

## Datasheet for Magnetic Beads-conjugated Mouse anti DDDDK-Tag mAb

**Catalog number:** mAb-02419-100

**Background:** FLAG-tag, or FLAG octapeptide, or FLAG epitope, is a polypeptide protein tag that can be added to a protein using recombinant DNA technology, having the sequence motif DYKDDDDK. It has been used for studying proteins in living cells and for protein purification by affinity chromatography. It has been used to separate recombinant, overexpressed protein from wild-type protein expressed by the host organism. It can also be used in the isolation of protein complexes with multiple subunits, because its mild purification procedure tends not to disrupt such complexes. It has been used to obtain proteins of sufficient purity and quality to carry out 3D structure determination by x-ray crystallography. A FLAG-tag can be used in many different assays that require recognition by an antibody. If there is no antibody against a given protein, adding a FLAG-tag to a protein allows the protein to be studied with an antibody against the FLAG sequence. Examples are cellular localization studies by immunofluorescence or detection by SDS PAGE protein electrophoresis and Western blotting.

### ■ Product Information

**Symbol:** DDDDK  
**Alias:** DDDDK;DDDDK tag;DDDDK-tag  
**Cross reactivity:** Species independent  
**Expected Protein Weight:**  
**Species:**  
**Source:** Mouse  
**Isotype:** IgG1  
**Purity:** Affinity purification  
**Storage:** Store at 4°C Avoid freeze / thaw cycles. | Buffer: PBS with 0.05% proclin300, pH7.3.

### ■ Immunogen Information

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

### ■ Recommended Dilutions

**IP** 30ul antibody (bead slurry) for 200µg-400µg extracts of whole cells

**Citation of product:** If use of this item results in a publication, please use this information:  
Magnetic Beads-conjugated Mouse anti DDDDK-Tag mAb(mAb-02419-100; 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.