

Anti-HA

Mouse Monoclonal Antibody

Catalog # H98-61M

Lot # L340-1

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the HA protein

Cross Reactivity

- All proteins with the HA tag

Host/Isotype/Clone#

Mouse, IgG3, C10

Immunogen

YPYDVPDYA peptide conjugated to KLH

Formulation

10mM PBS pH 7.2

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

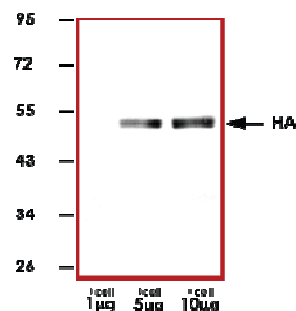
Scientific Background

The HA epitope tag (YPYDVPDYA) consists of 9 amino acid residues and comes from the Human influenza virus hemagglutinin [1]. Hemagglutinin is a viral cell surface glycoprotein responsible for the binding of viral particles to their target cell, and is named because of its ability to agglutinate erythrocytes in vitro [2].

References

1. Field, J. et al: Purification of a RAS-responsive adenyl cyclase complex from *Saccharomyces cerevisiae* by use of an epitope addition method. *Mol. Cell.* 1988 *Biol.* 8: 2159-2165
2. Medeiros, R, et al., Hemagglutinin residues of recent human A(H3N2) influenza viruses that contribute to the inability to agglutinate chicken erythrocytes. *Virology.* 2001; Oct. 10; 289(1): 74-85

Sample Data



Representative western blot with Anti-HA (1:2000) using 1 µg, 5 µg, and 10 µg of HEK-293 cell lysate transfected with an HA tagged fusion protein.

Anti-HA

Mouse Monoclonal Antibody

Catalog Number

H98-61M

Specific Lot Number

L340-1

Purification

Affinity purified using Protein A

Format

100 µg in 100 µl

Concentration

1.0 µg/µl

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.