

Datasheet for Rabbit anti His-Tag pAb

Catalog number: pAb-00009-100

Background: Consecutive histidine residues (usually 6 to 10 in length) are often inserted into the amino acid sequences of recombinant proteins. The resulting His-tagged proteins can be detected by using His antibodies.

■ Product Information

Symbol: His
Alias: His;His tag;His-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 His tag.

■ Recommended Dilutions

WB 1:500 - 1:1000
IF/ICC 1:50 - 1:200
IP 0.5µg-4µg antibody for 200µg-400µg extracts of whole cells

Citation of product: If use of this item results in a publication, please use this information: Rabbit anti His-Tag pAb(pAb-00009-100; 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.