

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; Web: www.genecopoeia.com

Copyright © 2024 GeneCopoeia, Inc.