

Datasheet for Myc-Tag Rabbit mAb

Catalog number: mAb-02420-020

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: Myc
Alias: Myc; Myc tag; Myc-tag; Myc-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 0.05% proclin300, 0.05% BSA, 50% glycerol, pH7.3.

■ Immunogen Information

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

■ Recommended Dilutions

WB 1:10000 - 1:120000
IF/ICC 1:50 - 1:200
IP 0.5µg-4µg antibody for 200µg-400µg extracts of whole cells
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: Myc-Tag Rabbit mAb(mAb-02420-020; 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.