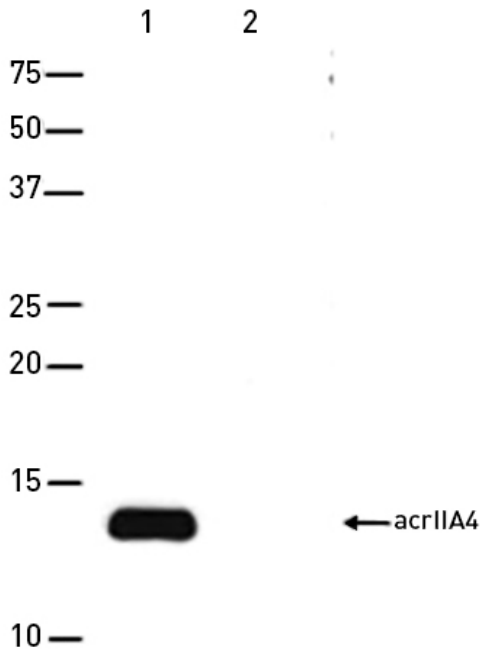


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against acrIIA4

Western blot was performed on protein extracts from U2OS cells spiked with the full length recombinant acrIIA4 protein (lane 1), using the Diagenode antibody against acrIIA4 (Cat. No. C15200248), diluted 1:1,000 in PBS-T containing 0.5% NFD. Lane 2 shows the result of U2OS protein extracts only, used as negative control. The marker is shown on the left, the position of the acrIIA4 protein is indicated on the right.