

Datasheet for Rabbit anti V5-Tag pAb

Catalog number: pAb-00010-100

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: V5
Alias: V5;V5 tag;V5-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 0.01% thimerosal, 50% glycerol, pH7.3.

■ Immunogen Information

Gene ID:
Swiss:
Immunogen: A synthetic peptide corresponding to V5 tag.

■ Recommended Dilutions

WB 1:1000 - 1:2000

Citation of product: If use of this item results in a publication, please use this information: Rabbit anti V5-Tag pAb(pAb-00010-100; 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.