

## HA tag monoclonal antibody - Classic

Cat. No. C15200190

Type: Monoclonal	Specificity: Specific for HA-tagged fusion proteins expressed in bacteria, insect, mammalian cells
Size: 50 µg	Isotype: NA
Concentration: 1.0 µg/µl	Source: Mouse
Lot No.: 003	Purity: Protein A purified
Storage buffer: NA	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.	

### 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.

### Applications

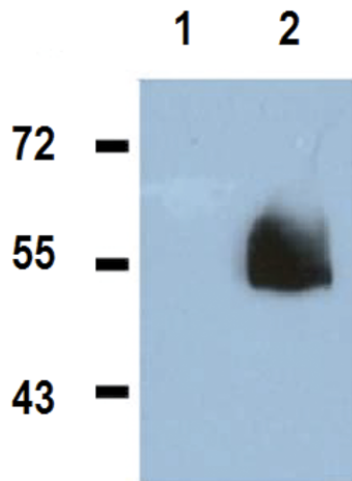
Applications	Suggested dilution *	References
Western Blotting	1:1,000 - 1:3,000	Fig 1
IP/ChIP	1:150 - 1:200	
IS	1:500 - 1:1,000	

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

### Target 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.

**Validation Data**



**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.