

## Anti-HA

Rabbit Polyclonal Antibody

**Catalog # H98-63R**

Lot # Q356-1

### Cited Applications

- Western blot (1:2000) - Immunostaining
- ELISA - Immunoprecipitation

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

### Specificity

Recognizes the HA protein tag

### Cross Reactivity

- All proteins with the HA tag

### Host

Rabbit

### Immunogen

YPYDVPDYA peptide conjugated to KLH

### Formulation

TBS pH 7.2, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

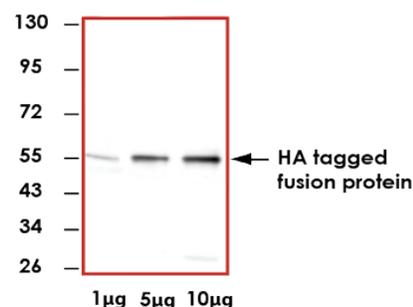
### Scientific Background

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

Rabbit Polyclonal Antibody

Catalog Number

H98-63R

Specific Lot Number

Q356-1

Purification

Affinity purified using the HA antigen

Concentration

0.25 µg/µL

Stability

1 yr 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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