

HA tag monoclonal antibody

Cat. No. C15200190 (MAb-190-050)

Type: Monoclonal

Isotype: IgG3

Source: Mouse

Lot #: 002

Size: 50 µg/ 25 µl

Concentration: 2 µg/µl

Specificity: The HA monoclonal is a multi-purpose antibody that recognizes HA epitopes located in the middle of protein sequences as well as at the N- or C-terminus.

Sensitivity: As low as 1 ng of HA-tagged protein can readily be detected In Western Blot.

Purity: Protein A purified monoclonal antibody in PBS (pH 7.2).

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
Western blotting	1:1,000 - 1:3,000	Fig 1
IP/ChIP	1:150 - 1:200	
IS	1:500 - 1:1,000	

*Please note that optimal antibody amount should be determined by the end-user for each specific assay condition.

Product description

Monoclonal antibody raised in mouse against the HA tag (amino acid sequence YPYDVPDYA) using a KLH-conjugated synthetic peptide. The HA antibody recognizes the influenza Hemagglutinin epitope which is extensively used as a protein tag in expression vectors. This antibody is extremely specific and allows unambiguous identification and quantitative analysis of HA tagged proteins.

Results

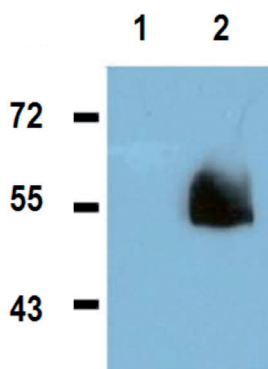


Figure 1. Western blot analysis using the Diagenode monoclonal antibody against the HA protein tag

Western blot was performed on whole cell extracts from HEK293 cells transfected with an HA-tagged expression vector using the Diagenode monoclonal antibody directed against HA (Cat. No. MAb-190-050). The antibody was used at a dilution of 1:1,000. Figure 1 shows the results for untransfected cells, used as a negative control (lane 1) and for cells transfected with a HA-tagged protein (lane 2). The MW marker (in kDa) is shown on the left. A single band is clearly visible in lane 2, but absent in lane 1 showing the high specificity of the HA antibody.

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: June 19, 2012