

## Datasheet for Rabbit anti HA-Tag pAb

**Catalog number:** pAb-00008-100

**Background:** Protein tags are peptide sequences genetically grafted onto a recombinant protein. Often these tags are removable by chemical agents or by enzymatic means, such as proteolysis or intein splicing. Tags are attached to proteins for various purposes. Epitope tags are short peptide sequences which are chosen because high-affinity antibodies can be reliably produced in many different species. These are usually derived from viral genes, which explain their high immunoreactivity. Epitope tags include V5-tag, Myc-tag, HA-tag and NE-tag. These tags are particularly useful for western blotting, immunofluorescence and immunoprecipitation experiments, although they also find use in antibody purification.

### ■ Product Information

**Symbol:** HA  
**Alias:** HA;HA tag;HA-tag  
**Cross reactivity:** Species independent  
**Expected Protein Weight:**  
**Speices:**  
**Source:** Rabbit  
**Isotype:** IgG  
**Purity:** Affinity purification  
**Storage:** Store at -20°C. Avoid freeze / thaw cycles. | Buffer: PBS with 0.09% Sodium azide, 50% glycerol, pH7.3.

### ■ Immunogen Information

**Gene ID:**  
**Swiss:**  
**Immunogen:** A synthetic peptide corresponding to HA-Tag.

### ■ Recommended Dilutions

**WB** 1:2000 - 1:6000  
**IF** 1:50 - 1:200

**Citation of product:** If use of this item results in a publication, please use this information: Rabbit anti HA-Tag pAb(pAb-00008-100; Genecopoeia, Inc, Rockville, MD).

GeneCopoeia Products are for Research Use Only  
9620 Medical Center Drive, #101 Rockville, MD 20850 USA  
Tel: 301-762-0888; Fax: 301-762-3888  
Email: [support@genecopoeia.com](mailto:support@genecopoeia.com); Web: [www.genecopoeia.com](http://www.genecopoeia.com)

Copyright © 2024 GeneCopoeia, Inc.