

CRISPR/Cas9-HRP antibody

Cat. No. C15200215

Type: Monoclonal

Source: Mouse

Lot: 001

Size: 500 µl

Concentration: not determined

Specificity: Streptococcus pyogenes

Purity: Protein G purified monoclonal antibody in Guardian Peroxidase Conjugate Stabilizer (Thermo Scientific cat. No.37548)

Storage: Store at 4°C.

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

Description: Monoclonal antibody raised in mouse against the Cas9 nuclease (CRISPR-associated protein 9) coupled to HRP.

Applications

Applications	Suggested dilution	References
Western blotting	1:100	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, UniProtKB/Swiss-Prot entry Q99ZW2) 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 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

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: December, 2014

Results

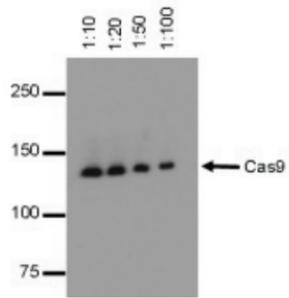


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against Cas9

Western blot was performed on protein extracts from HeLa cells transfected with a FLAG-tagged Cas9 using the Diagenode antibody against Cas9 (cat. No. C15200215). The antibody was used at different dilutions. The marker is shown on the left, the position of the FLAG-tagged Cas9 protein is indicated on the right.