

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; Web:www.genecopoeia.com

Copyright © 2024 GeneCopoeia, Inc.