

FLAG monoclonal antibody - Classic

Cat. No. C15200209

Type: Monoclonal

Source: Mouse

Lot #: 001

Size: 100 µg

Concentration: 1.0 µg/µl

Specificity: Carboxy and amino terminal linked FLAG™ tagged recombinant proteins

Purity: Purified by Protein A 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
ELISA	1:20,000 - 1:100,000	
Western blot	1:2,000 - 1:20,000	Figure 1, 2

Target description

Antibody for the detection of FLAG™ recognizes FLAG™ and is optimally suited for monitoring the expression of FLAG™ tagged fusion proteins. The antibody recognizes the epitope tag fused to either the amino- or carboxy- termini of targeted proteins. The epitope tag peptide sequence was first derived from the 11-amino-acid leader peptide of the gene-10 product from bacteriophage T7. DYKDDDDK is the most commonly used hydrophilic octapeptide tag.

Results

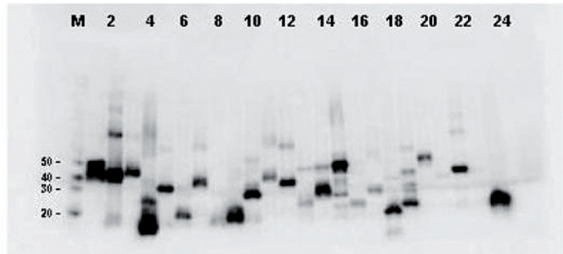


Figure 1. FLAG antibody western blot results

Twenty-four (24) clones were randomly selected and grown up from glycerol stocks by inoculating 0.5mL 2xYT medium. Expression of recombinant proteins was induced by the addition of IPTG. Proteins were purified by nickel affinity chromatography and eluted in 40 μ L. Samples were diluted 10-fold, transferred to nitrocellulose membrane and WB was performed with the FLAG monoclonal antibody.

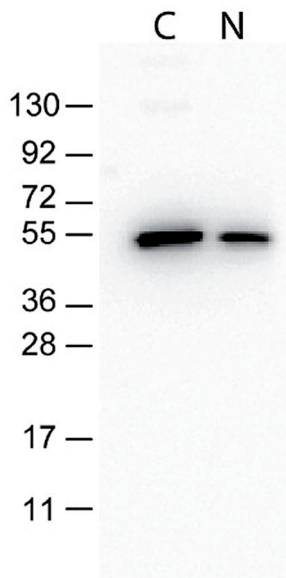


Figure 2. FLAG antibody western blot results

Monoclonal Antibody to detect FLAG™ conjugated proteins detects both C terminal linked and N terminal linked FLAG™ tagged recombinant proteins by western blot.

Diagenode sa. BELGIUM | EUROPE

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

Diagenode Inc. 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: September 10, 2014