

## Datasheet for S-Tag Rabbit pAb

**Catalog number:** pAb-00003-100

**Background:** S tag is an epitope tag composed of a 15 residue peptide, KETAAKFERQHMDs, derived from the pancreatic ribonuclease A.

### ■ Product Information

**Symbol:** S-Tag  
**Alias:** S-Tag; S  
**Cross reactivity:** Species independent  
**Expected Protein Weight:**  
**Species:**  
**Source:** Rabbit  
**Isotype:** IgG  
**Purity:** Affinity purification  
**Storage:** Store at -20°C. Avoid freeze / thaw cycles. | Buffer: PBS with 0.05% proclin300, 50% glycerol, pH7.3.

### ■ Immunogen Information

**Gene ID:**  
**Swiss:**  
**Immunogen:** A synthesized peptide derived from S-Tag.

### ■ Recommended Dilutions

**WB** 1:1000 - 1:5000  
**ELISA** Recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.

**Citation of product:** If use of this item results in a publication, please use this information: S-Tag Rabbit pAb(pAb-00003-100; 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.