

## Datasheet for Rabbit anti MBP-Tag pAb

**Catalog number:** pAb-00012-100

**Background:** Biofilm microarray result was validated with RT-PCR. [More information is available at EcoGene: EG10554]. In addition to being the periplasmic substrate-binding component of the maltose ABC transporter, MalE with bound substrate is also able to bind to the chemoreceptor Tar to induce chemotaxis toward maltose.

### ■ Product Information

**Symbol:**

**Alias:** ECK4026

**Cross reactivity:** Species independent

**Expected Protein Weight:** 43kDa

**Speices:**

**Source:** Rabbit

**Isotype:** IgG

**Purity:** Affinity purification

**Storage:** Store at -20°C. Avoid freeze / thaw cycles.[Buffer: PBS with 0.09% sodium azide,0.05% BSA,50% glycerol,pH7.3.

### ■ Immunogen Information

**Gene ID:** 948538

**Swiss:** P0AEX9

**Immunogen:** Recombinant protein containing a sequence corresponding to amino acids 27-396 of MBP protein (NP\_418458).

### ■ Recommended Dilutions

**WB** 1:10000 - 1:60000

**Citation of product:** If use of this item results in a publication, please use this information: Rabbit anti MBP-Tag pAb(pAb-00012-100; 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.